



2025 Price List

Products for DNA research exclusively
distributed by Eurogentec



Description	Reference	Scale/Quantity	Price List 2025				
			EUR	CHF	GBP	USD	
Glen Research products for oligonucleotide synthesis							EA1
GlenDNA							EA1A1Z
1c-dC-CE Phosphoramidite	10-1015-40	4 g	62.00	59.50	53.30	-	
1-Me-A-CE Phosphoramidite	10-3501-02	0.25 g	1,474.00	1,415.00	1,268.00	-	
1-Me-A-CE Phosphoramidite	10-3501-95	50 µmol	302.00	290.00	260.00	-	
3'-Phosphate CPG (High Load)	25-2900-46	4 x 2.5 µmol	173.00	166.00	149.00	-	
5% Phenoxyacetic Anhydride in Acetonitrile	40-4128-71	1000 mL	774.00	743.00	666.00	-	
5-I-U-CE Phosphoramidite	10-3091-95	50 µmol	194.00	186.00	167.00	-	
5-Me-dC Brancher Phosphoramidite	10-1018-02	0.25 g	726.00	697.00	624.00	-	
5-Me-dC Brancher Phosphoramidite	10-1018-90	100 µmol	295.00	283.00	254.00	-	
5-Me-dC Brancher Phosphoramidite, Expedite	10-1018-02E	0.25 g	726.00	697.00	624.00	-	
5-Me-U-CE Phosphoramidite	10-3050-95	50 µmol	137.00	132.00	118.00	-	
AAA Trimer Phosphoramidite	13-1000-90	100 µmol	1,294.00	1,242.00	1,113.00	-	
AAA Trimer Phosphoramidite	13-1000-90M	100 µmol	1,294.00	1,242.00	1,113.00	-	
AAA Trimer Phosphoramidite	13-1000-95	50 µmol	669.00	642.00	575.00	-	
AAA Trimer Phosphoramidite	13-1000-95M	50 µmol	669.00	642.00	575.00	-	
AAC Trimer Phosphoramidite	13-1001-90	100 µmol	1,294.00	1,242.00	1,113.00	-	
AAC Trimer Phosphoramidite	13-1001-90M	100 µmol	1,294.00	1,242.00	1,113.00	-	
AAC Trimer Phosphoramidite	13-1001-95	50 µmol	669.00	642.00	575.00	-	
AAC Trimer Phosphoramidite	13-1001-95M	50 µmol	669.00	642.00	575.00	-	
AB3900 Ac-dC-CPG 1000A, 1.0 µmol	20-2115-61	200 pce	968.00	929.00	832.00	-	
AB3900 dmf-dG-CPG 1000A, 1.0 µmol	20-2129-61	200 pce	968.00	929.00	832.00	-	
AB3900 dt-CPG 1000A, 1.0 µmol	20-2131-61	200 pce	968.00	929.00	832.00	-	
ACC Trimer Phosphoramidite	13-1011-95M	50 µmol	669.00	642.00	575.00	-	
ACC Trimer Phosphoramidite	20-2131-61	100 µmol	968.00	929.00	832.00	-	
ACC Trimer Phosphoramidite, Expedite	13-1011-95E	50 µmol	669.00	642.00	575.00	-	
Ac-dC-CE Phosphoramidite	10-1015-02	0.25 g	4.96	4.76	4.27	-	
Ac-dC-CE Phosphoramidite	10-1015-05	0.5 g	9.29	8.92	7.99	-	
Ac-dC-CE Phosphoramidite	10-1015-05M	0.5 g	9.29	8.92	7.99	-	
Ac-dC-CE Phosphoramidite	10-1015-10	1 g	15.50	14.90	13.30	-	
Ac-dC-CE Phosphoramidite	10-1015-10M	1 g	15.50	14.90	13.30	-	
Ac-dC-CE Phosphoramidite	10-1015-1C	1 g	15.50	14.90	13.30	-	
Ac-dC-CE Phosphoramidite	10-1015-1S	10 g	155.00	149.00	133.00	-	
Ac-dC-CE Phosphoramidite	10-1015-20	2 g	31.00	29.80	26.70	-	
Ac-dC-CE Phosphoramidite	10-1015-2C	2 g	31.00	29.80	26.70	-	
Ac-dC-CE Phosphoramidite	10-1015-50	5 g	77.40	74.30	66.60	-	
Ac-dC-CE Phosphoramidite	10-1015-C2	0.25 g	4.96	4.76	4.27	-	
Ac-dC-CE Phosphoramidite	10-1015-C5	0.5 g	9.29	8.92	7.99	-	
Ac-dC-CE Phosphoramidite,	10-1015-02M	0.25 g	4.96	4.76	4.27	-	
Ac-dC-CE Phosphoramidite, Expedite	10-1015-02E	0.25 g	4.96	4.76	4.27	-	
Ac-dC-CE Phosphoramidite, Pharmacia	10-1015-02P	0.25 g	4.96	4.76	4.27	-	
Ac-dC-CE Phosphoramidite, Pharmacia	10-1015-05P	0.5 g	9.29	8.92	7.99	-	
Ac-dC-CPG 1000	20-2015-01	0.1 g	13.30	12.80	11.40	-	
Ac-dC-CPG 1000	20-2015-02	0.25 g	17.70	17.00	15.20	-	
Ac-dC-CPG 1000	20-2015-10	1 g	66.40	63.70	57.10	-	
Ac-dC-CPG 1000	20-2015-61	48 x 1 µmol	265.00	254.00	228.00	-	
Ac-dC-CPG 1000	20-2015-62	200 x 0.2 µmol	664.00	637.00	571.00	-	
Ac-dC-CPG 1000	20-2115-13	1 x 10 µmol	88.50	85.00	76.10	-	
Ac-dC-CPG 1000	20-2115-41	4 x 1 µmol	31.00	29.80	26.70	-	
Ac-dC-CPG 1000	20-2115-42	4 x 0.2 µmol	31.00	29.80	26.70	-	
Ac-dC-CPG 1000	20-2115-45	4 x 40 nmol	31.00	29.80	26.70	-	
Ac-dC-CPG 1000	20-2115-62	200 x 200 nmol	830.00	797.00	714.00	-	
Ac-dC-CPG 1000	20-2215-14	1 x 15 µmol	133.00	128.00	114.00	-	
Ac-dC-CPG 1000	20-2215-41	4 x 1 µmol	22.10	21.20	19.00	-	
Ac-dC-CPG 1000	20-2215-42	4 x 0.2 µmol	15.50	14.90	13.30	-	
Ac-dC-CPG 1000	20-2215-45	4 x 40 nmol	15.50	14.90	13.30	-	
Ac-dC-CPG 1000 (1 g)	20-2015-10D	5 g	66.40	63.70	57.10	-	
Ac-dC-CPG 500	20-2013-01	0.1 g	13.30	12.80	11.40	-	
Ac-dC-CPG 500	20-2013-02	0.25 g	17.70	17.00	15.20	-	
Ac-dC-CPG 500	20-2113-13	1 x 10 µmol	88.50	85.00	76.10	-	
Ac-dC-CPG 500	20-2113-41	4 x 1 µmol	31.00	29.80	26.70	-	
Ac-dC-CPG 500	20-2113-42	4 x 0.2 µmol	31.00	29.80	26.70	-	
Ac-dC-CPG 500	20-2213-14	1 x 15 µmol	133.00	128.00	114.00	-	
Ac-dC-CPG 500	20-2213-41	4 x 1 µmol	22.10	21.20	19.00	-	
Ac-dC-CPG 500	20-2213-42	4 x 0.2 µmol	15.50	14.90	13.30	-	
Ac-G-CE Phosphoramidite	10-3025-10	1 g	99.60	95.60	85.70	-	
Ac-G-CE Phosphoramidite,	10-3025-02	0.25 g	35.40	34.00	30.40	-	
ACT Trimer Phosphoramidite	13-1013-90	100 µmol	1,294.00	1,242.00	1,113.00	-	
ACT Trimer Phosphoramidite	13-1013-90M	100 µmol	1,294.00	1,242.00	1,113.00	-	
ACT Trimer Phosphoramidite	13-1013-95	50 µmol	669.00	642.00	575.00	-	
ACT Trimer Phosphoramidite	13-1013-95M	50 µmol	669.00	642.00	575.00	-	
AGA Trimer Phosphoramidite	13-1020-90M	100 µmol	1,294.00	1,242.00	1,113.00	-	
AGA Trimer Phosphoramidite	13-1020-95M	50 µmol	669.00	642.00	575.00	-	
ATC Trimer Phosphoramidite	13-1031-90	100 µmol	1,294.00	1,242.00	1,113.00	-	
ATC Trimer Phosphoramidite	13-1031-90M	100 µmol	1,294.00	1,242.00	1,113.00	-	
ATC Trimer Phosphoramidite	13-1031-95	50 µmol	669.00	642.00	575.00	-	
ATC Trimer Phosphoramidite	13-1031-95M	50 µmol	669.00	642.00	575.00	-	
ATG Trimer Phosphoramidite	13-1032-90	100 µmol	1,294.00	1,242.00	1,113.00	-	
ATG Trimer Phosphoramidite	13-1032-90M	100 µmol	1,294.00	1,242.00	1,113.00	-	
ATG Trimer Phosphoramidite	13-1032-95	50 µmol	669.00	642.00	575.00	-	
ATG Trimer Phosphoramidite	13-1032-95M	50 µmol	669.00	642.00	575.00	-	
Br-U-CE Phosphoramidite	10-3090-95	50 µmol	180.00	173.00	155.00	-	
CAG Trimer Phosphoramidite	13-1102-90	100 µmol	1,294.00	1,242.00	1,113.00	-	
CAG Trimer Phosphoramidite	13-1102-90M	100 µmol	1,294.00	1,242.00	1,113.00	-	

Description	Reference	Scale/Quantity	Price List 2025			
			EUR	CHF	GBP	USD
CAG Trimer Phosphoramidite	13-1102-95	50 µmol	669.00	642.00	575.00	-
CAG Trimer Phosphoramidite	13-1102-95M	50 µmol	669.00	642.00	575.00	-
CAT Trimer Phosphoramidite	13-1103-90	100 µmol	1,294.00	1,242.00	1,113.00	-
CAT Trimer Phosphoramidite	13-1103-90M	100 µmol	1,294.00	1,242.00	1,113.00	-
CAT Trimer Phosphoramidite	13-1103-95	50 µmol	669.00	642.00	575.00	-
CAT Trimer Phosphoramidite	13-1103-95M	50 µmol	669.00	642.00	575.00	-
CAT Trimer Phosphoramidite, Expedite	13-1103-95E	50 µmol	669.00	642.00	575.00	-
CCA Trimer Phosphoramidite	13-1110-90M	100 µmol	1,294.00	1,242.00	1,113.00	-
CCA Trimer Phosphoramidite	13-1110-95M	50 µmol	669.00	642.00	575.00	-
CCG Trimer Phosphoramidite	13-1112-90	100 µmol	1,294.00	1,242.00	1,113.00	-
CCG Trimer Phosphoramidite	13-1112-90M	100 µmol	1,294.00	1,242.00	1,113.00	-
CCG Trimer Phosphoramidite	13-1112-95	50 µmol	669.00	642.00	575.00	-
CCG Trimer Phosphoramidite	13-1112-95M	50 µmol	669.00	642.00	575.00	-
CGG Trimer Phosphoramidite	13-1122-90M	100 µmol	1,294.00	1,242.00	1,113.00	-
CGG Trimer Phosphoramidite	13-1122-95M	50 µmol	669.00	642.00	575.00	-
CGT Trimer Phosphoramidite	13-1123-90	100 µmol	1,294.00	1,242.00	1,113.00	-
CGT Trimer Phosphoramidite	13-1123-90M	100 µmol	1,294.00	1,242.00	1,113.00	-
CGT Trimer Phosphoramidite	13-1123-95	50 µmol	669.00	642.00	575.00	-
CGT Trimer Phosphoramidite	13-1123-95M	50 µmol	669.00	642.00	575.00	-
Colonne dA - 10 µmol	20-2100-13F	10 nmol	88.50	85.00	76.10	-
Colonne dA - 2.5 µmol	20-2113-13F	2.5 nmol	88.50	85.00	76.10	-
Colonne dG - 10 µmol	20-2120-13F	10 nmol	88.50	85.00	76.10	-
Colonne dT - 10 µmol	20-2130-13F	10 nmol	88.50	85.00	76.10	-
CTG Trimer Phosphoramidite	13-1132-90	100 µmol	1,294.00	1,242.00	1,113.00	-
CTG Trimer Phosphoramidite	13-1132-90M	100 µmol	1,294.00	1,242.00	1,113.00	-
CTG Trimer Phosphoramidite	13-1132-95	50 µmol	669.00	642.00	575.00	-
CTG Trimer Phosphoramidite	13-1132-95M	50 µmol	669.00	642.00	575.00	-
dA LV-PS	26-2100-42	4 x 0.2 µmol	43.60	43.60	39.20	-
dA LV-PS	26-2100-45	4 x 40 nmol	43.60	43.60	39.20	-
dA PS	26-2600-62	200 x 200 nmol	653.00	627.00	562.00	-
dA PS	26-2600-65	200 x 40 nmol	653.00	627.00	562.00	-
dA, dC, dG, dT-CPG 1000	20-2241-41	4 x 1 µmol	22.10	21.20	19.00	-
dA, dC, dG, dT-CPG 1000	20-2241-42	4 x 0.2 µmol	15.50	14.90	13.30	-
dA, dC, dG, dT-CPG 1000	20-2241-45	4 x 40 nmol	15.50	14.90	13.30	-
dA, dC, dG, dT CPG 1000	20-2141-41	4 x 1 µmol	31.00	29.80	26.70	-
dA, dC, dG, dT CPG 1000	20-2141-42	4 x 0.2 µmol	31.00	29.80	26.70	-
dA, dC, dG, dT CPG 1000	20-2141-45	4 x 40 nmol	31.00	29.80	26.70	-
dA, dC, dG, dT- LV-PS 0.2µm Columns	26-2140-42	4 x 0.2 µmol	43.60	43.60	39.20	-
dA, dC, dG, dT- LV-PS 40nm Columns	26-2140-45	4 x 40 nmol	43.60	43.60	39.20	-
dA, dC, dG, dT-CPG 2000	20-2242-42	4 x 0.2 µmol	15.50	14.90	13.30	-
dA, dC, dG, dT-CPG 500	20-2140-41	4 x 1 µmol	31.00	29.80	26.70	-
dA, dC, dG, dT-CPG 500	20-2140-42	4 x 0.2 µmol	31.00	29.80	26.70	-
dA, dC, dG, dT-CPG 500	20-2240-41	4 x 1 µmol	22.10	21.20	19.00	-
dA, dC, dG, dT-CPG 500	20-2240-42	4 x 0.2 µmol	15.50	14.90	13.30	-
dA-CE Phosphoramidite	10-1000-02	0.25 g	4.96	4.76	4.27	-
dA-CE Phosphoramidite	10-1000-05	0.5 g	9.29	8.92	7.99	-
dA-CE Phosphoramidite	10-1000-10	1 g	15.50	14.90	13.30	-
dA-CE Phosphoramidite	10-1000-1C	1 g	15.50	14.90	13.30	-
dA-CE Phosphoramidite	10-1000-1S	10 g	155.00	149.00	133.00	-
dA-CE Phosphoramidite	10-1000-20	2 g	31.00	29.80	26.70	-
dA-CE Phosphoramidite	10-1000-2C	2 g	31.00	29.80	26.70	-
dA-CE Phosphoramidite	10-1000-40	4 g	62.00	59.50	53.30	-
dA-CE Phosphoramidite	10-1000-50	5 g	77.40	74.30	66.60	-
dA-CE Phosphoramidite	10-1000-5S	5 g	77.40	74.30	66.60	-
dA-CE Phosphoramidite	10-1000-C2	0.25 g	4.96	4.76	4.27	-
dA-CE Phosphoramidite	10-1000-C5	0.5 g	9.29	8.92	7.99	-
dA-CE Phosphoramidite, Mermade	10-1000-02M	0.25 g	4.96	4.76	4.27	-
dA-CE Phosphoramidite, Mermade	10-1000-05M	0.5 g	9.29	8.92	7.99	-
dA-CE Phosphoramidite, Mermade	10-1000-10M	1 g	15.50	14.90	13.30	-
dA-CPG 1 µmol columns	20-2101-61	200 pce	968.00	929.00	832.00	-
dA-CPG 1000	20-2001-01	0.1 g	13.30	12.80	11.40	-
dA-CPG 1000	20-2001-02	0.25 g	17.70	17.00	15.20	-
dA-CPG 1000	20-2001-10	1 g	66.40	63.70	57.10	-
dA-CPG 1000	20-2001-61	48 x 1 µmol	265.00	254.00	228.00	-
dA-CPG 1000	20-2001-62	200 x 0.2 µmol	664.00	637.00	571.00	-
dA-CPG 1000	20-2101-13	1 x 10 µmol	88.50	85.00	76.10	-
dA-CPG 1000	20-2101-41	4 x 1 µmol	31.00	29.80	26.70	-
dA-CPG 1000	20-2101-42	4 x 0.2 µmol	31.00	29.80	26.70	-
dA-CPG 1000	20-2101-45	4 x 40 nmol	31.00	29.80	26.70	-
dA-CPG 1000	20-2101-62	200 x 200 nm	830.00	797.00	714.00	-
dA-CPG 1000	20-2201-14	1 x 15 µmol	133.00	128.00	114.00	-
dA-CPG 1000	20-2201-41	4 x 1 µmol	22.10	21.20	19.00	-
dA-CPG 1000	20-2201-42	4 x 0.2 µmol	15.50	14.90	13.30	-
dA-CPG 1000	20-2201-45	4 x 40 nmol	15.50	14.90	13.30	-
dA-CPG 1000, Expedite	20-2201-42E	4 x 0.2 µmol	15.50	14.90	13.30	-
dA-CPG 2000	20-2002-01	0.1 g	16.60	15.90	14.30	-
dA-CPG 2000	20-2002-02	0.25 g	33.20	31.90	28.60	-
dA-CPG 2000	20-2002-10	1 g	166.00	159.00	143.00	-
dA-CPG 2000	20-2102-42	4 x 0.2 µmol	31.00	29.80	26.70	-
dA-CPG 2000	20-2202-42	4 x 0.2 µmol	15.50	14.90	13.30	-
dA-CPG 500	20-2000-01	0.1 g	13.30	12.80	11.40	-
dA-CPG 500	20-2000-02	0.25 g	17.70	17.00	15.20	-
dA-CPG 500	20-2000-10	1 g	66.40	63.70	57.10	-
dA-CPG 500	20-2100-13	1 x 10 µmol	88.50	85.00	76.10	-

Description	Reference	Scale/Quantity	Price List 2025			
			EUR	CHF	GBP	USD
dA-CPG 500	20-2100-41	4 x 1 µmol	31.00	29.80	26.70	-
dA-CPG 500	20-2100-42	4 x 0.2 µmol	31.00	29.80	26.70	-
dA-CPG 500	20-2200-14	1 x 15 µmol	133.00	128.00	114.00	-
dA-CPG 500	20-2200-41	4 x 1 µmol	22.10	21.20	19.00	-
dA-CPG 500	20-2200-42	4 x 0.2 µmol	15.50	14.90	13.30	-
dA-High Load-CPG	25-2000-02	0.25 g	22.10	21.20	19.00	-
dA-High Load-CPG	25-2000-10	1 g	79.70	76.50	68.50	-
dA-High Load-CPG	25-2100-17	1 x 25 µmol	111.00	107.00	95.50	-
dA-High Load-CPG	25-2100-46	4 x 2.5 µmol	66.40	63.70	57.10	-
dA-High Load-CPG	25-2200-18	1 x 35 µmol	164.00	157.00	141.00	-
dA-High Load-CPG	25-2200-46	4 x 2.5 µmol	66.40	63.70	57.10	-
dA-Icaa-CPG 500 Å col.(96pcs)	20-2200-42SP	0.2 µmol	15.50	14.90	13.30	-
dA-Thiophosphoramidite	10-1700-02	0.25 g	604.00	580.00	519.00	-
dA-Thiophosphoramidite	10-1700-90	100 µmol	252.00	242.00	217.00	-
dC LV-PS	26-2110-42	4 x 0.2 µmol	43.60	43.60	39.20	-
dC LV-PS	26-2110-45	4 x 40 nmol	43.60	43.60	39.20	-
dC PS	26-2610-62	200 x 200 nmol	653.00	627.00	562.00	-
dC PS	26-2610-65	200 x 40 nmol	653.00	627.00	562.00	-
dC Thiophosphoramidite	10-1710-02	0.25 g	604.00	580.00	519.00	-
dC-CE Phosphoramidite	10-1010-02	0.25 g	4.96	4.76	4.27	-
dC-CE Phosphoramidite	10-1010-02M	0.25 g	4.96	4.76	4.27	-
dC-CE Phosphoramidite	10-1010-05	0.5 g	9.29	8.92	7.99	-
dC-CE Phosphoramidite	10-1010-05M	0.5 g	9.29	8.92	7.99	-
dC-CE Phosphoramidite	10-1010-10	1 g	15.50	14.90	13.30	-
dC-CE Phosphoramidite	10-1010-10M	1 g	15.50	14.90	13.30	-
dC-CE Phosphoramidite	10-1010-1C	1 g	15.50	14.90	13.30	-
dC-CE Phosphoramidite	10-1010-1S	10 g	155.00	149.00	133.00	-
dC-CE Phosphoramidite	10-1010-20	2 g	31.00	29.80	26.70	-
dC-CE Phosphoramidite	10-1010-2C	2 g	31.00	29.80	26.70	-
dC-CE Phosphoramidite	10-1010-40	4 g	62.00	59.50	53.30	-
dC-CE Phosphoramidite	10-1010-50	5 g	77.40	74.30	66.60	-
dC-CE Phosphoramidite	10-1010-5S	5 g	77.40	74.30	66.60	-
dC-CE Phosphoramidite	10-1010-C2	0.25 g	4.96	4.76	4.27	-
dC-CE Phosphoramidite	10-1010-C5	0.5 g	9.29	8.92	7.99	-
dC-CPG 1000	20-2011-01	0.1 g	13.30	12.80	11.40	-
dC-CPG 1000	20-2011-02	0.25 g	17.70	17.00	15.20	-
dC-CPG 1000	20-2011-10	1 g	66.40	63.70	57.10	-
dC-CPG 1000	20-2111-13	1 x 10 µmol	88.50	85.00	76.10	-
dC-CPG 1000	20-2111-41	4 x 1 µmol	31.00	29.80	26.70	-
dC-CPG 1000	20-2111-42	4 x 0.2 µmol	31.00	29.80	26.70	-
dC-CPG 1000	20-2111-45	4 x 40 nmol	31.00	29.80	26.70	-
dC-CPG 1000	20-2211-14	1 x 15 µmol	133.00	128.00	114.00	-
dC-CPG 1000	20-2211-41	4 x 1 µmol	22.10	21.20	19.00	-
dC-CPG 1000	20-2211-42	4 x 0.2 µmol	15.50	14.90	13.30	-
dC-CPG 1000	20-2211-45	4 x 40 nmol	15.50	14.90	13.30	-
dC-CPG 1000, Expedite	20-2211-42E	4 x 0.2 µmol	15.50	14.90	13.30	-
dC-CPG 2000	20-2012-01	0.1 g	16.60	15.90	14.30	-
dC-CPG 2000	20-2012-02	0.25 g	33.20	31.90	28.60	-
dC-CPG 2000	20-2012-10	1 g	166.00	159.00	143.00	-
dC-CPG 2000	20-2112-42	4 x 0.2 µmol	31.00	29.80	26.70	-
dC-CPG 2000	20-2212-42	4 x 0.2 µmol	15.50	14.90	13.30	-
dC-CPG 500	20-2010-01	0.1 g	13.30	12.80	11.40	-
dC-CPG 500	20-2010-02	0.25 g	17.70	17.00	15.20	-
dC-CPG 500	20-2010-10	1 g	66.40	63.70	57.10	-
dC-CPG 500	20-2110-13	1 x 10 µmol	88.50	85.00	76.10	-
dC-CPG 500	20-2110-41	4 x 1 µmol	31.00	29.80	26.70	-
dC-CPG 500	20-2110-42	4 x 0.2 µmol	31.00	29.80	26.70	-
dC-CPG 500	20-2210-14	1 x 15 µmol	133.00	128.00	114.00	-
dC-CPG 500	20-2210-41	4 x 1 µmol	22.10	21.20	19.00	-
dC-CPG 500	20-2210-42	4 x 0.2 µmol	15.50	14.90	13.30	-
dC-High Load-CPG	25-2010-02	0.25 g	22.10	21.20	19.00	-
dC-High Load-CPG	25-2010-10	1 g	79.70	76.50	68.50	-
dC-High Load-CPG	25-2110-17	1 x 25 µmol	111.00	107.00	95.50	-
dC-High Load-CPG	25-2110-46	4 x 2.5 µmol	66.40	63.70	57.10	-
dC-High Load-CPG	25-2210-18	1 x 35 µmol	164.00	157.00	141.00	-
dC-High Load-CPG	25-2210-46	4 x 2.5 µmol	66.40	63.70	57.10	-
dC-Icaa-CPG 500 Å -cot.(96pcs)	20-2210-42SP	0.2 µmol	15.50	14.90	13.30	-
dC-Thiophosphoramidite	10-1710-90	100 µmol	252.00	242.00	217.00	-
dG LV-PS	26-2120-42	4 x 0.2 µmol	43.60	43.60	39.20	-
dG LV-PS	26-2120-45	4 x 40 nmol	43.60	43.60	39.20	-
dG-CE Phosphoramidite	10-1020-02	0.25 g	4.96	4.76	4.27	-
dG-CE Phosphoramidite	10-1020-05	0.5 g	9.29	8.92	7.99	-
dG-CE Phosphoramidite	10-1020-05M	0.5 g	9.29	8.92	7.99	-
dG-CE Phosphoramidite	10-1020-10	1 g	15.50	14.90	13.30	-
dG-CE Phosphoramidite	10-1020-10M	1 g	15.50	14.90	13.30	-
dG-CE Phosphoramidite	10-1020-1C	1 g	15.50	14.90	13.30	-
dG-CE Phosphoramidite	10-1020-1S	10 g	155.00	149.00	133.00	-
dG-CE Phosphoramidite	10-1020-20	2 g	31.00	29.80	26.70	-
dG-CE Phosphoramidite	10-1020-2C	2 g	31.00	29.80	26.70	-
dG-CE Phosphoramidite	10-1020-40	4 g	62.00	59.50	53.30	-
dG-CE Phosphoramidite	10-1020-50	5 g	77.40	74.30	66.60	-
dG-CE Phosphoramidite	10-1020-5S	5 g	77.40	74.30	66.60	-
dG-CE Phosphoramidite	10-1020-C2	0.25 g	4.96	4.76	4.27	-
dG-CE Phosphoramidite,	10-1020-02M	0.25 g	4.96	4.76	4.27	-

Description	Reference	Scale/Quantity	Price List 2025			
			EUR	CHF	GBP	USD
dG-CE Phosphoramidite, Expedite	10-1020-C5	0.5 g	9.29	8.92	7.99	-
dG-CPG 1000	20-2021-01	0.1 g	13.30	12.80	11.40	-
dG-CPG 1000	20-2021-02	0.25 g	17.70	17.00	15.20	-
dG-CPG 1000	20-2021-10	1 g	66.40	63.70	57.10	-
dG-CPG 1000	20-2021-61	48 x 1 µmol	265.00	254.00	228.00	-
dG-CPG 1000	20-2021-62	200 x 0.2 µmol	664.00	637.00	571.00	-
dG-CPG 1000	20-2121-13	1 x 10 µmol	88.50	85.00	76.10	-
dG-CPG 1000	20-2121-41	4 x 1 µmol	31.00	29.80	26.70	-
dG-CPG 1000	20-2121-42	4 x 0.2 µmol	31.00	29.80	26.70	-
dG-CPG 1000	20-2121-45	4 x 40 nmol	31.00	29.80	26.70	-
dG-CPG 1000	20-2221-14	1 x 15 µmol	133.00	128.00	114.00	-
dG-CPG 1000	20-2221-41	4 x 1 µmol	22.10	21.20	19.00	-
dG-CPG 1000	20-2221-42	4 x 0.2 µmol	15.50	14.90	13.30	-
dG-CPG 1000	20-2221-45	4 x 40 nmol	15.50	14.90	13.30	-
dG-CPG 1000, Expedite	20-2221-42E	4 x 0.2 µmol	15.50	14.90	13.30	-
dG-CPG 2000	20-2022-01	0.1 g	16.60	15.90	14.30	-
dG-CPG 2000	20-2022-02	0.25 g	33.20	31.90	28.60	-
dG-CPG 2000	20-2022-10	1 g	166.00	159.00	143.00	-
dG-CPG 2000	20-2122-42	4 x 0.2 µmol	31.00	29.80	26.70	-
dG-CPG 2000	20-2222-42	4 x 0.2 µmol	15.50	14.90	13.30	-
dG-CPG 500	20-2020-01	0.1 g	13.30	12.80	11.40	-
dG-CPG 500	20-2020-02	0.25 g	17.70	17.00	15.20	-
dG-CPG 500	20-2020-10	1 g	66.40	63.70	57.10	-
dG-CPG 500	20-2120-13	1 x 10 µmol	88.50	85.00	76.10	-
dG-CPG 500	20-2120-41	4 x 1 µmol	31.00	29.80	26.70	-
dG-CPG 500	20-2120-42	4 x 0.2 µmol	31.00	29.80	26.70	-
dG-CPG 500	20-2220-14	1 x 15 µmol	133.00	128.00	114.00	-
dG-CPG 500	20-2220-41	4 x 1 µmol	22.10	21.20	19.00	-
dG-CPG 500	20-2220-42	4 x 0.2 µmol	15.50	14.90	13.30	-
dG-High Load CPG	25-2020-02	0.25 g	22.10	21.20	19.00	-
dG-High Load CPG	25-2020-10	1 g	79.70	76.50	68.50	-
dG-High Load CPG	25-2120-17	1 x 25 µmol	111.00	107.00	95.50	-
dG-High Load CPG	25-2120-46	4 x 2.5 µmol	66.40	63.70	57.10	-
dG-High Load CPG	25-2120-46M	4 x 2.5 µmol	66.40	63.70	57.10	-
dG-High Load CPG	25-2220-18	1 x 35 µmol	164.00	157.00	141.00	-
dG-High Load CPG	25-2220-46	4 x 2.5 µmol	66.40	63.70	57.10	-
dG-lcaa-CPG 1000 Å	20-2121-41P	1 µmol	31.00	29.80	26.70	-
dG-lcaa-CPG 500 Å - col.(96pcs)	20-2220-42SP	0.2 µmol	15.50	14.90	13.30	-
dG-Thiophosphoramidite	10-1720-90	100 µmol	252.00	242.00	217.00	-
dG-Thiophosphoramidite,	10-1720-02	0.25 g	604.00	580.00	519.00	-
dmf-dG-CE Phosphoramidite	10-1029-05M	0.5 g	9.29	8.92	7.99	-
dmf-dG-CE Phosphoramidite	10-1029-10M	1 g	15.50	14.90	13.30	-
dmf-dG-CE Phosphoramidite	10-1029-2C	2 g	31.00	29.80	26.70	-
dmf-dG-CE Phosphoramidite	10-1029-40	4 g	62.00	59.50	53.30	-
dmf-dG-CE Phosphoramidite,	10-1029-02M	0.25 g	4.96	4.76	4.27	-
dmf-dG-CPG	20-2129-62	200 x 200 nm	830.00	797.00	714.00	-
dT LV-PS	26-2130-42	4 x 0.2 µmol	43.60	43.60	39.20	-
dT LV-PS	26-2130-45	4 x 40 nmol	43.60	43.60	39.20	-
dT-CE Phosphoramidite	10-1030-02	0.25 g	4.96	4.76	4.27	-
dT-CE Phosphoramidite	10-1030-05	0.5 g	9.29	8.92	7.99	-
dT-CE Phosphoramidite	10-1030-05M	0.5 g	9.29	8.92	7.99	-
dT-CE Phosphoramidite	10-1030-10	1 g	15.50	14.90	13.30	-
dT-CE Phosphoramidite	10-1030-10M	1 g	15.50	14.90	13.30	-
dT-CE Phosphoramidite	10-1030-1C	1 g	15.50	14.90	13.30	-
dT-CE Phosphoramidite	10-1030-1S	10 g	155.00	149.00	133.00	-
dT-CE Phosphoramidite	10-1030-20	2 g	31.00	29.80	26.70	-
dT-CE Phosphoramidite	10-1030-2C	2 g	31.00	29.80	26.70	-
dT-CE Phosphoramidite	10-1030-40	4 g	62.00	59.50	53.30	-
dT-CE Phosphoramidite	10-1030-50	5 g	77.40	74.30	66.60	-
dT-CE Phosphoramidite	10-1030-5S	5 g	77.40	74.30	66.60	-
dT-CE Phosphoramidite	10-1030-C2	0.25 g	4.96	4.76	4.27	-
dT-CE Phosphoramidite	10-1030-C5	0.5 g	9.29	8.92	7.99	-
dT-CE Phosphoramidite,	10-1030-02M	0.25 g	4.96	4.76	4.27	-
dT-CPG 1000	20-2031-01	0.1 g	13.30	12.80	11.40	-
dT-CPG 1000	20-2031-02	0.25 g	17.70	17.00	15.20	-
dT-CPG 1000	20-2031-10	1 g	66.40	63.70	57.10	-
dT-CPG 1000	20-2031-61	48 x 1 µmol	265.00	254.00	228.00	-
dT-CPG 1000	20-2031-62	200 x 0.2 µmol	664.00	637.00	571.00	-
dT-CPG 1000	20-2131-13	1 x 10 µmol	88.50	85.00	76.10	-
dT-CPG 1000	20-2131-41	4 x 1 µmol	31.00	29.80	26.70	-
dT-CPG 1000	20-2131-42	4 x 0.2 µmol	31.00	29.80	26.70	-
dT-CPG 1000	20-2131-45	4 x 40 nmol	31.00	29.80	26.70	-
dT-CPG 1000	20-2131-62	200 x 200 nm	830.00	797.00	714.00	-
dT-CPG 1000	20-2231-14	1 x 15 µmol	133.00	128.00	114.00	-
dT-CPG 1000	20-2231-41	4 x 1 µmol	22.10	21.20	19.00	-
dT-CPG 1000	20-2231-42	4 x 0.2 µmol	15.50	14.90	13.30	-
dT-CPG 1000	20-2231-45	4 x 40 nmol	15.50	14.90	13.30	-
dT-CPG 1000 (1 g)	20-2031-10D	1 g	66.40	63.70	57.10	-
dT-CPG 1000, Expedite	20-2231-42E	4 x 0.2 µmol	15.50	14.90	13.30	-
dT-CPG 2000	20-2032-01	0.1 g	16.60	15.90	14.30	-
dT-CPG 2000	20-2032-02	0.25 g	33.20	31.90	28.60	-
dT-CPG 2000	20-2032-10	1 g	166.00	159.00	143.00	-
dT-CPG 2000	20-2132-42	4 x 0.2 µmol	31.00	29.80	26.70	-
dT-CPG 2000	20-2232-42	4 x 0.2 µmol	15.50	14.90	13.30	-

Description	Reference	Scale/Quantity	Price List 2025			
			EUR	CHF	GBP	USD
dT-CPG 2000, ABI 3900	20-2132-42A	4 x 0.2 µmol	31.00	29.80	26.70	-
dT-CPG 500	20-2030-01	0.1 g	13.30	12.80	11.40	-
dT-CPG 500	20-2030-02	0.25 g	17.70	17.00	15.20	-
dT-CPG 500	20-2030-10	1 g	66.40	63.70	57.10	-
dT-CPG 500	20-2130-13	1 x 10 µmol	88.50	85.00	76.10	-
dT-CPG 500	20-2130-41	4 x 1 µmol	31.00	29.80	26.70	-
dT-CPG 500	20-2130-42	4 x 0.2 µmol	31.00	29.80	26.70	-
dT-CPG 500	20-2230-14	1 x 15 µmol	133.00	128.00	114.00	-
dT-CPG 500	20-2230-41	4 x 1 µmol	22.10	21.20	19.00	-
dT-CPG 500	20-2230-42	4 x 0.2 µmol	15.50	14.90	13.30	-
dT-High Load-CPG	25-2030-02	0.25 g	22.10	21.20	19.00	-
dT-High Load-CPG	25-2030-10	1 g	79.70	76.50	68.50	-
dT-High Load-CPG	25-2130-17	1 x 25 µmol	111.00	107.00	95.50	-
dT-High Load-CPG	25-2130-46	4 x 2.5 µmol	66.40	63.70	57.10	-
dT-High Load-CPG	25-2230-18	1 x 35 µmol	164.00	157.00	141.00	-
dT-High Load-CPG	25-2230-46	4 x 2.5 µmol	66.40	63.70	57.10	-
dT-lcaa-CPG 500 Å - col.(96pcs)	20-2230-42SP	0.2 µmol	15.50	14.90	13.30	-
dT-PS	26-2630-62	200 x 200 nmol	653.00	627.00	562.00	-
dT-PS	26-2630-65	200 x 40 nmol	653.00	627.00	562.00	-
dT-Q-CPG 500	21-2230-42	4 x 0.2 µmol	44.30	42.50	38.10	-
dT-Thiophosphoramidite	10-1730-90	100 µmol	252.00	242.00	217.00	-
dT-Thiophosphoramidite,	10-1730-02	0.25 g	604.00	580.00	519.00	-
Empty MerMade Columns - 50 nmol	20-0050-05	Pk / 48	177.00	170.00	152.00	-
Empty Synthesis Columns - TWIST 10um/15um	20-0040-00	Pk / 10	265.00	254.00	228.00	-
Empty Synthesis Columns, 1µmol Expedite Style	20-0021-01	Pk / 10	42.50	40.80	36.60	-
Empty Synthesis Columns, 40 nmol, 0.2 µmol Expedite Style	20-0021-02	Pk / 10	42.50	40.80	36.60	-
Empty Synthesis Columns-TWIST 40nm, 0.2um or 1um	20-0030-00	Pk / 10	53.10	51.00	45.70	-
GAA Trimer Phosphoramidite	13-1200-90	100 µmol	1,294.00	1,242.00	1,113.00	-
GAA Trimer Phosphoramidite	13-1200-90M	100 µmol	1,294.00	1,242.00	1,113.00	-
GAA Trimer Phosphoramidite	13-1200-95	50 µmol	669.00	642.00	575.00	-
GAA Trimer Phosphoramidite	13-1200-95M	50 µmol	669.00	642.00	575.00	-
GAC Trimer Phosphoramidite	13-1201-90	100 µmol	1,294.00	1,242.00	1,113.00	-
GAC Trimer Phosphoramidite	13-1201-90M	100 µmol	1,294.00	1,242.00	1,113.00	-
GAC Trimer Phosphoramidite	13-1201-95	50 µmol	669.00	642.00	575.00	-
GAC Trimer Phosphoramidite	13-1201-95M	50 µmol	669.00	642.00	575.00	-
GAT Trimer Phosphoramidite	13-1203-90M	100 µmol	1,294.00	1,242.00	1,113.00	-
GAT Trimer Phosphoramidite	13-1203-95M	50 µmol	669.00	642.00	575.00	-
GCA Trimer Phosphoramidite	13-1210-90M	100 µmol	1,294.00	1,242.00	1,113.00	-
GCA Trimer Phosphoramidite	13-1210-95M	50 µmol	669.00	642.00	575.00	-
GCA Trimer Phosphoramidite, Expedite	13-1210-95E	50 µmol	669.00	642.00	575.00	-
GCG Trimer Phosphoramidite	13-1212-90M	100 µmol	1,294.00	1,242.00	1,113.00	-
GCG Trimer Phosphoramidite	13-1212-95M	50 µmol	669.00	642.00	575.00	-
GCT Trimer Phosphoramidite	13-1213-90	100 µmol	1,294.00	1,242.00	1,113.00	-
GCT Trimer Phosphoramidite	13-1213-90M	100 µmol	1,294.00	1,242.00	1,113.00	-
GCT Trimer Phosphoramidite	13-1213-95	50 µmol	669.00	642.00	575.00	-
GCT Trimer Phosphoramidite	13-1213-95M	50 µmol	669.00	642.00	575.00	-
GGT Trimer Phosphoramidite	13-1223-90	100 µmol	1,294.00	1,242.00	1,113.00	-
GGT Trimer Phosphoramidite	13-1223-90M	100 µmol	1,294.00	1,242.00	1,113.00	-
GGT Trimer Phosphoramidite	13-1223-95	50 µmol	669.00	642.00	575.00	-
GGT Trimer Phosphoramidite	13-1223-95M	50 µmol	669.00	642.00	575.00	-
Glen Gel-Pak™ 2.5 Desalting Column	61-5025-05	Pk / 5	93.50	89.80	80.40	-
Glen Gel-Pak™ 2.5 Desalting Column	61-5025-25	Pk / 25	460.00	442.00	396.00	-
Glen UnySupport PS	26-5140-52	10 x 200 nmol	97.40	93.50	83.80	-
Glen UnySupport PS	26-5140-55	10 x 40 nmol	97.40	93.50	83.80	-
Glen UnySupport™ 1000	20-5041-01	0.1 g	16.60	15.90	14.30	-
Glen UnySupport™ 1000	20-5041-02	0.25 g	27.70	26.60	23.80	-
Glen UnySupport™ 1000	20-5041-10	1 g	105.00	101.00	90.30	-
Glen UnySupport™ 1000	20-5141-13	1 x 10 µmol	88.50	85.00	76.10	-
Glen UnySupport™ 1000	20-5141-41	4 x 1 µmol	53.10	51.00	45.70	-
Glen UnySupport™ 1000	20-5141-42	4 x 0.2 µmol	35.40	34.00	30.40	-
Glen UnySupport™ 1000	20-5141-95	96 x 40 nmol	387.00	372.00	333.00	-
Glen UnySupport™ 1000	20-5241-14	1 x 15 µmol	123.00	123.00	111.00	-
Glen UnySupport™ 1000	20-5241-42	4 x 0.2 µmol	35.40	34.00	30.40	-
Glen UnySupport™ 1000 - Mermade	20-5141-91	96 x 1.0µm	581.00	558.00	500.00	-
Glen UnySupport™ 2000	20-5042-10	1.0 g	105.00	101.00	90.30	-
Glen UnySupport™ 500	20-5040-01	0.1 g	16.60	15.90	14.30	-
Glen UnySupport™ 500	20-5040-02	0.25 g	27.70	26.60	23.80	-
Glen UnySupport™ 500	20-5040-10	1 g	105.00	101.00	90.30	-
GTA Trimer Phosphoramidite	13-1230-90M	100 µmol	1,294.00	1,242.00	1,113.00	-
GTA Trimer Phosphoramidite	13-1230-95M	50 µmol	669.00	642.00	575.00	-
GTT Trimer Phosphoramidite	13-1233-90	100 µmol	1,294.00	1,242.00	1,113.00	-
GTT Trimer Phosphoramidite	13-1233-90M	100 µmol	1,294.00	1,242.00	1,113.00	-
GTT Trimer Phosphoramidite	13-1233-95	50 µmol	669.00	642.00	575.00	-
GTT Trimer Phosphoramidite	13-1233-95M	50 µmol	669.00	642.00	575.00	-
I-CE Phosphoramidite	10-3040-90	100 µmol	360.00	346.00	310.00	-
I-CE Phosphoramidite	10-3040-95	50 µmol	194.00	186.00	167.00	-
N4-Ac-N4-Et-dC-CE Phosphoramidite	10-1513-90	100 µmol	338.00	324.00	291.00	-
pdU-CE Phosphoramidite	10-1054-90	100 µmol	101.00	97.00	86.90	-
TAC Trimer Phosphoramidite	13-1301-90	100 µmol	1,294.00	1,242.00	1,113.00	-
TAC Trimer Phosphoramidite	13-1301-90M	100 µmol	1,294.00	1,242.00	1,113.00	-
TAC Trimer Phosphoramidite	13-1301-95	50 µmol	669.00	642.00	575.00	-
TAC Trimer Phosphoramidite	13-1301-95M	50 µmol	669.00	642.00	575.00	-
TCT Trimer Phosphoramidite	13-1313-90	100 µmol	1,294.00	1,242.00	1,113.00	-
TCT Trimer Phosphoramidite	13-1313-90M	100 µmol	1,294.00	1,242.00	1,113.00	-

Description	Reference	Scale/Quantity	Price List 2025				
			EUR	CHF	GBP	USD	
TCT Trimer Phosphoramidite	13-1313-95	50 µmol	669.00	642.00	575.00	-	
TCT Trimer Phosphoramidite	13-1313-95M	50 µmol	669.00	642.00	575.00	-	
TCT Trimer Phosphoramidite, Expedite	13-1313-90E	100 µmol	1,294.00	1,242.00	1,113.00	-	
TCT Trimer Phosphoramidite, Expedite	13-1313-95E	50 µmol	669.00	642.00	575.00	-	
TGC Trimer Phosphoramidite	13-1321-90	100 µmol	1,294.00	1,242.00	1,113.00	-	
TGC Trimer Phosphoramidite	13-1321-90M	100 µmol	1,294.00	1,242.00	1,113.00	-	
TGC Trimer Phosphoramidite	13-1321-95	50 µmol	669.00	642.00	575.00	-	
TGC Trimer Phosphoramidite	13-1321-95M	50 µmol	669.00	642.00	575.00	-	
TGG Trimer Phosphoramidite	13-1322-90	100 µmol	1,294.00	1,242.00	1,113.00	-	
TGG Trimer Phosphoramidite	13-1322-90M	100 µmol	1,294.00	1,242.00	1,113.00	-	
TGG Trimer Phosphoramidite	13-1322-95	50 µmol	669.00	642.00	575.00	-	
TGG Trimer Phosphoramidite	13-1322-95M	50 µmol	669.00	642.00	575.00	-	
Trimer Phosphoramidite Mix 1	13-1991-90	100 µmol	1,869.00	1,794.00	1,607.00	-	
Trimer Phosphoramidite Mix 1	13-1991-95	50 µmol	956.00	918.00	822.00	-	
TTC Trimer Phosphoramidite	13-1331-90	100 µmol	1,294.00	1,242.00	1,113.00	-	
TTC Trimer Phosphoramidite	13-1331-90M	100 µmol	1,294.00	1,242.00	1,113.00	-	
TTC Trimer Phosphoramidite	13-1331-95	50 µmol	669.00	642.00	575.00	-	
TTC Trimer Phosphoramidite	13-1331-95M	50 µmol	669.00	642.00	575.00	-	
TTT Trimer Phosphoramidite	13-1333-90M	100 µmol	1,294.00	1,242.00	1,113.00	-	
TTT Trimer Phosphoramidite	13-1333-95M	50 µmol	669.00	642.00	575.00	-	
Universal Support III	26-5010-10	1 g	252.00	242.00	217.00	-	
Glen Solvant							EA1B1Z
0.02M Iodine in Tetrahydrof./Pyrid./Wat.	40-4032-57		55.80	53.60	48.00	-	
0.02M Iodine in Tetrahydrof./Pyrid./Wat.	40-4032-62		252.00	242.00	217.00	-	
0.02M Iodine in Tetrahydrofuran/ Water/ Pyridine	40-4132-52	200 mL	34.90	33.50	30.00	-	
0.02M Iodine in Tetrahydrofuran/ Water/ Pyridine	40-4132-57	450 mL	74.30	71.30	63.90	-	
0.02M Iodine in Tetrahydrofuran/ Water/ Pyridine	40-4132-62	2000 mL	252.00	242.00	217.00	-	
0.02M Iodine in Tetrahydrofuran/Pyridine/Water	40-4330-52	200 mL	31.00	29.80	26.70	-	
0.02M Iodine in Tetrahydrofuran/Pyridine/Water	40-4330-57	450 mL	55.80	53.60	48.00	-	
0.02M Iodine in Tetrahydrofuran/Pyridine/Water	40-4330-61	960 mL	119.00	114.00	102.00	-	
0.02M Iodine in Tetrahydrofuran/Pyridine/Water	40-4330-62	2000 mL	252.00	242.00	217.00	-	
0.02M Iodine in Tetrahydrofuran/Pyridine/Water, Mermade	40-4330-57M	450 ml	55.80	53.60	48.00	-	
0.05 M Potassium Carbonate in Methanol	60-4600-30	30 mL	23.20	22.30	20.00	-	
0.05 M Sulfurizing Reagent II in pyridine/acetonitrile, Expedite	40-4137-52E	200 ml	155.00	149.00	133.00	-	
0.05M I2 in Pyridine/H2O	40-4035-71	1000 mL	182.00	175.00	157.00	-	
0.25 M 5-Benzylthio-1H-Tetrazole in Acetonitrile	30-3172-52	200 mL	85.20	81.80	73.30	-	
0.25 M 5-Benzylthio-1H-tetrazole in Acetonitrile	30-3170-45	45 mL	34.90	33.50	30.00	-	
0.25 M 5-Benzylthio-1H-tetrazole in Acetonitrile	30-3170-52	200 mL	85.20	81.80	73.30	-	
0.25 M 5-Benzylthio-1H-tetrazole in Acetonitrile	30-3170-57	450 mL	166.00	159.00	143.00	-	
0.25 M 5-Benzylthio-1H-tetrazole in Acetonitrile	30-3170-62	2000 mL	600.00	576.00	516.00	-	
0.25 M 5-Ethylthio-1H-Tetrazole in Anhydrous Acetonitrile	30-3140-45	45 mL	34.90	33.50	30.00	-	
0.25 M 5-Ethylthio-1H-Tetrazole in Anhydrous Acetonitrile	30-3140-52	200 mL	85.20	81.80	73.30	-	
0.25 M 5-Ethylthio-1H-Tetrazole in Anhydrous Acetonitrile	30-3140-57	450 mL	166.00	159.00	143.00	-	
0.25 M 5-Ethylthio-1H-Tetrazole in Anhydrous Acetonitrile	30-3140-61	960 mL	290.00	278.00	249.00	-	
0.25 M 5-Ethylthio-1H-Tetrazole in Anhydrous Acetonitrile	30-3140-62	2000 mL	600.00	576.00	516.00	-	
0.25 M 5-Ethylthio-1H-Tetrazole in Anhydrous Acetonitrile	30-3142-52	200 mL	85.20	81.80	73.30	-	
0.25 M DCI in Anhydrous Acetonitrile	30-3150-45	45 mL	34.90	33.50	30.00	-	
0.25 M DCI in Anhydrous Acetonitrile	30-3150-52	200 mL	85.20	81.80	73.30	-	
0.25 M DCI in Anhydrous Acetonitrile	30-3150-57	450 mL	166.00	159.00	143.00	-	
0.25 M DCI in Anhydrous Acetonitrile	30-3150-62	2000 mL	600.00	576.00	516.00	-	
0.25 M DCI in Anhydrous Acetonitrile	30-3152-52	200 mL	85.20	81.80	73.30	-	
0.25M 5-Benzylthio-1H-Tetrazole in Acetonitrile, Expedite	30-3172-52E	200 ml	85.20	81.80	73.30	-	
0.5 M CSO in Anhydr Acetonitrile 450 ml	40-4632-57	450 ml	441.00	423.00	379.00	-	
0.5 M CSO in Anhydrous Acetonitrile	40-4632-62	2 L	1,955.00	1,877.00	1,681.00	-	
0.5 M CSO in Anhydrous Acetonitrile (Expedite)	40-4632-52E	200 mL	201.00	193.00	173.00	-	
0.5 M CSO in Anhydrous Acetonitrile 1 l	40-4632-71	1 l	968.00	929.00	832.00	-	
0.5M CSO in Anhydrous Acetonitrile	40-4632-52	200 mL	201.00	193.00	173.00	-	
10 % 1-Methylimidazole in Tetrahydrofuran/Pyridine, Expedite	40-4122-52E	200 ml	34.90	33.50	30.00	-	
10% 1-Methylimidazole in Tetrahydrofuran	40-4120-52	200 mL	31.00	29.80	26.70	-	
10% 1-Methylimidazole in Tetrahydrofuran	40-4120-57	450 mL	69.70	66.90	59.90	-	
10% 1-Methylimidazole in Tetrahydrofuran	40-4120-62	2000 mL	252.00	242.00	217.00	-	
10% 1-Methylimidazole in Tetrahydrofuran, Mermade	40-4120-57M	450 ml	69.70	66.90	59.90	-	
10% 1-Methylimidazole in Tetrahydrofuran/Pyridine	40-4122-52	200 mL	34.90	33.50	30.00	-	
10% 1-Methylimidazole in Tetrahydrofuran/Pyridine	40-4122-57	450 mL	74.30	71.30	63.90	-	
10% 1-Methylimidazole in Tetrahydrofuran/Pyridine	40-4122-62	2000 mL	252.00	242.00	217.00	-	
16% 1-Methylimidazole in Tetrahydrofuran	40-4220-52	200 mL	34.90	33.50	30.00	-	
16% 1-Methylimidazole in Tetrahydrofuran	40-4220-57	450 mL	74.30	71.30	63.90	-	
16% 1-Methylimidazole in Tetrahydrofuran	40-4220-61	960 mL	158.00	152.00	136.00	-	
16% 1-Methylimidazole in Tetrahydrofuran	40-4220-62	2000 mL	329.00	316.00	283.00	-	
2% Aqueous Trifluoroacetic Acid	60-4040-57	450 mL	44.30	42.50	38.10	-	
2.0M Triethylamine Acetate, HPLC grade, pH=7	60-4110-52	200 mL	53.10	51.00	45.70	-	
2.0M Triethylamine Acetate, HPLC grade, pH=7	60-4110-57	450 mL	106.00	102.00	91.20	-	
2.0M Triethylamine Acetate, HPLC grade, pH=7	60-4110-60	960 mL	177.00	170.00	152.00	-	
2.0M Triethylamine Acetate, HPLC grade, pH=7	60-4110-62	2000 mL	354.00	340.00	304.00	-	
2.0M Triethylamine Acetate, HPLC grade, pH=7, Mermade	60-4110-62M	2000 ml	354.00	340.00	304.00	-	
2.5% Dichloroacetic acid in Dichloromethane	40-4042-57	450 mL	54.20	52.00	46.60	-	
2.5% Dichloroacetic acid in Dichloromethane	40-4042-62	2000 mL	159.00	153.00	137.00	-	
20 % 1-Methylimidazole in Acetonitrile / 2,6-Lutidine	40-4115-71	1000 ml	271.00	260.00	233.00	-	
20% Acetic Anhydride in Acetonitrile 1l	40-4224-71	1 l	112.00	108.00	96.30	-	
3 % TCA/DCM, Expedite	40-4140-61E	960 mL	73.60	70.70	63.30	-	
3% DCA in Toluene	40-4240-71	1000 mL	155.00	149.00	133.00	-	
3% Dichloroacetic acid in Dichloromethane	40-4040-57	450 mL	54.20	52.00	46.60	-	
3% Dichloroacetic acid in Dichloromethane	40-4040-61	960 mL	89.10	85.50	76.60	-	
3% Dichloroacetic acid in Dichloromethane	40-4040-62	2000 mL	174.00	167.00	150.00	-	

Description	Reference	Scale/Quantity	Price List 2025				
			EUR	CHF	GBP	USD	
3% Dichloroacetic acid in Dichloromethane	40-4040-71	1000 mL	96.80	92.90	83.20	-	
3% Dichloroacetic acid in Dichloromethane, Mermade	40-4040-62M	2000 ml	174.00	167.00	150.00	-	
3% TCA/DCM	40-4140-57	450 mL	38.70	37.20	33.30	-	
3% TCA/DCM	40-4140-61	960 mL	73.60	70.70	63.30	-	
3% TCA/DCM	40-4140-62	2000 mL	136.00	131.00	117.00	-	
3% TCA/DCM	40-4140-71	1000 mL	77.40	74.30	66.60	-	
4,5-Dicyanoimidazole (DCI), 5 g	30-3060-50	5 g	54.20	52.00	46.60	-	
4,5-Dicyanoimidazole (DCI), Crystalline	30-3050-10	1 g	27.10	26.00	23.30	-	
4,5-Dicyanoimidazole (DCI), Crystalline	30-3050-25	25 g	387.00	372.00	333.00	-	
5% Phenoxyacetic anhydride in THF	40-4212-52	200 mL	155.00	149.00	133.00	-	
5% Phenoxyacetic anhydride in THF	40-4212-57	450 mL	348.00	334.00	299.00	-	
5% Phenoxyacetic anhydride in THF, Expedite	40-4212-52E	200 mL	155.00	149.00	133.00	-	
5% Phenoxyacetic anhydride in THF/Pyridine	40-4210-52	200 mL	163.00	156.00	140.00	-	
5% Phenoxyacetic anhydride in THF/Pyridine	40-4210-57	450 mL	356.00	342.00	306.00	-	
5-Benzylthio-1H-tetrazole (BTT)	30-3070-10	1 g	27.10	26.00	23.30	-	
5-Benzylthio-1H-tetrazole (BTT)	30-3070-20	2 g	46.50	44.60	40.00	-	
5-Benzylthio-1H-tetrazole (BTT)	30-3070-25	25 g	387.00	372.00	333.00	-	
5-Benzylthio-1H-tetrazole (BTT), Mermade	30-3070-25M	25 g	387.00	372.00	333.00	-	
6.5% Dimethylaminopyridine in Tetrahydrofuran	40-4020-52	200 mL	50.30	48.30	43.30	-	
Acetonitrile, anhydrous	40-4050-45	60 mL	15.50	14.90	13.30	-	
Acetonitrile, anhydrous	40-4050-50	100 mL	19.40	18.60	16.70	-	
Acetonitrile, anhydrous	40-4050-53	300 mL	34.90	33.50	30.00	-	
Acetonitrile, anhydrous	40-4050-57	450 mL	38.70	37.20	33.30	-	
Acetonitrile, anhydrous, Mermade	40-4050-57M	450 mL	38.70	37.20	33.30	-	
Acetonitrile/Melm	40-4015-71	1000 mL	163.00	156.00	140.00	-	
Cap Mix (40-4026-75+40-4027-75ref)	40-4028-71	500 mL	201.00	193.00	173.00	-	
Sublimed 1H-Tetrazole in Anhydrous Acetonitrile	30-3100-45	45 mL	34.90	33.50	30.00	-	
Sublimed 1H-Tetrazole in Anhydrous Acetonitrile	30-3100-52	200 mL	85.20	81.80	73.30	-	
Sublimed 1H-Tetrazole in Anhydrous Acetonitrile	30-3100-57	450 mL	166.00	159.00	143.00	-	
Sublimed 1H-Tetrazole in Anhydrous Acetonitrile	30-3100-62	2000 mL	600.00	576.00	516.00	-	
Sublimed 1H-Tetrazole in Anhydrous Acetonitrile	30-3102-52	200 mL	85.20	81.80	73.30	-	
Sublimed 1H-Tetrazole in Anhydrous Acetonitrile, Expedite	30-3102-52E	200 mL	85.20	81.80	73.30	-	
Sulfurizing Reagent II	40-4037-10	1 g	38.70	37.20	33.30	-	
Sulfurizing Reagent II	40-4037-20	2 g	77.40	74.30	66.60	-	
Sulfurizing Reagent II, Mermade	40-4037-20M	2 g	77.40	74.30	66.60	-	
Sulfurizing Solution	40-4137-51	150 mL	77.40	74.30	66.60	-	
Sulfurizing Solution	40-4137-52	200 mL	155.00	149.00	133.00	-	
Sulfurizing Solution	40-4137-57	450 mL	348.00	334.00	299.00	-	
Tetrahydrofuran/ 2,6-Lutidine/ Acetic Anhydride	40-4010-52	200 mL	31.00	29.80	26.70	-	
Tetrahydrofuran/ 2,6-Lutidine/ Acetic Anhydride	40-4010-57	450 mL	69.70	66.90	59.90	-	
Tetrahydrofuran/ 2,6-Lutidine/ Acetic Anhydride	40-4010-61	960 mL	119.00	114.00	102.00	-	
Tetrahydrofuran/ 2,6-Lutidine/ Acetic Anhydride	40-4010-62	2000 mL	252.00	242.00	217.00	-	
Tetrahydrofuran/ Acetic Anhydride	40-4012-52	200 mL	31.00	29.80	26.70	-	
Tetrahydrofuran/ Acetic Anhydride	40-4012-57	450 mL	55.80	53.60	48.00	-	
Tetrahydrofuran/ Acetic Anhydride	40-4012-62	2000 mL	213.00	204.00	183.00	-	
Tetrahydrofuran/ Pyridine/ Acetic Anhydride	40-4110-52	200 mL	31.00	29.80	26.70	-	
Tetrahydrofuran/ Pyridine/ Acetic Anhydride	40-4110-57	450 mL	55.80	53.60	48.00	-	
Tetrahydrofuran/ Pyridine/ Acetic Anhydride	40-4110-62	2000 mL	252.00	242.00	217.00	-	
Tetrahydrofuran/ Pyridine/ Acetic Anhydride, Mermade	40-4110-57M	450 ml	55.80	53.60	48.00	-	
Glen Purifications							EA1C1Z
2.0M Hexylammonium Acetate, HPLC grade, pH=7 - 200 ml	60-4210-52	200 ml	66.40	63.70	57.10	-	
Adapter Rack for 96 well manifolds, Pk/1	60-0010-01		22.10	21.20	19.00	-	
Glen Gel-Pak™ 1.0 Desalting Column - Pk / 50	61-5010-50	Pk / 50	690.00	662.00	593.00	-	
Glen UnySupport™ 1000	20-5241-41	4 x 1 µmol	53.10	51.00	45.70	-	
Glen-Pak 50 mg DNA Purification Cartridge	60-5000-96	1 pack of 96 pce	459.00	441.00	395.00	-	
Glen-Pak™ DNA 3mg 384-Well Plate	60-5500-01	Pk / 1	597.00	573.00	513.00	-	
Glen-Pak™ DNA 30 mg 96-Well Plate	60-5400-01	Pk / 1	525.00	504.00	452.00	-	
Glen-Pak™ DNA Cartridge 3 g	60-5300-01	Pk / 1	133.00	128.00	114.00	-	
Glen-Pak™ DNA purification cartridge	60-5100-30	Pk / 30	177.00	170.00	152.00	-	
Glen-Pak™ DNA purification cartridge (for use with disposable syringes)	60-5200-01	Pk / 1	22.10	21.20	19.00	-	
Seal for unused wells on Adapter Rack, Pk/1	60-0020-01		31.00	29.80	26.70	-	
Glen Research reagents for cartridge purification							EA1D1Z
1-Ethynyl-dSpacer CE Phosphoramidite	10-1910-90	100 µmol	561.00	539.00	482.00	-	
1-Me-A-CE Phosphoramidite	10-3501-90	100 µmol	575.00	552.00	495.00	-	
1-Me-A-CE Phosphoramidite, Exp	10-3501-90E	100 µmol	575.00	552.00	495.00	-	
1-Me-dA-CE Phosphoramidite	10-1501-02	0.25 g	1,079.00	1,036.00	928.00	-	
1-Me-dA-CE Phosphoramidite	10-1501-90	100 µmol	360.00	346.00	310.00	-	
1-Me-dA-CE Phosphoramidite	10-1501-95	50 µmol	180.00	173.00	155.00	-	
1-Methyl-PseudoUridine Phosphoramidite	10-3056-02	0.25 g	4,034.00	3,873.00	3,469.00	-	
1-Methyl-PseudoUridine Phosphoramidite	10-3056-95	50 µmol	726.00	697.00	624.00	-	
1-Methyl-PseudoUridine Phosphoramidite - 100 µmol	10-3056-90	100 µmol	1,424.00	1,367.00	1,225.00	-	
1-MT-LA-CE Phosphoramidite, Expedite	10-2030-05E	0.5 g	108.00	104.00	92.90	-	
2'-F-5-Me-U-ANA-CE Phosphoramidite	10-3850-02	0.25 g	439.00	421.00	378.00	-	
2'-F-5-Me-U-ANA-CE Phosphoramidite	10-3850-05	0.5 g	863.00	828.00	742.00	-	
2'-F-G-CE Phosphoramidite, Expedite	10-3420-02E	0.25 g	108.00	104.00	92.90	-	
2',3'-ddA-CE Phosphoramidite	10-7001-02	0.25 g	964.00	925.00	829.00	-	
2',3'-ddA-CE Phosphoramidite	10-7001-90	100 µmol	230.00	221.00	198.00	-	
2',3'-ddC-CE Phosphoramidite	10-7101-02	0.25 g	964.00	925.00	829.00	-	
2',3'-ddC-CE Phosphoramidite	10-7101-90	100 µmol	230.00	221.00	198.00	-	
2',3'-ddC-CE Phosphoramidite - Mermade	10-7101-02M	0.25 g	964.00	925.00	829.00	-	
2',3'-ddC-CPG	20-2017-01	0.1 g	332.00	319.00	286.00	-	
2',3'-ddC-CPG	20-2017-10	1 g	3,319.00	3,186.00	2,854.00	-	
2',3'-ddC-CPG	20-2117-41	4 x 1 µmol	664.00	637.00	571.00	-	
2',3'-ddC-CPG	20-2117-42	4 x 0.2 µmol	221.00	212.00	190.00	-	

Description	Reference	Scale/Quantity	Price List 2025			
			EUR	CHF	GBP	USD
2',3'-ddC-CPG, Mermade	20-2117-42M	4 x 0.2 µmol	221.00	212.00	190.00	-
2',3'-ddG-CE Phosphoramidite	10-7201-02	0.25 g	1,122.00	1,077.00	965.00	-
2',3'-ddG-CE Phosphoramidite	10-7201-90	100 µmol	230.00	221.00	198.00	-
2',3'-ddT-CE Phosphoramidite	10-7301-02	0.25 g	964.00	925.00	829.00	-
2',3'-ddT-CE Phosphoramidite	10-7301-90	100 µmol	230.00	221.00	198.00	-
2,6-Diaminopurine-TOM-CE Phosphoramidite	10-3085-90	100 µmol	611.00	587.00	525.00	-
2,6-Diaminopurine-TOM-CE Phosphoramidite	10-3085-95	50 µmol	306.00	294.00	263.00	-
2.0 M Hexylammonium Acetate, HPLC grade, pH = 7	60-4210-57	450 ml	133.00	128.00	114.00	-
2',3'-ddC-Icaa-CPG 500 Å, Beckman - 0.2 µmol	20-2117-42B	0.2 µmol	221.00	212.00	190.00	-
2',3'-ddC-Icaa-CPG 500 Å - 4 x 1 µmol	20-2117-41E	4 x 1 µmol	664.00	637.00	571.00	-
2'-OMe-C-Thiophosphoramidite	10-3171-02	0.25 g	1,035.00	994.00	890.00	-
2-Amino-dA-CE Phosphoramidite	10-1085-02	0.25 g	417.00	400.00	359.00	-
2-Amino-dA-CE Phosphoramidite	10-1085-90	100 µmol	187.00	180.00	161.00	-
2-Amino-dA-CE Phosphoramidite	10-1085-95	50 µmol	108.00	104.00	92.90	-
2-Aminopurine-CE Phosphoramidite	10-1046-90E	100 µmol	209.00	201.00	180.00	-
2-Aminopurine-CE Phosphoramidite	10-1046-02	0.25 g	597.00	573.00	513.00	-
2-Aminopurine-CE Phosphoramidite	10-1046-90	100 µmol	209.00	201.00	180.00	-
2-Aminopurine-CE Phosphoramidite, Expedite - 0.25 g	10-1046-02E	0.25 g	597.00	573.00	513.00	-
2'-DeoxyNebularine-CE Phosphoramidite (Purine)	10-1041-02	0.25 g	453.00	435.00	390.00	-
2'-DeoxyNebularine-CE Phosphoramidite (Purine)	10-1041-90	100 µmol	165.00	158.00	142.00	-
2'-DeoxyNebularine-CE Phosphoramidite (Purine), Expedite	10-1041-02E	0.25 g	453.00	435.00	390.00	-
2'-deoxypseudoU-CE Phosphora.	10-1055-95P	50 µmol	288.00	276.00	248.00	-
2'-deoxypseudoU-CE Phosphoramidite	10-1055-02	0.25 g	1,848.00	1,774.00	1,589.00	-
2'-deoxypseudoU-CE Phosphoramidite	10-1055-90	100 µmol	546.00	524.00	470.00	-
2'-deoxypseudoU-CE Phosphoramidite	10-1055-95	50 µmol	288.00	276.00	248.00	-
2'-deoxypseudoU-CE Phosphoramidite, Expedite - 0.25 g	10-1055-02E	0.25 g	1,848.00	1,774.00	1,589.00	-
2'-F-A-ANA-CE Phosphoramidite	10-3800-02	0.25 g	769.00	738.00	661.00	-
2'-F-A-ANA-CE Phosphoramidite, Mermade	10-3800-02M	0.25 g	769.00	738.00	661.00	-
2'-F-Ac-C-ANA-CE Phosphoramidite	10-3815-02	0.25 g	762.00	732.00	655.00	-
2'-F-Ac-C-ANA-CE Phosphoramidite	10-3815-05	0.5 g	1,510.00	1,450.00	1,299.00	-
2'-F-Ac-C-ANA-CE Phosphoramidite, Mermade	10-3815-02M	0.25 g	762.00	732.00	655.00	-
2'-F-A-CE Phosphoramidite	10-3400-02	0.25 g	108.00	104.00	92.90	-
2'-F-A-CE Phosphoramidite	10-3400-05	0.5 g	216.00	207.00	186.00	-
2'-F-A-CE Phosphoramidite, Expedite	10-3400-05E	0.5 g	216.00	207.00	186.00	-
2'-F-dl-CE Phosphoramidite-50 Exp.	10-1085-95E	50 µmol	108.00	104.00	92.90	-
2'-F-dl-CE Phosphoramidite	10-1082-02	0.25 g	1,539.00	1,477.00	1,324.00	-
2'-F-dl-CE Phosphoramidite	10-1082-90	100 µmol	582.00	559.00	501.00	-
2'-F-dl-CE Phosphoramidite	10-1082-95	50 µmol	316.00	303.00	272.00	-
2'-F-dl-CE Phosphoramidite- Expedite	10-1082-95E	50 µmol	316.00	303.00	272.00	-
2'-F-G-ANA-CE Phosphoramidite	10-3820-02	0.25 g	748.00	718.00	643.00	-
2'-F-G-ANA-CE Phosphoramidite	10-3820-90	100 µmol	273.00	262.00	235.00	-
2'-F-G-CE Phosphoramidite	10-3420-02	0.25 g	108.00	104.00	92.90	-
2'-F-G-CE Phosphoramidite	10-3420-05	0.5 g	216.00	207.00	186.00	-
2'-F-G-CE Phosphoramidite, Expedite	10-3420-05E	0.5 g	216.00	207.00	186.00	-
2'-F-I-CE Phosphoramidite, Mermade	10-3440-02M	0.25 g	273.00	262.00	235.00	-
2'-F-U-CE Phosphoramidite, Expedite	10-3430-05E	144.00	138.00	124.00	-	
2'-OMe-Ac-C-CE Phosphoramidite, Expedite	10-3115-10E	1 g	122.00	117.00	105.00	-
2'-OMe-Ac-C-RNA-CPG, Expedite	20-3715-41E	4 x 1 µmol	144.00	138.00	124.00	-
2'-OMe-A-CE Phosphoramidite, Expedite	10-3100-10E	1 g	122.00	117.00	105.00	-
2'-OMe-A-RNA-CPG, Expedite	20-3700-41E	4 x 1 µmol	144.00	138.00	124.00	-
2'-OMe-C-RNA-CPG, Expedite	20-3710-41E	4 x 1 µmol	144.00	138.00	124.00	-
2'-OMe-G-CE Phosphoramidite, Expedite	10-3121-10E	1 g	199.00	191.00	171.00	-
2'-OMe-G-RNA-CPG, Expedite	20-3721-41E	4 x 1 µmol	144.00	138.00	124.00	-
2'-OMe-ibu-G-CE Phosphoramidite	10-3120-10	1.0 g	122.00	117.00	105.00	-
2'-OMe-ibu-G-CE Phosphoramidite, Expedite	10-3120-10E	1.0 g	122.00	117.00	105.00	-
2'-OMe-iPr-Pac-G-CE Phosphoramidite, Expedite	10-3621-10E	1 g	360.00	346.00	310.00	-
2'-OMe-Pac-A-CE Phosphoramidite	10-3601-10	1 g	360.00	346.00	310.00	-
2'-OMe-Pac-A-CE Phosphoramidite, Expedite	10-3601-10E	1 g	360.00	346.00	310.00	-
2'-OMe-U-CE Phosphoramidite, Expedite	10-3130-10E	1 g	122.00	117.00	105.00	-
2'-OMe-U-RNA-CPG	20-3630-01	0.1 g	57.50	55.20	49.50	-
2'-OMe-U-RNA-CPG, Expedite	20-3730-41E	4 x 1 µmol	144.00	138.00	124.00	-
2-Thio-dT-CE Phosphoramidite	10-1036-02	0.25 g	1,136.00	1,091.00	977.00	-
2-Thio-dT-CE Phosphoramidite	10-1036-90	100 µmol	446.00	428.00	384.00	-
2-Thio-dT-CE Phosphoramidite	10-1036-95	50 µmol	244.00	234.00	210.00	-
3' - (6 - FAM)CPG -0.2 µm(1pk/4)Mermade	20-2961-42M	4 x 0.2 µmol	144.00	138.00	124.00	-
3' - Amino-ModifierC7 CPG1000-(1pk/4)Mer	20-2958-41M	4 x 1 µmol	170.00	163.00	146.00	-
3' - CPR II CPG	20-2903-01	0.1 g	101.00	97.00	86.90	-
3' -Amino-ModifierC7 CPG1000-(1pk/4)Mer.	20-2958-42M	4 x 0.2 µmol	94.10	90.30	80.90	-
3' Thiol-Modifier 6 S-S CPG	20-2938-01	0.1 g	129.00	124.00	111.00	-
3' Thiol-Modifier 6 S-S CPG	20-2938-10	1 g	1,186.00	1,139.00	1,020.00	-
3'-(6-FAM) CPG	20-2961-01	0.1 g	180.00	173.00	155.00	-
3'-(6-FAM) CPG	20-2961-10	1 g	1,510.00	1,450.00	1,299.00	-
3'-(6-FAM) CPG	20-2961-13	1 x 10 µmol	475.00	456.00	409.00	-
3'-(6-FAM) CPG	20-2961-14	1 x 15 µmol	712.00	684.00	612.00	-
3'-(6-FAM) CPG	20-2961-41	4 x 1 µmol	256.00	246.00	220.00	-
3'-(6-FAM) CPG	20-2961-42	4 x 0.2 µmol	144.00	138.00	124.00	-
3'-(6-FAM) PS	26-2961-01	0.1 g	216.00	207.00	186.00	-
3'-(6-FAM) PS	26-2961-10	1 g	1,977.00	1,898.00	1,700.00	-
3'-(6-FAM) PS	26-2961-52	10 x 200 nmol	475.00	456.00	409.00	-
3'-(6-FAM) PS	26-2961-55	10 x 40 nmol	475.00	456.00	409.00	-
3'-(6-Fluorescein) CPG	20-2964-01	0.1 g	173.00	166.00	149.00	-
3'-(6-Fluorescein) CPG	20-2964-10	1 g	1,431.00	1,374.00	1,231.00	-
3'-(6-Fluorescein) CPG	20-2964-13	1 x 10 µmol	431.00	414.00	371.00	-
3'-(6-Fluorescein) CPG	20-2964-14	1 x 15 µmol	647.00	621.00	556.00	-

Description	Reference	Scale/Quantity	Price List 2025			
			EUR	CHF	GBP	USD
3 ⁺ -(6-Fluorescein) CPG	20-2964-41	4 x 1 µmol	288.00	276.00	248.00	-
3 ⁺ -(6-Fluorescein) CPG	20-2964-42	4 x 0.2 µmol	173.00	166.00	149.00	-
3 ⁺ -Acridine CPG	20-2973-01	0.1 g	173.00	166.00	149.00	-
3 ⁺ -Acridine CPG	20-2973-10	1 g	1,431.00	1,374.00	1,231.00	-
3 ⁺ -Acridine CPG	20-2973-13	1 x 10 µmol	431.00	414.00	371.00	-
3 ⁺ -Acridine CPG	20-2973-14	1 x 15 µmol	647.00	621.00	556.00	-
3 ⁺ -Acridine CPG	20-2973-41	4 x 1 µmol	288.00	276.00	248.00	-
3 ⁺ -Acridine CPG	20-2973-42	4 x 0.2 µmol	173.00	166.00	149.00	-
3 ⁺ -Alkyne-Modifier Serinol CPG	20-2992-01	0.1 g	151.00	145.00	130.00	-
3 ⁺ -Alkyne-Modifier Serinol CPG	20-2992-10	1 g	1,150.00	1,104.00	989.00	-
3 ⁺ -Alkyne-Modifier Serinol CPG	20-2992-41	4 x 1 µmol	252.00	242.00	217.00	-
3 ⁺ -Alkyne-Modifier Serinol CPG	20-2992-42	4 x 0.2 µmol	144.00	138.00	124.00	-
3 ⁺ -Alkyne-Modifier Serinol CPG	20-2992-42	4 x 0.2 µmol	144.00	138.00	124.00	-
3 ⁺ -Amino-dT CPG	20-2981-01	0.1 g	173.00	166.00	149.00	-
3 ⁺ -Amino-Modifier C6 dT CPG	20-2039-01	0.1 g	144.00	138.00	124.00	-
3 ⁺ -Amino-Modifier C6 dT CPG	20-2039-10	1 g	1,208.00	1,160.00	1,039.00	-
3 ⁺ -Amino-Modifier C6 dT CPG	20-2039-13	1 x 10 µmol	367.00	352.00	316.00	-
3 ⁺ -Amino-Modifier C6 dT CPG	20-2039-14	1 x 15 µmol	546.00	524.00	470.00	-
3 ⁺ -Amino-Modifier C6 dT CPG	20-2039-41	4 x 1 µmol	237.00	228.00	204.00	-
3 ⁺ -Amino-Modifier C6 dT CPG	20-2039-42	4 x 0.2 µmol	144.00	138.00	124.00	-
3 ⁺ -Amino-Modifier C6 dT CPG, Expedite	20-2039-41E	4 x 1 µmol	237.00	228.00	204.00	-
3 ⁺ -Amino-Modifier C7 CPG 1000	20-2958-01	0.1 g	137.00	132.00	118.00	-
3 ⁺ -Amino-Modifier C7 CPG 1000	20-2958-10	1 g	971.00	932.00	835.00	-
3 ⁺ -Amino-Modifier C7 CPG 1000	20-2958-13	1 x 10 µmol	360.00	346.00	310.00	-
3 ⁺ -Amino-Modifier C7 CPG 1000	20-2958-14	1 x 15 µmol	539.00	517.00	464.00	-
3 ⁺ -Amino-Modifier C7 CPG 1000	20-2958-41	4 x 1 µmol	170.00	163.00	146.00	-
3 ⁺ -Amino-Modifier C7 CPG 1000	20-2958-42	4 x 0.2 µmol	94.10	90.30	80.90	-
3 ⁺ -Amino-Modifier C7 CPG 1000, Expedite	20-2958-41E	4 x 1 µmol	170.00	163.00	146.00	-
3 ⁺ -Amino-Modifier Serinol CPG	20-2997-01	0.1 g	137.00	132.00	118.00	-
3 ⁺ -Amino-Modifier Serinol CPG	20-2997-10	1 g	971.00	932.00	835.00	-
3 ⁺ -Amino-Modifier Serinol CPG	20-2997-13	1 x 10 µmol	360.00	346.00	310.00	-
3 ⁺ -Amino-Modifier Serinol CPG	20-2997-14	1 x 15 µmol	539.00	517.00	464.00	-
3 ⁺ -Amino-Modifier Serinol CPG	20-2997-41	4 x 1 µmol	201.00	193.00	173.00	-
3 ⁺ -Amino-Modifier Serinol CPG	20-2997-41M	4 x 1.0 µmol	201.00	193.00	173.00	-
3 ⁺ -Amino-Modifier Serinol CPG	20-2997-42	4 x 0.2 µmol	122.00	117.00	105.00	-
3 ⁺ -Azido-Modifier Serinol CPG	20-2999-01	0.1 g	173.00	166.00	149.00	-
3 ⁺ -Azido-Modifier Serinol CPG	20-2999-10	1 g	1,431.00	1,374.00	1,231.00	-
3 ⁺ -Azido-Modifier Serinol CPG, Mermade	20-2999-10M	1 g	1,431.00	1,374.00	1,231.00	-
3 ⁺ -BBQ 650 CPG	20-5934-42	4 x 0.2 µmol	165.00	158.00	142.00	-
3 ⁺ -BHQ-1 CPG	20-5931-01	0.1 g	395.00	379.00	340.00	-
3 ⁺ -BHQ-1 CPG	20-5931-10	1 g	3,308.00	3,176.00	2,845.00	-
3 ⁺ -BHQ-1 CPG	20-5931-13	1 x 10 µmol	1,035.00	994.00	890.00	-
3 ⁺ -BHQ-1 CPG	20-5931-14	1 x 15 µmol	1,546.00	1,484.00	1,330.00	-
3 ⁺ -BHQ-1 CPG	20-5931-41	4 x 1 µmol	365.00	350.00	314.00	-
3 ⁺ -BHQ-1 CPG	20-5931-42	4 x 0.2 µmol	127.00	122.00	109.00	-
3 ⁺ -BHQ-1 CPG, Mermade	20-5931-42M	4 x 0.2 µmol	127.00	122.00	109.00	-
3 ⁺ -BHQ-1 CPG, ABI 3900	20-5931-42A	4 x 0.2 µmol	127.00	122.00	109.00	-
3 ⁺ -BHQ-1 CPG-0.2um(1pk/4) Expedite	20-5931-42E	4 x 0.2 µmol	127.00	122.00	109.00	-
3 ⁺ -BHQ-1 CPG-1.0um(1pk/4)Expedite	20-5931-41E	4 x 0.2 µmol	365.00	350.00	314.00	-
3 ⁺ -BHQ-1 CPG-1.0um(1pk/4)Mermade	20-5931-41M	4 x 0.2 µmol	365.00	350.00	314.00	-
3 ⁺ -BHQ-2 CPG	20-5932-01	0.1 g	395.00	379.00	340.00	-
3 ⁺ -BHQ-2 CPG	20-5932-10	1 g	3,308.00	3,176.00	2,845.00	-
3 ⁺ -BHQ-2 CPG	20-5932-13	1 x 10 µmol	1,035.00	994.00	890.00	-
3 ⁺ -BHQ-2 CPG	20-5932-14	1 x 15 µmol	1,546.00	1,484.00	1,330.00	-
3 ⁺ -BHQ-2 CPG	20-5932-41	4 x 1 µmol	365.00	350.00	314.00	-
3 ⁺ -BHQ-2 CPG	20-5932-42	4 x 0.2 µmol	115.00	110.00	98.90	-
3 ⁺ -BHQ-2 CPG, Mermade	20-5932-42M	4 x 0.2 µmol	115.00	110.00	98.90	-
3 ⁺ -BHQ-2 CPG-0.2um(1pk/4)Expedite	20-5932-42E	4 x 0.2 µmol	115.00	110.00	98.90	-
3 ⁺ -BHQ-2 CPG-1 µmol (1 pk/4) Mermade	20-5932-41M	4 x 0.2 µmol	365.00	350.00	314.00	-
3 ⁺ -BHQ-2 CPG-1.0um(1pk/4)Expedite	20-5932-41E	4 x 0.2 µmol	365.00	350.00	314.00	-
3 ⁺ -BHQ-3 CPG	20-5933-01	0.1 g	395.00	379.00	340.00	-
3 ⁺ -BHQ-3 CPG	20-5933-10	1 g	3,308.00	3,176.00	2,845.00	-
3 ⁺ -BHQ-3 CPG	20-5933-13	1 x 10 µmol	1,186.00	1,139.00	1,020.00	-
3 ⁺ -BHQ-3 CPG	20-5933-14	1 x 15 µmol	1,783.00	1,712.00	1,533.00	-
3 ⁺ -BHQ-3 CPG	20-5933-41	4 x 1 µmol	539.00	517.00	464.00	-
3 ⁺ -BHQ-3 CPG	20-5933-42	4 x 0.2 µmol	165.00	158.00	142.00	-
3 ⁺ -BHQ-3 CPG - ABI 3900	20-5933-42A	4 x 0.2 µmol	165.00	158.00	142.00	-
3 ⁺ -BHQ-3 CPG, Mermade	20-5933-42M	4 x 0.2 µmol	165.00	158.00	142.00	-
3 ⁺ -BHQ-3 CPG-4x0.2um Expedite	20-5933-42E	4 x 0.2 µmol	165.00	158.00	142.00	-
3 ⁺ -BHQ-3 CPG-4x1.0um	20-5933-41E	4 x 0.2 µmol	539.00	517.00	464.00	-
3 ⁺ -BiotinTEG CPG	20-2955-01	0.1 g	173.00	166.00	149.00	-
3 ⁺ -BiotinTEG CPG	20-2955-10	1 g	1,431.00	1,374.00	1,231.00	-
3 ⁺ -BiotinTEG CPG	20-2955-13	1 x 10 µmol	431.00	414.00	371.00	-
3 ⁺ -BiotinTEG CPG	20-2955-14	1 x 15 µmol	647.00	621.00	556.00	-
3 ⁺ -BiotinTEG CPG	20-2955-41	4 x 1 µmol	243.00	233.00	209.00	-
3 ⁺ -BiotinTEG CPG	20-2955-42	4 x 0.2 µmol	133.00	128.00	114.00	-
3 ⁺ -BiotinTEG CPG - ABI 3900	20-2955-42A	4 x 0.2 µmol	133.00	128.00	114.00	-
3 ⁺ -BiotinTEG PS	26-2955-52	10 x 200 nmol	431.00	414.00	371.00	-
3 ⁺ -BiotinTEG PS	26-2955-55	10 x 40 nmol	431.00	414.00	371.00	-
3 ⁺ -Cholesteryl-TEG CPG	20-2975-01	0.1 g	122.00	117.00	105.00	-
3 ⁺ -Cholesteryl-TEG CPG	20-2975-10	1 g	1,007.00	967.00	866.00	-
3 ⁺ -Cholesteryl-TEG CPG	20-2975-10	1 g	1,007.00	967.00	866.00	-
3 ⁺ -Cholesteryl-TEG CPG	20-2975-13	1 x 10 µmol	302.00	290.00	260.00	-
3 ⁺ -Cholesteryl-TEG CPG	20-2975-14	1 x 15 µmol	453.00	435.00	390.00	-

Description	Reference	Scale/Quantity	Price List 2025			
			EUR	CHF	GBP	USD
3 ⁻ -Cholesteryl-TEG CPG	20-2975-41	4 x 1 µmol	201.00	193.00	173.00	-
3 ⁻ -Cholesteryl-TEG CPG	20-2975-42	4 x 0.2 µmol	121.00	116.00	104.00	-
3 ⁻ -Cholesteryl-TEG CPG - Mermade	20-2975-41M	4 x 1 µmol	201.00	193.00	173.00	-
3 ⁻ -CPR II CPG	20-2903-10	1 g	690.00	662.00	593.00	-
3 ⁻ -CPR II CPG, Mermade	20-2903-41M	4 x 0.1 µmol	144.00	138.00	124.00	-
3 ⁻ -CPR II CPG, Mermade	20-2903-42M	4 x 0.2 µmol	86.30	82.80	74.20	-
3-Cyanovinylcarbazole Phosphoramidite (CNVK)	10-4960-02	0.25 g	1,618.00	1,553.00	1,391.00	-
3-Cyanovinylcarbazole Phosphoramidite (CNVK)	10-4960-90	100 µmol	561.00	539.00	482.00	-
3-Cyanovinylcarbazole Phosphoramidite (CNVK)	10-4960-95	50 µmol	288.00	276.00	248.00	-
3-Cyanovinylcarbazole Phosphoramidite (CNVK), Expedite	10-4960-90E	100 µmol	561.00	539.00	482.00	-
3 ⁻ -DabcyI CPG	20-5912-01	0.1 g	159.00	153.00	137.00	-
3 ⁻ -DabcyI CPG	20-5912-10	1 g	1,294.00	1,242.00	1,113.00	-
3 ⁻ -DabcyI CPG	20-5912-13	1 x 10 µmol	465.00	446.00	400.00	-
3 ⁻ -DabcyI CPG	20-5912-14	1 x 15 µmol	664.00	637.00	571.00	-
3 ⁻ -DabcyI CPG	20-5912-41	4 x 1 µmol	243.00	233.00	209.00	-
3 ⁻ -DabcyI CPG	20-5912-42	4 x 0.2 µmol	133.00	128.00	114.00	-
3 ⁻ -DabcyI PS	26-5912-52	10 x 200 nmol	475.00	456.00	409.00	-
3 ⁻ -DabcyI PS	26-5912-55	10 x 40 nmol	475.00	456.00	409.00	-
3 ⁻ -DabcyI-CPG-0.2µmol col.(1pk/4)Mermade	20-5912-42M	4 x 0.2 µmol	133.00	128.00	114.00	-
3 ⁻ -DabcyI-CPG-1000 col.	20-5912-41E	4 x 1 µmol	243.00	233.00	209.00	-
3 ⁻ -DabcyI-CPG1000 col1µM(1pk/4)Mermade	20-5912-41M	4 x 1 µmol	243.00	233.00	209.00	-
3 ⁻ -Dabsyl CPG	20-5911-01	0.1 g	187.00	180.00	161.00	-
3 ⁻ -Dabsyl CPG	20-5911-10	1 g	1,474.00	1,415.00	1,268.00	-
3 ⁻ -Dabsyl CPG	20-5911-13	1 x 10 µmol	554.00	532.00	476.00	-
3 ⁻ -Dabsyl CPG	20-5911-14	1 x 15 µmol	791.00	759.00	680.00	-
3 ⁻ -Dabsyl CPG	20-5911-41	4 x 1 µmol	268.00	257.00	230.00	-
3 ⁻ -Dabsyl CPG	20-5911-42	4 x 0.2 µmol	149.00	143.00	128.00	-
3 ⁻ -Dabsyl-CPG Col 0.2 µmol pk/4 Mermade	20-5911-42M	0.2 µmol	149.00	143.00	128.00	-
3 ⁻ -Dabsyl-CPG col.	20-5911-41E	4 x 1 µmol	268.00	257.00	230.00	-
3 ⁻ -Dabsyl-CPG Col. (Exp)	20-5911-13E	10 µmol	554.00	532.00	476.00	-
3 ⁻ -Dabsyl-CPG col. Expedite	20-5911-42E	0.2 µmol	149.00	143.00	128.00	-
3 ⁻ -Dabsyl-CPG-Col-1µM (4/pack)Mermade	20-5911-41M	4 x 1 µmol	268.00	257.00	230.00	-
3 ⁻ -dA-CE Phosphoramidite	10-1004-02	0.25 g	1,898.00	1,822.00	1,632.00	-
3 ⁻ -dA-CE Phosphoramidite	10-1004-90	100 µmol	683.00	656.00	587.00	-
3 ⁻ -dA-CE Phosphoramidite	10-1004-95	50 µmol	360.00	346.00	310.00	-
3 ⁻ -dA-CE Phosphoramidite - Expedite	10-1004-02E	0.25 g	1,898.00	1,822.00	1,632.00	-
3 ⁻ -dA-CPG	20-2004-01	0.1 g	813.00	780.00	699.00	-
3 ⁻ -dA-CPG	20-2104-41	4 x 1 µmol	1,014.00	973.00	872.00	-
3 ⁻ -dA-CPG	20-2104-42	4 x 0.2 µmol	360.00	346.00	310.00	-
3 ⁻ -dA-CPG 500 Å, column, Expedite - 0.2 µmol	20-2104-42E	0.2 µmol	360.00	346.00	310.00	-
3 ⁻ -dA-lcaa-CPG 500 Å - 1 µmol	20-2104-41E	1 µmol	1,014.00	973.00	872.00	-
3 ⁻ -DBCO-Serinol CPG - Expedite	20-2998-42E	4 x 0.2 µmol	129.00	124.00	111.00	-
3 ⁻ -dC-CE Phosphoramidite	10-1064-02	0.25 g	3,387.00	3,252.00	2,913.00	-
3 ⁻ -dC-CE Phosphoramidite	10-1064-90	100 µmol	1,136.00	1,091.00	977.00	-
3 ⁻ -dC-CE Phosphoramidite	10-1064-95	50 µmol	582.00	559.00	501.00	-
3 ⁻ -dC-CE Phosphoramidite - Expedite	10-1064-02E	0.25 g	3,387.00	3,252.00	2,913.00	-
3 ⁻ -dC-CPG	20-2064-01	0.1 g	431.00	414.00	371.00	-
3 ⁻ -dC-CPG	20-2164-41	4 x 1 µmol	863.00	828.00	742.00	-
3 ⁻ -dC-CPG	20-2164-42	4 x 0.2 µmol	288.00	276.00	248.00	-
3 ⁻ -dC-lcaa-CPG 500 Å	20-2164-41E	1 µmol	863.00	828.00	742.00	-
3 ⁻ -dC-lcaa-CPG 500 Å - 0.2 µmol	20-2164-42E	0.2 µmol	288.00	276.00	248.00	-
3 ⁻ -dG-CE Phosphoramidite	10-1074-02	0.25 g	2,020.00	1,939.00	1,737.00	-
3 ⁻ -dG-CE Phosphoramidite	10-1074-90	100 µmol	690.00	662.00	593.00	-
3 ⁻ -dG-CE Phosphoramidite	10-1074-95	50 µmol	360.00	346.00	310.00	-
3 ⁻ -dG-CE Phosphoramidite - Expedite	10-1074-02E	0.25 g	2,020.00	1,939.00	1,737.00	-
3 ⁻ -dG-CPG	20-2074-01	0.1 g	431.00	414.00	371.00	-
3 ⁻ -dG-CPG	20-2174-41	4 x 1 µmol	863.00	828.00	742.00	-
3 ⁻ -dG-CPG	20-2174-42	4 x 0.2 µmol	288.00	276.00	248.00	-
3 ⁻ -dG-lcaa-CPG 500 Å	20-2174-42E	0.2 µmol	288.00	276.00	248.00	-
3 ⁻ -dG-lcaa-CPG 500 Å - 1 µmol	20-2174-41E	1 µmol	863.00	828.00	742.00	-
3 ⁻ -Dithiol Serinol CPG	20-2991-01	0.1 g	173.00	166.00	149.00	-
3 ⁻ -dT-CE Phosphoramidite	10-1084-02	0.25 g	2,423.00	2,326.00	2,084.00	-
3 ⁻ -dT-CE Phosphoramidite	10-1084-90	100 µmol	762.00	732.00	655.00	-
3 ⁻ -dT-CE Phosphoramidite	10-1084-95	50 µmol	403.00	387.00	347.00	-
3 ⁻ -dT-CE Phosphoramidite - Expedite	10-1084-02E	0.25 g	2,423.00	2,326.00	2,084.00	-
3 ⁻ -dT-CPG	20-2084-01	0.1 g	431.00	414.00	371.00	-
3 ⁻ -dT-CPG	20-2184-41	4 x 1 µmol	863.00	828.00	742.00	-
3 ⁻ -dT-CPG	20-2184-42	4 x 0.2 µmol	288.00	276.00	248.00	-
3 ⁻ -dT-lcaa-CPG 500 Å	20-2184-42E	0.2 µmol	288.00	276.00	248.00	-
3 ⁻ -dT-lcaa-CPG 500 Å - 1 µmol	20-2184-41E	1 µmol	863.00	828.00	742.00	-
3 ⁻ -Fluorescein CPG	20-2963-01	0.1 g	173.00	166.00	149.00	-
3 ⁻ -Fluorescein CPG	20-2963-10	1 g	1,431.00	1,374.00	1,231.00	-
3 ⁻ -Fluorescein CPG	20-2963-13	1 x 10 µmol	431.00	414.00	371.00	-
3 ⁻ -Fluorescein CPG	20-2963-14	1 x 15 µmol	647.00	621.00	556.00	-
3 ⁻ -Fluorescein CPG	20-2963-41	4 x 1 µmol	288.00	276.00	248.00	-
3 ⁻ -Fluorescein CPG	20-2963-42	4 x 0.2 µmol	173.00	166.00	149.00	-
3 ⁻ -Fluorescein CPG, Expedite	20-2963-41E	4 x 1 µmol	288.00	276.00	248.00	-
3 ⁻ Fluorescein dT CPG, column - 1 µmol	20-2056-41E	1 µmol	302.00	290.00	260.00	-
3 ⁻ -Fluorescein-dT CPG	20-2056-01	0.1 g	180.00	173.00	155.00	-
3 ⁻ -Fluorescein-dT CPG	20-2056-10	1 g	1,503.00	1,443.00	1,293.00	-
3 ⁻ -Fluorescein-dT CPG	20-2056-13	1 x 10 µmol	453.00	435.00	390.00	-
3 ⁻ -Fluorescein-dT CPG	20-2056-14	1 x 15 µmol	676.00	649.00	581.00	-
3 ⁻ -Fluorescein-dT CPG	20-2056-41	4 x 1 µmol	302.00	290.00	260.00	-
3 ⁻ -Fluorescein-dT CPG	20-2056-42	4 x 0.2 µmol	180.00	173.00	155.00	-

Description	Reference	Scale/Quantity	Price List 2025			
			EUR	CHF	GBP	USD
3'-Glyceryl CPG	20-2902-01	0.1 g	122.00	117.00	105.00	-
3'-Glyceryl CPG	20-2902-10	1 g	863.00	828.00	742.00	-
3'-Glyceryl CPG	20-2902-13	1 x 10 µmol	324.00	311.00	279.00	-
3'-Glyceryl CPG	20-2902-14	1 x 15 µmol	503.00	483.00	433.00	-
3'-Glyceryl CPG	20-2902-41	4 x 1 µmol	180.00	173.00	155.00	-
3'-Glyceryl CPG	20-2902-42	4 x 0.2 µmol	108.00	104.00	92.90	-
3'-Glyceryl CPG, column, Pharmacia	20-2902-41P	1 µmol (4 x 1 pce)	180.00	173.00	155.00	-
3'-Phosphate CPG	20-2900-01	0.1 g	93.50	89.80	80.40	-
3'-Phosphate CPG	20-2900-10	1 g	662.00	636.00	569.00	-
3'-Phosphate CPG	20-2900-13	1 x 10 µmol	244.00	234.00	210.00	-
3'-Phosphate CPG	20-2900-14	1 x 15 µmol	388.00	372.00	334.00	-
3'-Phosphate CPG	20-2900-41	4 x 1 µmol	122.00	117.00	105.00	-
3'-Phosphate CPG	20-2900-41E	1 µmol	122.00	117.00	105.00	-
3'-Phosphate CPG	20-2900-42	4 x 0.2 µmol	66.40	63.70	57.10	-
3'-Phosphate CPG (High Load)	25-2900-01	0.1 g	122.00	117.00	105.00	-
3'-Phosphate CPG (High Load)	25-2900-10	1 g	863.00	828.00	742.00	-
3'-Phosphate CPG, column	20-2900-41P	4 x 1 µmol	122.00	117.00	105.00	-
3'-Phosphate CPG, Exp.	20-2900-42E	4 x 0.2 µmol	66.40	63.70	57.10	-
3'-Phosphate PS	26-2900-52	10 x 200 nmol	216.00	207.00	186.00	-
3'-Phosphate PS	26-2900-55	10 x 40 nmol	216.00	207.00	186.00	-
3'-PhosphateCPG-0.2µm.col(1pk/4)Mermade	20-2900-42M	0.2 µmol	66.40	63.70	57.10	-
3'-PhosphateCPG-1 µm.col.(1pk/4)Mermade	20-2900-41M	1 µmol	122.00	117.00	105.00	-
3'-Propargyl-5-Me-dC CPG	20-2982-01	0.1 g	273.00	262.00	235.00	-
3'-Propargyl-5-Me-dC CPG	20-2982-10	1 g	2,682.00	2,575.00	2,307.00	-
3'-Propargyl-5-Me-dC CPG, column	20-2982-41	4 x 1 µmol	546.00	524.00	470.00	-
3'-Propargyl-5-Me-dC CPG, column	20-2982-42	4 x 0.2 µmol	216.00	207.00	186.00	-
3'-Protected Biotin Serinol CPG	20-2993-01	0.1 g	173.00	166.00	149.00	-
3'-Protected Biotin Serinol CPG	20-2993-10	1 g	1,431.00	1,374.00	1,231.00	-
3'-Protected Biotin Serinol CPG	20-2993-13	1 x 10 µmol	431.00	414.00	371.00	-
3'-Protected Biotin Serinol CPG	20-2993-14	1 x 15 µmol	647.00	621.00	556.00	-
3'-Protected Biotin Serinol CPG	20-2993-41	4 x 1 µmol	288.00	276.00	248.00	-
3'-Protected Biotin Serinol CPG	20-2993-42	4 x 0.2 µmol	173.00	166.00	149.00	-
3'-Protected BiotinLC Serinol CPG	20-2995-01	0.1 g	177.00	177.00	159.00	-
3'-Protected BiotinLC Serinol CPG	20-2995-10	1 g	1,470.00	1,470.00	1,323.00	-
3'-Protected BiotinLC Serinol CPG	20-2995-13	1 x 10 µmol	443.00	443.00	399.00	-
3'-Protected BiotinLC Serinol CPG	20-2995-14	1 x 15 µmol	665.00	665.00	599.00	-
3'-Protected BiotinLC Serinol CPG	20-2995-41	4 x 1 µmol	295.00	295.00	266.00	-
3'-Protected BiotinLC Serinol CPG	20-2995-42	4 x 0.2 µmol	177.00	177.00	159.00	-
3'-Protected BiotinLC Serinol CPG, Expedite	20-2995-42E	4 x 0.2 µmol	177.00	177.00	159.00	-
3'-PT-Amino-Modifier C3 CPG	20-2954-01	0.1 g	137.00	132.00	118.00	-
3'-PT-Amino-Modifier C3 CPG	20-2954-10	1 g	971.00	932.00	835.00	-
3'-PT-Amino-Modifier C3 CPG	20-2954-13	1 x 10 µmol	360.00	346.00	310.00	-
3'-PT-Amino-Modifier C3 CPG	20-2954-14	1 x 15 µmol	539.00	517.00	464.00	-
3'-PT-Amino-Modifier C3 CPG	20-2954-41	4 x 1 µmol	201.00	193.00	173.00	-
3'-PT-Amino-Modifier C3 CPG	20-2954-42	4 x 0.2 µmol	122.00	117.00	105.00	-
3'-PT-Amino-Modifier C3 CPG - 4 x 1 µmol	20-2954-41P	4 x 1 µmol	201.00	193.00	173.00	-
3'-PT-Amino-Modifier C6 CPG	20-2956-01	0.1 g	137.00	132.00	118.00	-
3'-PT-Amino-Modifier C6 CPG	20-2956-10	1 g	971.00	932.00	835.00	-
3'-PT-Amino-Modifier C6 CPG	20-2956-13	1 x 10 µmol	360.00	346.00	310.00	-
3'-PT-Amino-Modifier C6 CPG	20-2956-14	1 x 15 µmol	539.00	517.00	464.00	-
3'-PT-Amino-Modifier C6 CPG	20-2956-41	4 x 1 µmol	201.00	193.00	173.00	-
3'-PT-Amino-Modifier C6 CPG	20-2956-42	4 x 0.2 µmol	122.00	117.00	105.00	-
3'-PT-Amino-Modifier C6 CPG, Expedite	20-2956-41E	4 x 1 µmol	201.00	193.00	173.00	-
3'-PT-Amino-Modifier C6 CPG,Expedite	20-2956-42E	4 x 0.2 µmol	122.00	117.00	105.00	-
3'-PT-Amino-Modifier C6 PS	26-2956-10	1 g	1,474.00	1,415.00	1,268.00	-
3'-PT-Amino-Modifier C6 PS	26-2956-52	10 x 200 nmol	316.00	303.00	272.00	-
3'-PT-Amino-Modifier C6 PS	26-2956-55	10 x 40 nmol	316.00	303.00	272.00	-
3'-Spacer C3 CPG	20-2913-01	0.1 g	101.00	97.00	86.90	-
3'-Spacer C3 CPG	20-2913-10	1 g	690.00	662.00	593.00	-
3'-Spacer C3 CPG	20-2913-13	1 x 10 µmol	259.00	249.00	223.00	-
3'-Spacer C3 CPG	20-2913-14	1 x 15 µmol	403.00	387.00	347.00	-
3'-Spacer C3 CPG	20-2913-41	4 x 1 µmol	144.00	138.00	124.00	-
3'-Spacer C3 CPG	20-2913-41E	1 µmol	144.00	138.00	124.00	-
3'-Spacer C3 CPG	20-2913-42	4 x 0.2 µmol	86.30	82.80	74.20	-
3'-Spacer C3 CPG, Exp.	20-2913-42E	4 x 0.2 µmol	86.30	82.80	74.20	-
3'-TAMRA CPG	20-5910-01	0.1 g	159.00	153.00	137.00	-
3'-TAMRA CPG	20-5910-10	1 g	1,321.00	1,268.00	1,136.00	-
3'-TAMRA CPG	20-5910-41	4 x 1 µmol	221.00	212.00	190.00	-
3'-TAMRA CPG	20-5910-42	4 x 0.2 µmol	106.00	102.00	91.20	-
3'-TAMRA CPG 500	20-5910-42E	4 x 0.2 µmol	106.00	102.00	91.20	-
3'-TAMRA CPG 500 - 4 x 1 µmol	20-5910-41E	4 x 1 µmol	221.00	212.00	190.00	-
3'-TAMRA CPG 500 0.2µM (1pk/4)Mermade	20-5910-42M	106.00	102.00	91.20	-	
3'-TAMRA CPG 500 1µM (1pk/4)Mermade	20-5910-41M	221.00	212.00	190.00	-	
3'-TAMRA PS	26-5910-52	10 x 200 nmol	475.00	456.00	409.00	-
3'-TAMRA PS	26-5910-55	10 x 40 nmol	475.00	456.00	409.00	-
3'-Thiol-Modifier 6 S-S CPG	20-2938-41E	4 x 1 µmol	187.00	180.00	161.00	-
3'-Thiol-Modifier 6 S-S CPG	20-2938-42	4 x 0.2 µmol	115.00	110.00	98.90	-
3'-Thiol-Modifier 6 S-S CPG	20-2938-42E	4 x 0.2 µmol	115.00	110.00	98.90	-
3'-Thiol-Modifier 6 S-S CPG - ABI 3900	20-2938-42A	4 x 0.2 µmol	115.00	110.00	98.90	-
3'-Thiol-Modifier C3 S-S CPG	20-2933-01	0.1 g	113.00	108.00	97.20	-
3'-Thiol-Modifier C3 S-S CPG	20-2933-10	1 g	796.00	764.00	685.00	-
3'-Thiol-Modifier C3 S-S CPG	20-2933-13	1 x 10 µmol	299.00	287.00	257.00	-
3'-Thiol-Modifier C3 S-S CPG	20-2933-14	1 x 15 µmol	465.00	446.00	400.00	-
3'-Thiol-Modifier C3 S-S CPG	20-2933-41	4 x 1 µmol	152.00	146.00	131.00	-

Description	Reference	Scale/Quantity	Price List 2025			
			EUR	CHF	GBP	USD
3'-Thiol-Modifier C3 S-S CPG	20-2933-41E		152.00	146.00	131.00	-
3'-Thiol-Modifier C3 S-S CPG	20-2933-42	4 x 0.2 µmol	83.00	79.70	71.40	-
3'-Thiol-Modifier C3 S-S CPG, ABI	20-2933-42A	4 x 0.2 µmol	83.00	79.70	71.40	-
3'-Thiol-Modifier C3 S-S CPG, Beckman	20-2933-41B		152.00	146.00	131.00	-
3'-Thiol-Modifier C3 S-S CPG, Pharmacia	20-2933-41P		152.00	146.00	131.00	-
3'-Thiol-Modifier C3 S-SCPG 1µmol 4 pc M	20-2933-41M		152.00	146.00	131.00	-
3'-Thiol-Modifier C6 S-S CPG	20-2938-41	4x1.0µm	187.00	180.00	161.00	-
3'-Thiol-Modifier C3 S-SCPG-0.2µm(1pk/4)M	20-2933-42M	4 x 0.2 µmol	83.00	79.70	71.40	-
3'-Uaq Cap CPG	20-2980-01	0.1 g	259.00	249.00	223.00	-
4,5-Dicyanoimidazole (DCI)	30-3060-1K	1000 g	5,531.00	5,310.00	4,757.00	-
4,5-Dicyanoimidazole (DCI)	30-3060-K5	500 g	3,319.00	3,186.00	2,854.00	-
4-Thio-dT CE Phosphoramidite, Expedite	10-1034-02E	0.25 g	1,136.00	1,091.00	977.00	-
4-Thio-dT-CE Phosphoramidite	10-1034-02	0.25 g	1,136.00	1,091.00	977.00	-
4-Thio-dT-CE Phosphoramidite	10-1034-90	100 µmol	460.00	442.00	396.00	-
4-Thio-dT-CE Phosphoramidite	10-1034-95	50 µmol	252.00	242.00	217.00	-
4-Thio-U-TOM-CE Phosphoramidite - Expedite	10-3052-02E	0.25 g	1,402.00	1,346.00	1,206.00	-
5'- DMS(O)MT-Amino-Modifier C6, Expedite	10-1907-02E	0.25 g	324.00	311.00	279.00	-
5' Stearyl Phosphoramidite	10-1979-90	100 µmol	64.70	62.10	55.60	-
5,6-Dihydro-dU-CE Phosphoramidite	10-1550-02	0.25 g	1,071.00	1,028.00	921.00	-
5,6-Dihydro-dU-CE Phosphoramidite	10-1550-90	100 µmol	316.00	303.00	272.00	-
5'-Aldehyde-Modifier C2 Phosphoramidite	10-1933-02	0.25 g	604.00	580.00	519.00	-
5'-Aldehyde-Modifier C2 Phosphoramidite	10-1933-90	100 µmol	137.00	132.00	118.00	-
5'-Aldehyde-Modifier C2 Phosphoramidite, Expedite	10-1933-90E	100 µmol	137.00	132.00	118.00	-
5'-Amino-dT-CE Phosphoramidite	10-1932-02	0.25 g	676.00	649.00	581.00	-
5'-Amino-dT-CE Phosphoramidite	10-1932-90	100 µmol	237.00	228.00	204.00	-
5'-Amino-Modifier 5	10-1905-02	0.25 g	288.00	276.00	248.00	-
5'-Amino-Modifier 5	10-1905-90	100 µmol	86.30	82.80	74.20	-
5'-Amino-Modifier 5	10-1905-90E	100 µmol	86.30	82.80	74.20	-
5'-Amino-Modifier C12	10-1912-02	0.25 g	431.00	414.00	371.00	-
5'-Amino-Modifier C12	10-1912-02E	0.25 g	431.00	414.00	371.00	-
5'-Amino-Modifier C12	10-1912-90	100 µmol	129.00	124.00	111.00	-
5'-Amino-Modifier C12 - Mermade	10-1912-02M	0.25 g	431.00	414.00	371.00	-
5'-Amino-Modifier C12, Expedite - 100 µmol	10-1912-90E	100 µmol	129.00	124.00	111.00	-
5'-Amino-Modifier C12-PDA, Mermade	10-1948-02M	0.25 g	512.00	492.00	440.00	-
5'-Amino-Modifier C3-TFA	10-1923-02	0.25 g	252.00	242.00	217.00	-
5'-Amino-Modifier C3-TFA	10-1923-90	100 µmol	71.90	69.00	61.80	-
5'-Amino-Modifier C3-TFA, Exped.	10-1923-02E	0.25 g	252.00	242.00	217.00	-
5'-Amino-Modifier C6	10-1906-02	0.25 g	177.00	170.00	152.00	-
5'-Amino-Modifier C6	10-1906-90	100 µmol	53.10	51.00	45.70	-
5'-Amino-Modifier C6 - Mermade	10-1906-02M	0.25 g	177.00	170.00	152.00	-
5'-Amino-Modifier C6, Beckman	10-1906-90B	100 µmol	53.10	51.00	45.70	-
5'-Amino-Modifier C6, Expedite - 0.25 g	10-1906-02E	0.25 g	177.00	170.00	152.00	-
5'-Amino-Modifier C6, Expedite - 100 µmol	10-1906-90E	100 µmol	53.10	51.00	45.70	-
5'-Amino-Modifier C6, Pharmacia - 100 µmol	10-1906-90P	100 µmol	53.10	51.00	45.70	-
5'-Amino-Modifier C6-PDA	10-1947-02	0.25 g	144.00	138.00	124.00	-
5'-Amino-Modifier C6-PDA	10-1947-02	0.25 g	144.00	138.00	124.00	-
5'-Amino-Modifier C6-PDA	10-1947-90	100 µmol	43.10	41.40	37.10	-
5'-Amino-Modifier C6-TFA	10-1916-02	0.25 g	88.50	85.00	76.10	-
5'-Amino-Modifier C6-TFA	10-1916-90	100 µmol	26.60	25.50	22.90	-
5'-Amino-Modifier C6-TFA, Exped.	10-1916-02E	0.25 g	88.50	85.00	76.10	-
5'-Amino-Modifier C6-TFA, Mermade	10-1916-02M	0.25 g	88.50	85.00	76.10	-
5'-Amino-Modifier TEG CE-Phosphoramidite	10-1917-02	0.25 g	719.00	690.00	618.00	-
5'-Amino-Modifier TEG CE-Phosphoramidite	10-1917-90	100 µmol	165.00	158.00	142.00	-
5'-Amino-Modifier TEG CE-Phosphoramidite, Expedite	10-1917-02E	0.25 g	719.00	690.00	618.00	-
5'-Amino-Modifier TEG CE-Phosphoramidite, Expedite	10-1917-90E	100 µmol	165.00	158.00	142.00	-
5'-AminoOxy - Modifier 11 - CE - Phosphoramidite	10-1919-95	50 µmol	216.00	207.00	186.00	-
5'-AminoOxy-Modifier 11-CE-Phosphoramidite	10-1919-02	0.25 g	1,337.00	1,284.00	1,150.00	-
5'-AminoOxy-Modifier 11-CE-Phosphoramidite	10-1919-90	100 µmol	403.00	387.00	347.00	-
5'-AquaPhluor® 593 CE Phosphor.	10-5923-95	50µmol	647.00	621.00	556.00	-
5'-AquaPhluor® 593 CE Phosphoramidite	10-5923-90	100 µmol	1,294.00	1,242.00	1,113.00	-
5'-AquaPhluor® 639 CE Phosphoramidite	10-5926-90		805.00	773.00	692.00	-
5'-BBQ-650®-CE Phosphoramidite, Expedite	10-5934-02E	0.25 g	1,956.00	1,878.00	1,682.00	-
5'-BBQ-650®-CE Phosphoramidite, Expedite	10-5934-90E	100 µmol	647.00	621.00	556.00	-
5'-BBQ-650®-CE Phosphoramidite, Expedite	10-5934-95E	50 µmol	345.00	331.00	297.00	-
5'-BBQ-650-CE Phosphoramidite	10-5934-90	100 µmol	647.00	621.00	556.00	-
5'-BHQ-1 Phosphoramidite	10-5931-02	0.25 g	1,373.00	1,318.00	1,181.00	-
5'-BHQ-1 Phosphoramidite	10-5931-90	100 µmol	388.00	372.00	334.00	-
5'-BHQ-1 Phosphoramidite	10-5931-95	50 µmol	209.00	201.00	180.00	-
5'-BHQ-2 Phosphoramidite	10-5932-02	0.25 g	1,373.00	1,318.00	1,181.00	-
5'-BHQ-2 Phosphoramidite	10-5932-90	100 µmol	388.00	372.00	334.00	-
5'-BHQ-2 Phosphoramidite	10-5932-95	50 µmol	209.00	201.00	180.00	-
5'-Biotin Phosphoramidite	10-5950-02	0.25 g	277.00	266.00	238.00	-
5'-Biotin Phosphoramidite	10-5950-90	100 µmol	99.60	95.60	85.70	-
5'-Biotin Phosphoramidite	10-5950-95	50 µmol	66.40	63.70	57.10	-
5'-Biotin Phosphoramidite, Exp.	10-5950-02E	0.25 g	277.00	266.00	238.00	-
5'-Biotin Phosphoramidite, Exp.	10-5950-90E	100 µmol	99.60	95.60	85.70	-
5'-Biotin Phosphoramidite, Expedite	10-5950-95E	50 µmol	66.40	63.70	57.10	-
5'-Biotin Phosphoramidite, Pharmacia - 50 µmol	10-5950-95P	50 µmol	66.40	63.70	57.10	-
5-Br-dC-CE Phosphoramidite	10-1080-02	0.25 g	288.00	276.00	248.00	-
5-Br-dC-CE Phosphoramidite	10-1080-90	100 µmol	115.00	110.00	98.90	-
5-Br-dC-CE Phosphoramidite, Expedite - 100 µmol	10-1080-90E	100 µmol	115.00	110.00	98.90	-
5-Br-dU-CE Phosphoramidite	10-1090-02	0.25 g	177.00	170.00	152.00	-
5-Br-dU-CE Phosphoramidite	10-1090-02M	0.25 g	182.00	182.00	164.00	-
5-Br-dU-CE Phosphoramidite	10-1090-90	100 µmol	66.40	63.70	57.10	-

Description	Reference	Scale/Quantity	Price List 2025			
			EUR	CHF	GBP	USD
5-Br-dU-CE Phosphoramidite	10-1090-90M	100 µmol	68.00	68.00	61.00	-
5-Br-dU-CE Phosphoramidite, Expedite	10-1090-02E	0.25 g	177.00	170.00	152.00	-
5 ¹ -Bromoethyl Phosphoramidite, Mermade	10-1946-02M	0.25 g	288.00	276.00	248.00	-
5 ¹ -Bromoethyl Phosphoramidite	10-1946-02	0.25 g	288.00	276.00	248.00	-
5 ¹ -Bromoethyl Phosphoramidite	10-1946-90	100 µmol	86.30	82.80	74.20	-
5 ¹ -Bromoethyl Phosphoramidite, Expedite	10-1946-02E	0.25 g	288.00	276.00	248.00	-
5-Carboxy-dC-CE Phosphoramidite	10-1066-02	0.25 g	2,668.00	2,561.00	2,294.00	-
5-Carboxy-dC-CE Phosphoramidite	10-1066-90	100 µmol	978.00	939.00	841.00	-
5-Carboxy-dC-CE Phosphoramidite	10-1066-95	50 µmol	503.00	483.00	433.00	-
5 ¹ -Carboxy-Modifier C10	10-1935-02	0.25 g	381.00	366.00	328.00	-
5 ¹ -Carboxy-Modifier C10	10-1935-90	100 µmol	93.50	89.80	80.40	-
5 ¹ -Carboxy-Modifier C5	10-1945-02	0.25 g	554.00	532.00	476.00	-
5 ¹ -Carboxy-Modifier C5	10-1945-90	100 µmol	158.00	152.00	136.00	-
5 ¹ -Carboxy-Modifier C5 - Mermade	10-1945-02M	0.25 g	554.00	532.00	476.00	-
5 ¹ -Carboxy-Modifier C5 - Mermade	10-1945-90M	100 µmol	158.00	152.00	136.00	-
5 ¹ -CDPI3 MGB™ Phosphoramidite	10-5924-90	100 µmol	2,085.00	2,002.00	1,793.00	-
5 ¹ -Cholesteryl-TEG Phospho.	10-1976-02	0.25 g	856.00	822.00	736.00	-
5 ¹ -Cholesteryl-TEG Phospho.	10-1976-02	0.25 g	856.00	822.00	736.00	-
5 ¹ -Cholesteryl-TEG Phospho.	10-1976-90	100 µmol	288.00	276.00	248.00	-
5 ¹ -Cholesteryl-TEG Phospho.	10-1976-95	50 µmol	158.00	152.00	136.00	-
5 ¹ -Cholesteryl-TEG Phosphoramidite, Mermade	10-1976-90M	100 µmol	288.00	276.00	248.00	-
5 ¹ -Dabcytl Phosphoramidite	10-5912-02	0.25 g	935.00	898.00	804.00	-
5 ¹ -Dabcytl Phosphoramidite	10-5912-90	100 µmol	324.00	311.00	279.00	-
5 ¹ -Dabcytl Phosphoramidite	10-5912-95	50 µmol	180.00	173.00	155.00	-
5 ¹ -Dabcytl Phosphoramidite, Expedite - 50 µmol	10-5912-95E	50 µmol	180.00	173.00	155.00	-
5 ¹ -Dabcytl Phosphoramidite, Pharmacia - 100 µmol	10-5912-90P	100 µmol	324.00	311.00	279.00	-
5 ¹ -Dabsyl-dT	10-1532-02	0.25 g	1,352.00	1,298.00	1,163.00	-
5 ¹ -Dabsyl-dT	10-1532-90	100 µmol	410.00	394.00	353.00	-
5 ¹ -DBCO-TEG Phosphoramidite	10-1941-95	50 µmol	201.00	193.00	173.00	-
5 ¹ -DBCO-TEG Phosphoramidite, Expedite	10-1941-02E	0.25 g	1,309.00	1,257.00	1,126.00	-
5 ¹ -DBCO-TEG Phosphoramidite, Mermade	10-1941-02M	0.25 g	1,309.00	1,257.00	1,126.00	-
5 ¹ -DBCO-TEG Phosphoramidite, Mermade	10-1941-90M	100 µmol	374.00	359.00	322.00	-
5 ¹ -Dichloro-dimethoxy-Fluorescein Phosphoramidite II (JOE II)	10-5906-02	0.25 g	712.00	684.00	612.00	-
5 ¹ -Dichloro-dimethoxy-Fluorescein Phosphoramidite II (JOE II)	10-5906-90	100 µmol	285.00	274.00	245.00	-
5 ¹ -Dimethoxytrityl-N-benzoyl-5-cyanoethoxy- methyl-2'-deoxyCytidine,3'-[(2-cyanoethyl)- (N,N-diisopropyl)]-phospho	10-1062-02	0.25 g	2,790.00	2,678.00	2,399.00	-
5 ¹ -Dimethoxytrityl-N-benzoyl-5-cyanoethoxy- methyl-2'-deoxyCytidine,3'-[(2-cyanoethyl)- (N,N-diisopropyl)]-phospho	10-1062-90	100 µmol	1,035.00	994.00	890.00	-
5 ¹ -Dimethoxytrityl-N-benzoyl-5-cyanoethoxy- methyl-2'-deoxyCytidine,3'-[(2-cyanoethyl)- (N,N-diisopropyl)]-phospho	10-1062-95	50 µmol	532.00	511.00	458.00	-
5 ¹ -DMS(O)MT-Amino-Modifier C6	10-1907-02	0.25 g	324.00	311.00	279.00	-
5 ¹ -DMS(O)MT-Amino-Modifier C6	10-1907-90	100 µmol	115.00	110.00	98.90	-
5 ¹ -DMS(O)MT-Amino-Modifier C6, Expedite - 100 µmol	10-1907-90E	100 µmol	115.00	110.00	98.90	-
5-Ethynyl-dU-CE Phosphoramidite	10-1554-02	0.25 g	1,345.00	1,291.00	1,157.00	-
5-Ethynyl-dU-CE Phosphoramidite	10-1554-90	100 µmol	410.00	394.00	353.00	-
5-F-dU-CE Phosphoramidite	10-1092-02	0.25 g	662.00	636.00	569.00	-
5-F-dU-CE Phosphoramidite	10-1092-90	100 µmol	216.00	207.00	186.00	-
5-F-dU-CE Phosphoramidite, Exp	10-1092-02E	0.25 g	662.00	636.00	569.00	-
5-F-dU-CE Phosphoramidite, Exp	10-1092-90E	100 µmol	216.00	207.00	186.00	-
5 ¹ -Fluorescein amidite (6-FAM)	10-5901-90E		83.00	79.70	71.40	-
5 ¹ -Fluorescein amidite (6-FAM)	10-5901-95E		55.30	53.10	47.60	-
5 ¹ -Fluorescein Phosphoramidite	10-5901-02	0.25 g	210.00	202.00	181.00	-
5 ¹ -Fluorescein Phosphoramidite	10-5901-90	100 µmol	83.00	79.70	71.40	-
5 ¹ -Fluorescein Phosphoramidite	10-5901-90P	100 µmol	83.00	79.70	71.40	-
5 ¹ -Fluorescein Phosphoramidite	10-5901-95	50 µmol	55.30	53.10	47.60	-
5 ¹ -Fluorescein Phosphoramidite, Expedite	10-5901-02E	0.25 g	210.00	202.00	181.00	-
5 ¹ -Fluorescein Phosphoramidite, Mermade	10-5901-02M	0.25 g	210.00	202.00	181.00	-
5-Formyl dC III CE Phosphoramidite	10-1564-90	100 µmol	1,956.00	1,878.00	1,682.00	-
5-Formyl-dC-CE Phosphoramidite	10-1514-02	0.25 g	4,638.00	4,452.00	3,989.00	-
5-Formyl-dC-CE Phosphoramidite	10-1514-95	50 µmol	877.00	842.00	754.00	-
5-Formylindole-CE Phosphoramidite	10-1934-02	0.25 g	575.00	552.00	495.00	-
5-Formylindole-CE Phosphoramidite	10-1934-90	100 µmol	180.00	173.00	155.00	-
5-Formylindole-CE Phosphoramidite, Expedite	10-1934-90E	100 µmol	180.00	173.00	155.00	-
5-Formylindole-CE Phosphoramidite, Mermade	10-1934-02M	0.25 g	575.00	552.00	495.00	-
5 ¹ -GalNAc C3 Phosphoramidite	10-1974-02	0.25 g	762.00	732.00	655.00	-
5 ¹ -GalNAc C3 Phosphoramidite	10-1974-90	100 µmol	367.00	352.00	316.00	-
5 ¹ -GalNAc C3 Phosphoramidite, Mermade	10-1974-02M	0.25 g	762.00	732.00	655.00	-
5 ¹ -GalNAc C3 Phosphoramidite, Mermade	10-1974-90M	100 µmol	367.00	352.00	316.00	-
5 ¹ -Hexachloro-Fluorescein Phosph.	10-5902-02E	0.25 g	830.00	797.00	714.00	-
5 ¹ -Hexachloro-Fluorescein Phosphoramidite	10-5902-02	0.25 g	830.00	797.00	714.00	-
5 ¹ -Hexachloro-Fluorescein Phosphoramidite	10-5902-90	100 µmol	365.00	350.00	314.00	-
5 ¹ -Hexachloro-Fluorescein Phosphoramidite	10-5902-95	50 µmol	188.00	180.00	162.00	-
5 ¹ -HEX-Fluo amidite	10-5902-90E	100 µmol	365.00	350.00	314.00	-
5 ¹ -HEX-Fluo amidite, Expedite - 50 µmol	10-5902-95E	50 µmol	188.00	180.00	162.00	-
5 ¹ -Hexynyl Phosphoramidite	10-1908-02	0.25 g	288.00	276.00	248.00	-
5 ¹ -Hexynyl Phosphoramidite	10-1908-90	100 µmol	86.30	82.80	74.20	-
5 ¹ -Hexynyl Phosphoramidite, Expedite	10-1908-02E	0.25 g	288.00	276.00	248.00	-
5 ¹ -Hexynyl Phosphoramidite, Mermade	10-1908-02M	0.25 g	288.00	276.00	248.00	-
5-Hydroxymethyl dC II CE Phosphoramidite	10-1510-90	100 µmol	964.00	925.00	829.00	-
5-Hydroxymethyl-dC II-CE Phosphoramidite	10-1510-95	50 µmol	496.00	476.00	427.00	-
5-Hydroxymethyl-dC II-CE Phosphoramidite – 0.25 g	10-1510-02	0.25 g	3,020.00	2,899.00	2,597.00	-
5-Hydroxymethyl-dU-CE Phosphor.	10-1093-90M	100 µmol	403.00	387.00	347.00	-
5-Hydroxymethyl-dU-CE Phosphoramidite	10-1093-02	0.25 g	1,230.00	1,181.00	1,058.00	-
5-Hydroxymethyl-dU-CE Phosphoramidite	10-1093-90	100 µmol	403.00	387.00	347.00	-
5-I-dC-CE Phosphoramidite	10-1081-02	0.25 g	561.00	539.00	482.00	-
5-I-dC-CE Phosphoramidite	10-1081-90	100 µmol	230.00	221.00	198.00	-
5-I-dC-CE Phosphoramidite, Expedite - 100 µmol	10-1081-90E	100 µmol	230.00	221.00	198.00	-

Description	Reference	Scale/Quantity	Price List 2025			
			EUR	CHF	GBP	USD
5'-I-dT-CE Phosphoramidite	10-1931-02	0.25 g	496.00	476.00	427.00	-
5'-I-dT-CE Phosphoramidite	10-1931-90	100 µmol	129.00	124.00	111.00	-
5-I-dU-CE Phosphoramidite	10-1091-02	0.25 g	266.00	255.00	229.00	-
5-I-dU-CE Phosphoramidite	10-1091-90	100 µmol	93.50	89.80	80.40	-
5-I-dU-CE Phosphoramidite, Expedite	10-1091-02E	0.25 g	266.00	255.00	229.00	-
5-I-dU-CE Phosphoramidite, Expedite	10-1091-90E	100 µmol	93.50	89.80	80.40	-
5'-Maleimide-Modifier Phosphoramidite	10-1938-02	0.25 g	482.00	463.00	415.00	-
5'-Maleimide-Modifier-Phosphoramidite	10-1938-90	100 µmol	101.00	97.00	86.90	-
5-Me-2'-deoxyZebularine-CE Phosphoramidite	10-1061-02	0.25 g	1,697.00	1,629.00	1,459.00	-
5-Me-2'-deoxyZebularine-CE Phosphoramidite	10-1061-90	100 µmol	618.00	593.00	531.00	-
5-Me-2'-deoxyZebularine-CE Phosphoramidite	10-1061-95	50 µmol	324.00	311.00	279.00	-
5-Me-C-2'-MOE-Phosphoramidite	10-3211-10	1.0 g	108.00	104.00	92.90	-
5-Me-C-2'-MOE-Phosphoramidite - Mermade	10-3211-10M	1.0 g	108.00	104.00	92.90	-
5-Me-C-TOM-CE Phosphoramidite - Expedite	10-3064-90E	100 µmol	273.00	262.00	235.00	-
5-Me-dC-CE Phosphoramidite	10-1060-02	0.25 g	133.00	128.00	114.00	-
5-Me-dC-CE Phosphoramidite	10-1060-90	100 µmol	55.30	53.10	47.60	-
5-Me-dC-CE Phosphoramidite	10-1060-90E	100 µmol	55.30	53.10	47.60	-
5-Me-dC-CE Phosphoramidite, Expedite - 0.25 g	10-1060-02E	0.25 g	133.00	128.00	114.00	-
5-Me-U-2'-MOE-Phosphoramidite	10-3231-10	1.0 g	108.00	104.00	92.90	-
5-Me-U-2'-MOE-Phosphoramidite, Mermade	10-3231-20M	2.0 g	216.00	207.00	186.00	-
5-Nitroindole-CE phosphoramidi	10-1044-90C	100 µmol	187.00	180.00	161.00	-
5-Nitroindole-CE Phosphoramidite	10-1044-02	0.25 g	518.00	497.00	445.00	-
5-Nitroindole-CE Phosphoramidite	10-1044-90	100 µmol	187.00	180.00	161.00	-
5-Nitroindole-CE Phosphoramidite, Expedite - 0.25 g	10-1044-02E	0.25 g	518.00	497.00	445.00	-
5-Nitroindole-CE Phosphoramidite, Expedite - 100 µmol	10-1044-90E	100 µmol	187.00	180.00	161.00	-
5-O-Dimethoxytrityl-1-O-tert-butylidimethylsilyl-2-deoxyribose-3-[(2-cyanoethyl)-(N,N-diisopropyl)]-phospho	10-1927-02	250 mg	733.00	704.00	630.00	-
5-O-Dimethoxytrityl-1-O-tert-butylidimethylsilyl-2-deoxyribose-3-[(2-cyanoethyl)-(N,N-diisopropyl)]-phospho	10-1927-90	100 µmol	244.00	234.00	210.00	-
5-O-Dimethoxytrityl-1-O-tert-butylidimethylsilyl-2-deoxyribose-3-[(2-cyanoethyl)-(N,N-diisopropyl)]-phospho	10-1927-95	50 µmol	137.00	132.00	118.00	-
5-OH-dC-CE Phosphoramidite	10-1063-02	0.25 g	1,431.00	1,374.00	1,231.00	-
5-OH-dC-CE Phosphoramidite	10-1063-90	100 µmol	554.00	532.00	476.00	-
5-OH-dU-CE Phosphoramidite	10-1053-02	0.25 g	1,445.00	1,387.00	1,243.00	-
5-OH-dU-CE Phosphoramidite	10-1053-90	100 µmol	467.00	448.00	402.00	-
5'-OMe-dT-CE Phosphoramidite	10-1031-02	0.25 g	590.00	566.00	507.00	-
5'-OMe-dT-CE Phosphoramidite	10-1031-90	100 µmol	201.00	193.00	173.00	-
5'-Pyrene Cap Phosphoramidite	10-1987-02	0.25 g	712.00	684.00	612.00	-
5'-Pyrene Cap Phosphoramidite	10-1987-90	100 µmol	280.00	269.00	241.00	-
5'-TCO C6 Phosphoramidite	10-1943-02	0.25 g	1,726.00	1,657.00	1,484.00	-
5'-TET-Fluo amidite	10-5903-90E	100 µmol	409.00	393.00	352.00	-
5'-TET-Fluo amidite, Expedite - 50 µmol	10-5903-95E	50 µmol	210.00	202.00	181.00	-
5'-Tetrachloro-Fluorescein Phosphoramidite	10-5903-02	0.25 g	985.00	946.00	847.00	-
5'-Tetrachloro-Fluorescein Phosphoramidite	10-5903-90	100 µmol	409.00	393.00	352.00	-
5'-Tetrachloro-Fluorescein Phosphoramidite	10-5903-95	50 µmol	210.00	202.00	181.00	-
5'-Thiol-Modifier C6	10-1926-02	0.25 g	288.00	276.00	248.00	-
5'-Thiol-Modifier C6	10-1926-90	100 µmol	86.30	82.80	74.20	-
5'-Thiol-Modifier C6	10-1926-90E	100 µmol	86.30	82.80	74.20	-
5'-Trimethoxystilbene Cap Phosphoramidite	10-1986-02	0.25 g	741.00	711.00	637.00	-
5'-Trimethoxystilbene Cap Phosphoramidite	10-1986-90	100 µmol	280.00	269.00	241.00	-
6-FAM CPG 1000 -1µm (1pk/4)Mermade	20-2961-41M	4 x 1 µmol	256.00	246.00	220.00	-
6-Fluorescein Phosphoramidite	10-1964-02	0.25 g	856.00	822.00	736.00	-
6-Fluorescein Phosphoramidite	10-1964-90	100 µmol	424.00	407.00	365.00	-
6-Fluorescein Phosphoramidite	10-1964-95	50 µmol	237.00	228.00	204.00	-
6-Fluorescein Phosphoramidite, Expedite	10-1964-95E	50 µmol	237.00	228.00	204.00	-
6-Fluorescein Serinol Phosphoramidite	10-1994-02	0.25 g	856.00	822.00	736.00	-
6-Fluorescein Serinol Phosphoramidite	10-1994-90	100 µmol	424.00	407.00	365.00	-
6-Fluorescein Serinol Phosphoramidite	10-1994-95	50 µmol	237.00	228.00	204.00	-
6-thio-dG-CE Phosphoramidite	10-1072-02	0.25 g	1,539.00	1,477.00	1,324.00	-
6-thio-dG-CE Phosphoramidite	10-1072-90	100 µmol	590.00	566.00	507.00	-
6-thio-dG-CE Phosphoramidite	10-1072-95	50 µmol	309.00	297.00	266.00	-
6-Thio-G-CE Phosphoramidite	10-3072-02	0.25 g	1,726.00	1,657.00	1,484.00	-
6-Thio-G-CE Phosphoramidite	10-3072-90	100 µmol	719.00	690.00	618.00	-
6-Thio-G-CE Phosphoramidite	10-3072-95	50 µmol	360.00	346.00	310.00	-
7-deaza-8-aza-dA-CE Phosphoramidite	10-1083-02	0.25 g	1,632.00	1,567.00	1,404.00	-
7-deaza-8-aza-dA-CE Phosphoramidite	10-1083-90	100 µmol	561.00	539.00	482.00	-
7-deaza-8-aza-dA-CE Phosphoramidite	10-1083-95	50 µmol	295.00	283.00	254.00	-
7-Deaza-8-aza-dG-CE Phosphoramidite (PPG)	10-1073-02	0.25 g	2,970.00	2,851.00	2,554.00	-
7-Deaza-8-aza-dG-CE Phosphoramidite (PPG)	10-1073-90	100 µmol	978.00	939.00	841.00	-
7-Deaza-8-aza-dG-CE Phosphoramidite (PPG)	10-1073-95	50 µmol	503.00	483.00	433.00	-
7-Deaza-dA-CE Phosphoramidite	10-1001-02	0.25 g	1,747.00	1,677.00	1,502.00	-
7-Deaza-dA-CE Phosphoramidite	10-1001-90	100 µmol	604.00	580.00	519.00	-
7-Deaza-dA-CE Phosphoramidite	10-1001-95	50 µmol	316.00	303.00	272.00	-
7-deaza-dG-CE Phosphoramidite	10-1021-02	0.25 g	1,510.00	1,450.00	1,299.00	-
7-deaza-dG-CE Phosphoramidite	10-1021-90	100 µmol	532.00	511.00	458.00	-
7-deaza-dG-CE Phosphoramidite	10-1021-95	50 µmol	280.00	269.00	241.00	-
7-Deaza-dG-CE Phosphoramidite	10-1021-95C	50 µmol	280.00	269.00	241.00	-
8-Amino-dA-CE Phosphoramidite	10-1086-02	0.25 g	2,229.00	2,140.00	1,917.00	-
8-Amino-dA-CE Phosphoramidite	10-1086-90	100 µmol	791.00	759.00	680.00	-
8-Amino-dA-CE Phosphoramidite	10-1086-95	50 µmol	410.00	394.00	353.00	-
8-Amino-dG-CE Phosphoramidite	10-1079-02	0.25 g	1,603.00	1,539.00	1,379.00	-
8-Amino-dG-CE Phosphoramidite	10-1079-90	100 µmol	575.00	552.00	495.00	-
8-Amino-dG-CE Phosphoramidite	10-1079-95	50 µmol	309.00	297.00	266.00	-
8-Br-dA-CE Phosphoramidite	10-1007-02	0.25 g	611.00	587.00	525.00	-
8-Br-dA-CE Phosphoramidite	10-1007-90	100 µmol	252.00	242.00	217.00	-
8-Br-dG-CE Phosphoramidite	10-1027-02	0.25 g	467.00	448.00	402.00	-
8-Br-dG-CE Phosphoramidite	10-1027-90	100 µmol	173.00	166.00	149.00	-

Description	Reference	Scale/Quantity	Price List 2025			
			EUR	CHF	GBP	USD
8-Br-dG-CE Phosphoramidite, Expedite	10-1027-02E	0.25 g	467.00	448.00	402.00	-
8-Br-dG-CE Phosphoramidite, Expedite - 100 µmol	10-1027-90E	100 µmol	173.00	166.00	149.00	-
8-oxo-dA-CE Phosphoramidite	10-1008-02	0.25 g	640.00	614.00	550.00	-
8-oxo-dA-CE Phosphoramidite	10-1008-90	100 µmol	230.00	221.00	198.00	-
8-oxo-dG-CE Phosphoramidite	10-1028-02	0.25 g	1,402.00	1,346.00	1,206.00	-
8-oxo-dG-CE Phosphoramidite	10-1028-90	100 µmol	511.00	491.00	439.00	-
8-oxo-dG-CE Phosphoramidite	10-1028-95	50 µmol	255.00	245.00	219.00	-
8-oxo-dG-CE Phosphoramidite, Expedite - 100 µmol	10-1028-90E	100 µmol	511.00	491.00	439.00	-
A-2'-MOE-Phosphoramidite	10-3200-10	1.0 g	108.00	104.00	92.90	-
A-2'-MOE-Phosphoramidite - Mermade	10-3200-10M	1.0 g	108.00	104.00	92.90	-
Abasic II Phosphoramidite - Mermade	10-1927-02M	250 mg	733.00	704.00	630.00	-
Ac-5-Me-dC-CE Phosphoramidite	10-1560-02	0.25 g	173.00	166.00	149.00	-
Ac-5-Me-dC-CE Phosphoramidite	10-1560-90	100 µmol	71.90	69.00	61.80	-
Ac-A-RNA-CPG	20-3404-42A	4 x 0.2 µmol	108.00	104.00	92.90	-
Ac-A-RNA-CPG, Expedite	20-3404-41E	4 x 1.0 µmol	144.00	138.00	124.00	-
ACC Trimer Phosphoramidite	13-1011-90	100 µmol	1,294.00	1,242.00	1,113.00	-
Ac-C-CE Phosphoramidite - Expedite	10-3015-10E	1 g	99.60	95.60	85.70	-
Ac-C-CE Phosphoramidite, Pharmacia	10-3015-05P	0.5 g	53.10	51.00	45.70	-
Ac-dC-5'-CE Phosphoramidite.	10-5101-10	1.0 g	374.00	359.00	322.00	-
Ac-dC-CE Phosphoramidite	10-1015-5S	5 g	77.40	74.30	66.60	-
Ac-dC-CPG 500	20-2013-10	1 g	66.40	63.70	57.10	-
Ac-dC-Me Phosphoramidite	10-1115-02	0.25 g	395.00	379.00	340.00	-
Ac-dC-Me Phosphoramidite	10-1115-05	0.5 g	791.00	759.00	680.00	-
Ac-dC-Me Phosphoramidite	10-1115-05E	0.5 g	791.00	759.00	680.00	-
Ac-dC-Q-CPG 500	21-2013-01	0.1 g	33.20	31.90	28.60	-
Ac-dC-Q-CPG 500	21-2013-10	1 g	277.00	266.00	238.00	-
Ac-dC-Q-CPG 500	21-2113-41	4 x 1 µmol	66.40	63.70	57.10	-
Ac-dC-Q-CPG 500	21-2113-42	4 x 0.2 µmol	44.30	42.50	38.10	-
Ac-dC-Q-CPG 500	21-2213-41	4 x 1 µmol	66.40	63.70	57.10	-
Ac-dC-Q-CPG 500	21-2213-42	4 x 0.2 µmol	44.30	42.50	38.10	-
Ac-dC-Q-CPG 500,	21-2013-02	0.25 g	66.40	63.70	57.10	-
Ac-G-CE Phosphoramidite	10-3025-02E	0.25 g	35.40	34.00	30.40	-
Ac-G-CE Phosphoramidite	10-3025-05	0.5 g	53.10	51.00	45.70	-
Ac-G-CE Phosphoramidite - Expedite - 0.5 g	10-3025-05E	0.5 g	53.10	51.00	45.70	-
Acridine CPG	20-2973-41E	1 µmol	288.00	276.00	248.00	-
Acridine CPG	20-2973-42E	0.2 µmol	173.00	166.00	149.00	-
Acridine Phosphoramidite	10-1973-02	0.25 g	971.00	932.00	835.00	-
Acridine Phosphoramidite	10-1973-90	100 µmol	424.00	407.00	365.00	-
Acridine Phosphoramidite	10-1973-95	50 µmol	237.00	228.00	204.00	-
AGA Trimer Phosphoramidite	13-1020-90	100 µmol	1,294.00	1,242.00	1,113.00	-
Alkyne-Modifier Serinol Phosphor.	10-1992-90M		266.00	255.00	229.00	-
Alkyne-Modifier Serinol Phosphoramidite	10-1992-02	0.25 g	827.00	794.00	711.00	-
Alkyne-Modifier Serinol Phosphoramidite	10-1992-90	100 µmol	266.00	255.00	229.00	-
Alkyne-Modifier Serinol Phosphoramidite	10-1992-95	50 µmol	144.00	138.00	124.00	-
Alkyne-Modifier Serinol Phosphoramidite, Mermade	10-1992-02M	0.25 g	827.00	794.00	711.00	-
Alkyne-NHS Ester	50-1905-24	23 mg	431.00	414.00	371.00	-
Alkyne-NHS Ester, 2.3mg	50-1905-23		86.30	82.80	74.20	-
Amino-Modifier C2 dT	10-1037-02	0.25 g	604.00	580.00	519.00	-
Amino-Modifier C2 dT	10-1037-05	0.5 g	1,208.00	1,160.00	1,039.00	-
Amino-Modifier C2 dT	10-1037-90	100 µmol	273.00	262.00	235.00	-
Amino-Modifier C6 dA	10-1089-02	0.25 g	762.00	732.00	655.00	-
Amino-Modifier C6 dA	10-1089-90	100 µmol	331.00	318.00	285.00	-
Amino-Modifier C6 dA - Mermade	10-1089-02M	0.25 g	762.00	732.00	655.00	-
Amino-Modifier C6 dC	10-1019-02	0.25 g	928.00	891.00	798.00	-
Amino-Modifier C6 dC	10-1019-90	100 µmol	395.00	379.00	340.00	-
Amino-Modifier C6 dC, Expedite - 100 µmol	10-1019-90E	100 µmol	395.00	379.00	340.00	-
Amino-Modifier C6 dC, Mermade	10-1019-02M	0.25 g	928.00	891.00	798.00	-
Amino-Modifier C6 dT	10-1039-02	0.25 g	358.00	344.00	308.00	-
Amino-Modifier C6 dT	10-1039-02P	0.25 g	358.00	344.00	308.00	-
Amino-Modifier C6 dT	10-1039-05	0.5 g	717.00	688.00	617.00	-
Amino-Modifier C6 dT	10-1039-90	100 µmol	179.00	172.00	154.00	-
Amino-Modifier C6 dT, Expedite	10-1039-90E	100 µmol	179.00	172.00	154.00	-
Amino-Modifier C6-U Phosphoramidite	10-3039-02	0.25 g	2,747.00	2,637.00	2,362.00	-
Amino-Modifier C6-U Phosphoramidite	10-3039-90	100 µmol	1,323.00	1,270.00	1,138.00	-
Amino-Modifier C6-U Phosphoramidite	10-3039-95	50 µmol	683.00	656.00	587.00	-
Amino-Modifier Serinol amidite	10-1997-02E	0.25 g	856.00	822.00	736.00	-
Amino-Modifier Serinol Phosphoramidite	10-1997-02	0.25 g	856.00	822.00	736.00	-
Amino-Modifier Serinol Phosphoramidite	10-1997-90	100 µmol	324.00	311.00	279.00	-
Amino-Modifier Serinol Phosphoramidite	10-1997-95	50 µmol	180.00	173.00	155.00	-
AP-dC-CE Phosphoramidite	10-1097-02	0.25 g	1,855.00	1,781.00	1,595.00	-
AP-dC-CE Phosphoramidite	10-1097-90	100 µmol	733.00	704.00	630.00	-
AP-dC-CE Phosphoramidite	10-1097-95	50 µmol	388.00	372.00	334.00	-
AP-dC-CE Phosphoramidite, Mermade	10-1097-02M	0.25 g	1,855.00	1,781.00	1,595.00	-
Asymmetric Doubler (Lev) Phosphoramidite, Expedite	10-1981-90E	100 µmol	151.00	145.00	130.00	-
a-Tocopherol-TEG Phosphoramidite	10-1977-02	0.25 g	949.00	911.00	816.00	-
a-Tocopherol-TEG Phosphoramidite	10-1977-90	100 µmol	475.00	456.00	409.00	-
a-Tocopherol-TEG Phosphoramidite	10-1977-95	50 µmol	252.00	242.00	217.00	-
Azidobutyrate NHS Ester	50-1904-23	2.3 mg	86.30	82.80	74.20	-
Azidobutyrate NHS Ester	50-1904-24	23 mg	431.00	414.00	371.00	-
Azobenzene Phosphoramidite	10-5800-02	0.25 g	791.00	759.00	680.00	-
Azobenzene Phosphoramidite	10-5800-90	100 µmol	288.00	276.00	248.00	-
Azobenzene Phosphoramidite	10-5800-95	50 µmol	151.00	145.00	130.00	-
Azobenzene Phosphoramidite, Expedite	10-5800-02E	0.25 g	791.00	759.00	680.00	-
BBQ-650° CPG	20-5934-01	0.1 g	410.00	394.00	353.00	-

Description	Reference	Scale/Quantity	Price List 2025			
			EUR	CHF	GBP	USD
BBQ-650® CPG	20-5934-10	1 g	3,473.00	3,334.00	2,987.00	-
BBQ-650® CPG, Expedite	20-5934-41E	4 x 1 µmol	489.00	469.00	421.00	-
BBQ-650®-dT-CE Phosphoramidite	10-5944-02	0.25 g	2,150.00	2,064.00	1,849.00	-
BBQ-650®-dT-CE Phosphoramidite	10-5944-90	100 µmol	1,251.00	1,201.00	1,076.00	-
BBQ-650®-dT-CE Phosphoramidite	10-5944-95	50 µmol	640.00	614.00	550.00	-
BBQ-650®-dT-CE Phosphoramidite - Expedite	10-5944-02E	0.25 g	2,150.00	2,064.00	1,849.00	-
beta-L-Pac-dA-CE Phosphoramidite	10-2101-02	0.25 g	216.00	207.00	186.00	-
BHQ-1-dT	10-5941-02	0.25 g	1,920.00	1,843.00	1,651.00	-
BHQ-1-dT	10-5941-90	100 µmol	1,064.00	1,021.00	915.00	-
BHQ-1-dT	10-5941-95	50 µmol	532.00	511.00	458.00	-
BHQ-1-dT, Mermade	10-5941-02M	0.25 g	1,920.00	1,843.00	1,651.00	-
BHQ-2 Phosphoramidite	10-1962-90	100 µmol	762.00	732.00	655.00	-
BHQ-2-dT	10-5942-02	0.25 g	1,920.00	1,843.00	1,651.00	-
BHQ-2-dT	10-5942-90	100 µmol	1,064.00	1,021.00	915.00	-
BHQ-2-dT	10-5942-95	50 µmol	532.00	511.00	458.00	-
BHQ-2-dT, Expedite	10-5942-95E	50 µmol	532.00	511.00	458.00	-
BHQ-2-dT, Mermade	10-5942-02M	0.25 g	1,920.00	1,843.00	1,651.00	-
BioTEG-CPG	20-2955-42E	0.2 µmol	133.00	128.00	114.00	-
BioTEG-CPG -0.2µmol col.(1pk/4)Mermade	20-2955-42M	0.2µmol	133.00	128.00	114.00	-
BioTEG-CPG, column	20-2955-41E	1 µmol	243.00	233.00	209.00	-
BioTEG-CPG-1.0 µmol col.(1pk/4)Mermade	20-2955-41M	1 µmol	243.00	233.00	209.00	-
Biotin Phosphoramidite	10-1953-02	0.25 g	971.00	932.00	835.00	-
Biotin Phosphoramidite	10-1953-90	100 µmol	424.00	407.00	365.00	-
Biotin Phosphoramidite	10-1953-95	50 µmol	237.00	228.00	204.00	-
Biotin Phosphoramidite - Expedite	10-1953-90E	100 µmol	424.00	407.00	365.00	-
Biotin Phosphoramidite, Expedite.	10-1953-02E	0.25 g	971.00	932.00	835.00	-
Biotin Phosphoramidite, Mermade	10-1953-02M	0.25 g	971.00	932.00	835.00	-
Biotin-dT	10-1038-02	0.25 g	830.00	797.00	714.00	-
Biotin-dT	10-1038-90	100 µmol	431.00	414.00	371.00	-
Biotin-dT	10-1038-95	50 µmol	222.00	213.00	191.00	-
Biotin-dT - Expedite - 100 µmol	10-1038-90E	100 µmol	431.00	414.00	371.00	-
BiotinTEG Azide	50-2000-90	100 µmol	647.00	621.00	556.00	-
BiotinTEG Azide	50-2000-92	25 µmol	216.00	207.00	186.00	-
BiotinTEG Phosphoramidite	10-1955-02	0.25 g	523.00	502.00	450.00	-
BiotinTEG Phosphoramidite	10-1955-90	100 µmol	228.00	219.00	196.00	-
BiotinTEG Phosphoramidite	10-1955-95	50 µmol	128.00	123.00	110.00	-
BiotinTEG Phosphoramidite - Expedite	10-1955-90E	100 µmol	228.00	219.00	196.00	-
BiotinTEG Phosphoramidite, Exp.	10-1955-02E	0.25 g	523.00	502.00	450.00	-
BiotinTEG Phosphoramidite, Mermade	10-1955-02M	0.25 g	523.00	502.00	450.00	-
Bz-5-Me-C-L-A-CE Phosphoramidite, Expedite	10-2011-10E	1 g	216.00	207.00	186.00	-
Bz-5-Me-C-L-A-CE Phosphoramidite	10-2011-10	1 g	216.00	207.00	186.00	-
Bz-5-Me-C-L-A-CE Phosphoramidite - Mermade	10-2011-05M	0.5 g	108.00	104.00	92.90	-
Bz-5-Me-C-L-A-CE Phosphoramidite - Mermade	10-2011-10M	1 g	216.00	207.00	186.00	-
Bz-5-Me-C-L-A-CPG	20-2511-41E	4 x 1.0 µmol	93.50	89.80	80.40	-
Bz-A-CE Phosphoramidite	10-3003-02E	0.25g	35.40	34.00	30.40	-
Bz-A-LA Phosphoramidite	10-2000-05	0.5 g	108.00	104.00	92.90	-
Bz-A-LA-CE Phosphoramidite	10-2000-10	1 g	216.00	207.00	186.00	-
Bz-A-LA-CE Phosphoramidite - Mermade	10-2000-05M	0.5 g	108.00	104.00	92.90	-
Bz-A-LA-CE Phosphoramidite - Mermade	10-2000-10M	1 g	216.00	207.00	186.00	-
Bz-A-LA-CE Phosphoramidite, Expedite	10-2000-10E	1 g	216.00	207.00	186.00	-
Bz-A-LA-CPG	20-2501-41E	4 x 1.0 µmol	93.50	89.80	80.40	-
C8-Alkyne-dC-CE Phosphoramidite	10-1543-02E		1,798.00	1,726.00	1,546.00	-
C8-Alkyne-dC-CE Phosphoramidite	10-1543-90	100 µmol	683.00	656.00	587.00	-
C8-Alkyne-dC-CE Phosphoramidite	10-1543-90E	100 µmol	683.00	656.00	587.00	-
C8-Alkyne-dT-CE Phosphoramidite	10-1540-02E	0.25 g	1,294.00	1,242.00	1,113.00	-
C8-Alkyne-dT-CE Phosphoramidite	10-1540-90	100 µmol	453.00	435.00	390.00	-
C8-Alkyne-dT-CE Phosphoramidite	10-1540-90E	100 µmol	453.00	435.00	390.00	-
Carboxy-dT	10-1035-02	0.25 g	604.00	580.00	519.00	-
Carboxy-dT	10-1035-90	100 µmol	288.00	276.00	248.00	-
CCA Trimer Phosphoramidite	13-1110-90	100 µmol	1,294.00	1,242.00	1,113.00	-
CDPI3 MGB™ CPG – 1 g	20-5924-10	1 g	3,221.00	3,092.00	2,770.00	-
CDPI3 MGB™ CPG, Expedite	20-5924-01	0.1 g	338.00	324.00	291.00	-
CGG Trimer Phosphoramidite	13-1122-90	100 µmol	1,294.00	1,242.00	1,113.00	-
Chemical Phosphoryl. Reagent II (CPR II)	10-1901-02E	0.25 g	221.00	212.00	190.00	-
Chemical Phosphorylation Reagent	10-1900-02	0.25 g	177.00	170.00	152.00	-
Chemical Phosphorylation Reagent	10-1900-90	100 µmol	49.80	47.80	42.80	-
Chemical Phosphorylation Reagent II	10-1901-02	0.25 g	221.00	212.00	190.00	-
Chemical Phosphorylation Reagent II	10-1901-90	100 µmol	66.40	63.70	57.10	-
Chemical Phosphorylation Reagent II (CPR II)	10-1901-90E		66.40	63.70	57.10	-
Cholesteryl-TEG CPG	20-2975-42E	0.2 µmol	121.00	116.00	104.00	-
Cholesteryl-TEG Phosphoramidite	10-1975-02	0.25 g	856.00	822.00	736.00	-
Cholesteryl-TEG Phosphoramidite	10-1975-02E	0.25 g	856.00	822.00	736.00	-
Cholesteryl-TEG Phosphoramidite	10-1975-90	100 µmol	446.00	428.00	384.00	-
Cholesteryl-TEG Phosphoramidite	10-1975-95	50 µmol	237.00	228.00	204.00	-
Cholesteryl-TEG Phosphoramidite, Mermade	10-1975-02M	0.25 g	856.00	822.00	736.00	-
Cis-syn Thymine Dimer Phosphoramidite	11-1330-02	0.25 g	20,017.00	20,017.00	18,015.00	-
Cis-syn Thymine Dimer Phosphoramidite	11-1330-90	100 µmol	8,017.00	7,696.00	6,895.00	-
Cis-syn Thymine Dimer Phosphoramidite	11-1330-95	50 µmol	4,027.00	3,866.00	3,463.00	-
Coumarin Azide	50-2004-90	100 µmol	503.00	483.00	433.00	-
Coumarin Azide	50-2004-92	25 µmol	165.00	158.00	142.00	-
C-TOM-CE Phosphoramidite - Expedite	10-3014-02E	0.25 g	83.00	79.70	71.40	-
Cy3 Phosphoramidite, Cyclone	10-5913-95C	50 µmol	295.00	283.00	254.00	-
Cy3 Phosphoramidite, Expedite	10-5913-90E	100 µmol	539.00	517.00	464.00	-
Cy3 Phosphoramidite, Expedite	10-5913-95E	50 µmol	295.00	283.00	254.00	-

Description	Reference	Scale/Quantity	Price List 2025			
			EUR	CHF	GBP	USD
Cy3.5* Phosphoramidite	10-5914-02	0.25 g	1,433.00	1,376.00	1,232.00	-
Cy3.5* Phosphoramidite	10-5914-90	100 µmol	619.00	594.00	532.00	-
Cy3.5* Phosphoramidite	10-5914-95	50 µmol	341.00	327.00	293.00	-
Cy3* Phosphoramidite	10-5913-02	0.25 g	1,330.00	1,277.00	1,144.00	-
Cy3* Phosphoramidite	10-5913-90	100 µmol	539.00	517.00	464.00	-
Cy3* Phosphoramidite	10-5913-95	50 µmol	295.00	283.00	254.00	-
Cy5 Phosphoramidite	10-5915-90E	100 µmol	539.00	517.00	464.00	-
Cy5 Phosphoramidite, Pharmacia	10-5915-95P	50 µmol	295.00	283.00	254.00	-
Cy5.5* Phosphoramidite	10-5916-02	0.25 g	1,433.00	1,376.00	1,232.00	-
Cy5.5* Phosphoramidite	10-5916-90	100 µmol	697.00	669.00	599.00	-
Cy5.5* Phosphoramidite	10-5916-90	100 µmol	697.00	669.00	599.00	-
Cy5.5* Phosphoramidite	10-5916-95	50 µmol	379.00	364.00	326.00	-
Cy5* Phosphoramidite	10-5915-02	0.25 g	1,330.00	1,277.00	1,144.00	-
Cy5* Phosphoramidite	10-5915-90	100 µmol	539.00	517.00	464.00	-
Cy5* Phosphoramidite	10-5915-95	50 µmol	295.00	283.00	254.00	-
Cyanine 3 CPG	20-5913-01	0.1 g	259.00	249.00	223.00	-
Cyanine 3 CPG	20-5913-10	1 g	2,236.00	2,147.00	1,923.00	-
Cyanine 3 CPG	20-5913-42	4 x 0.2 µmol	129.00	124.00	111.00	-
Cyanine 3 CPG - Expedite	20-5913-41E	4 x 1.0 µmol	446.00	428.00	384.00	-
Cyanine 3 CPG - Mermade	20-5913-41M	4 x 1 µmol	446.00	428.00	384.00	-
Cyanine 3 Phosphoramidite - Mermade	10-5913-90M	100 µmol	539.00	517.00	464.00	-
Cyanine 3 Phosphoramidite, M.	10-5913-02M	0.25 g	1,330.00	1,277.00	1,144.00	-
Cyanine 3.5 Phosphoramidite, M.	10-5914-02M	0.25 g	1,433.00	1,376.00	1,232.00	-
Cyanine 5 CPG - Expedite	20-5915-42E	4 x 0.2 µmol	129.00	124.00	111.00	-
Cyanine 5 Phosphoramidite	10-5915-02E	0.25 g	1,330.00	1,277.00	1,144.00	-
Cyanine 5 Phosphoramidite, Expedite	10-5915-95E	50 µmol	295.00	283.00	254.00	-
Cyanine 5 Phosphoramidite, Merm.	10-5915-02M	0.25 g	1,330.00	1,277.00	1,144.00	-
Cyanine 5.5 Phosphoramidite	10-5916-02E	0.25 g	1,433.00	1,376.00	1,232.00	-
Cyanine 5.5 Phosphoramidite, Mer.	10-5916-02M	0.25 g	1,433.00	1,376.00	1,232.00	-
dA-5'-CE Phosphoramidite	10-0001-02	0.25 g	85.20	81.80	73.30	-
dA-5'-CE Phosphoramidite	10-0001-05	0.5 g	164.00	157.00	141.00	-
dA-5'-CE Phosphoramidite	10-0001-10	1 g	316.00	303.00	272.00	-
dA-5'-CE Phosphoramidite, Expedite	10-0001-05E	0.5 g	164.00	157.00	141.00	-
dA-5'-CE Phosphoramidite, Expedite	10-0001-10E	1 g	316.00	303.00	272.00	-
dA-5'-CPG	20-0002-01	0.1 g	71.90	69.00	61.80	-
dA-5'-CPG	20-0002-10	1 g	633.00	608.00	544.00	-
dA-5'-CPG	20-0012-13	1 x 10 µmol	324.00	311.00	279.00	-
dA-5'-CPG	20-0012-14	1 x 15 µmol	431.00	414.00	371.00	-
dA-5'-CPG	20-0012-41	4 x 1 µmol	144.00	138.00	124.00	-
dA-5'-CPG	20-0012-42	4 x 0.2 µmol	108.00	104.00	92.90	-
dA-5'-lcaa-CPG 500 -	20-0012-42E	0.2 µmol	108.00	104.00	92.90	-
dA-5'-lcaa-CPG 500, column - 1 µmol	20-0012-41E	1 µmol	144.00	138.00	124.00	-
Dabcyl-dT	10-1058-02	0.25 g	971.00	932.00	835.00	-
Dabcyl-dT	10-1058-90	100 µmol	467.00	448.00	402.00	-
Dabcyl-dT	10-1058-95	50 µmol	259.00	249.00	223.00	-
Dabcyl-dT - Mermade	10-1058-02M	0.25 g	971.00	932.00	835.00	-
dA-CE Phosphoramidite, Expedite	10-1000-C5E	0.5 g	9.29	8.92	7.99	-
dA-Me Phosphoramidite	10-1100-02	0.25 g	395.00	379.00	340.00	-
dA-Me Phosphoramidite	10-1100-05	0.5 g	791.00	759.00	680.00	-
dA-Me Phosphoramidite	10-1100-05E	0.5 g	791.00	759.00	680.00	-
dA-Q-CPG 500	21-2000-01	0.1 g	33.20	31.90	28.60	-
dA-Q-CPG 500	21-2000-10	1 g	277.00	266.00	238.00	-
dA-Q-CPG 500	21-2100-41	4 x 1 µmol	66.40	63.70	57.10	-
dA-Q-CPG 500	21-2100-42	4 x 0.2 µmol	44.30	42.50	38.10	-
dA-Q-CPG 500	21-2200-41	4 x 1 µmol	66.40	63.70	57.10	-
dA-Q-CPG 500	21-2200-42	4 x 0.2 µmol	44.30	42.50	38.10	-
dA-Q-CPG 500,	21-2000-02	0.25 g	66.40	63.70	57.10	-
DBCO TEG	10-1941-02		1,309.00	1,257.00	1,126.00	-
dbco nhs ester	50-1941-23		86.30	82.80	74.20	-
DBCO-dT-CE Phosphoramidite	10-1539-90	100 µMol	733.00	704.00	630.00	-
DBCO-dT-CE phosphoramidite	10-1539-95	50 µmol	388.00	372.00	334.00	-
DBCO-dT-CE Phosphoramidite, Expedite	10-1539-95E	50 µmol	388.00	372.00	334.00	-
DBCO-PEG4-phosphoramidite	10-1941-90	100 µmol	374.00	359.00	322.00	-
DBCO-Serinol Phosphoramidite	10-1998-02	0.25 g	1,366.00	1,311.00	1,175.00	-
DBCO-Serinol Phosphoramidite	10-1998-02E		1,366.00	1,311.00	1,175.00	-
DBCO-Serinol Phosphoramidite	10-1998-95	50 µmol	259.00	249.00	223.00	-
DBCO-Serinol Phosphoramidite	10-1998-95E		259.00	249.00	223.00	-
DBCO-Serinol Phosphoramidite, Mermade	10-1998-02M	0.25 g	1,366.00	1,311.00	1,175.00	-
DBCO-sulfh-NHS Ester	50-1941-24	52 mg	431.00	414.00	371.00	-
dC-5'-CE Phosphoramidite	10-0101-02	0.25 g	85.20	81.80	73.30	-
dC-5'-CE Phosphoramidite	10-0101-05	0.5 g	164.00	157.00	141.00	-
dC-5'-CE Phosphoramidite	10-0101-10	1 g	316.00	303.00	272.00	-
dC-5'-CE Phosphoramidite, Expedite	10-0101-02E	0.25 g	85.20	81.80	73.30	-
dC-5'-CE Phosphoramidite, Expedite	10-0101-05E	0.5 g	164.00	157.00	141.00	-
dC-5'-CE Phosphoramidite, Expedite	10-0101-10E	1 g	316.00	303.00	272.00	-
dC-5'-CPG	20-0102-01	0.1 g	71.90	69.00	61.80	-
dC-5'-CPG	20-0102-10	1 g	633.00	608.00	544.00	-
dC-5'-CPG	20-0112-13	1 x 10 µmol	324.00	311.00	279.00	-
dC-5'-CPG	20-0112-14	1 x 15 µmol	431.00	414.00	371.00	-
dC-5'-CPG	20-0112-41	4 x 1 µmol	144.00	138.00	124.00	-
dC-5'-CPG	20-0112-42	4 x 0.2 µmol	108.00	104.00	92.90	-
dC-5'-lcaa-CPG 500	20-0112-42C	0.2 µmol	108.00	104.00	92.90	-
dC-5'-lcaa-CPG 500, column - 1 µmol	20-0112-41C	1 µmol	144.00	138.00	124.00	-
dC-Q-CPG 500	21-2010-01	0.1 g	33.20	31.90	28.60	-

Description	Reference	Scale/Quantity	Price List 2025			
			EUR	CHF	GBP	USD
dC-Q-CPG 500	21-2010-10	1 g	277.00	266.00	238.00	-
dC-Q-CPG 500	21-2110-41	4 x 1 µmol	66.40	63.70	57.10	-
dC-Q-CPG 500	21-2110-42	4 x 0.2 µmol	44.30	42.50	38.10	-
dC-Q-CPG 500	21-2210-41	4 x 1 µmol	66.40	63.70	57.10	-
dC-Q-CPG 500	21-2210-42	4 x 0.2 µmol	44.30	42.50	38.10	-
dC-Q-CPG 500,	21-2010-02	0.25 g	66.40	63.70	57.10	-
DEACM Caged-dG-CE Phosph.	10-1533-90E	100 µmol	647.00	621.00	556.00	-
DEACM Caged-dG-CE Phosphoramidite	10-1533-90	100 µmol	647.00	621.00	556.00	-
def-dA-CE Phosphoramidite	10-1504-10	1 g	101.00	97.00	86.90	-
def-dA-CE Phosphoramidite, Expedite	10-1504-10E	1 g	101.00	97.00	86.90	-
DesthiobiotinTEG Phosphoramidite	10-1952-02	0.25 g	1,301.00	1,249.00	1,119.00	-
DesthiobiotinTEG Phosphoramidite	10-1952-90	100 µmol	532.00	511.00	458.00	-
DesthiobiotinTEG Phosphoramidite	10-1952-95	50 µmol	280.00	269.00	241.00	-
DesthiobiotinTEG-CPG	20-2952-01	0.1 g	216.00	207.00	186.00	-
DesthiobiotinTEG-CPG	20-2952-10	1 g	1,726.00	1,657.00	1,484.00	-
DesthiobiotinTEG-CPG	20-2952-13	1 x 10 µmol	525.00	504.00	452.00	-
DesthiobiotinTEG-CPG	20-2952-14	1 x 15 µmol	791.00	759.00	680.00	-
DesthiobiotinTEG-CPG	20-2952-41	4 x 1 µmol	352.00	338.00	303.00	-
DesthiobiotinTEG-CPG	20-2952-42	4 x 0.2 µmol	216.00	207.00	186.00	-
dG-5'-CPG	20-0202-01	0.1 g	73.90	73.90	66.50	-
dG-5'-CPG	20-0202-10	1 g	591.00	591.00	532.00	-
dG-5'-CPG	20-0212-13	1 x 10 µmol	332.00	332.00	299.00	-
dG-5'-CPG	20-0212-14	1 x 15 µmol	443.00	443.00	399.00	-
dG-5'-CPG	20-0212-41	4 x 1 µmol	148.00	148.00	133.00	-
dG-5'-CPG	20-0212-42	4 x 0.2 µmol	111.00	111.00	100.00	-
dG-CE Phosphoramidite, Expedite	10-1020-C5E	0.5 g	9.29	8.92	7.99	-
dG-Me Phosphoramidite	10-1120-02	0.25 g	395.00	379.00	340.00	-
dG-Me Phosphoramidite	10-1120-05	0.5 g	791.00	759.00	680.00	-
dG-Me Phosphoramidite	10-1120-05E	0.5 g	791.00	759.00	680.00	-
dI-CE Phosphoramidite	10-1040-02	0.25 g	98.30	94.40	84.50	-
dI-CE Phosphoramidite	10-1040-90	100 µmol	42.20	40.50	36.30	-
dI-CE Phosphoramidite, Expedite - 100 µmol	10-1040-90E	100 µmol	42.20	40.50	36.30	-
dI-CPG 1000	20-2041-01	0.1 g	43.10	41.40	37.10	-
dI-CPG 1000	20-2191-41	4 x 1 µmol	173.00	166.00	149.00	-
dI-CPG 500	20-2040-01	0.1 g	57.50	55.20	49.50	-
dI-CPG 500	20-2190-41	4 x 1 µmol	180.00	173.00	155.00	-
dI-CPG 500	20-2190-42	4 x 0.2 µmol	108.00	104.00	92.90	-
dI-lcaa-CPG 500 Å, column - 0.2 µmol	20-2190-42E	0.2 µmol	108.00	104.00	92.90	-
dI-lcaa-CPG 500 Å, column - 1 µmol	20-2190-41E	1 µmol	180.00	173.00	155.00	-
Dithiol Serinol Phosphoramidite	10-1991-02	0.25 g	841.00	807.00	723.00	-
Dithiol Serinol Phosphoramidite	10-1991-90	100 µmol	309.00	297.00	266.00	-
dK-CE Phosphoramidite	10-1048-02	0.25 g	985.00	946.00	847.00	-
dK-CE Phosphoramidite	10-1048-90	100 µmol	338.00	324.00	291.00	-
dmf-5-Me-isodC-CE Phosphoramidite	10-1065-02	0.25 g	498.00	478.00	428.00	-
dmf-5-Me-isodC-CE Phosphoramidite	10-1065-90	100 µmol	159.00	153.00	137.00	-
dmf-5-Me-isodC-CE Phosphoramidite, Expedite	10-1065-02E	0.25 g	498.00	478.00	428.00	-
dmf-dG PS	26-2629-62	200 x 200 nmol	653.00	627.00	562.00	-
dmf-dG PS	26-2629-65	200 x 40 nmol	653.00	627.00	562.00	-
dmf-dG-5'-CE Phosphoramidite	10-9201-02	0.25 g	85.20	81.80	73.30	-
dmf-dG-5'-CE Phosphoramidite	10-9201-05	0.5 g	164.00	157.00	141.00	-
dmf-dG-5'-CE Phosphoramidite	10-9201-10	1 g	316.00	303.00	272.00	-
dmf-dG-5'-CE Phosphoramidite, Exp.	10-9201-10E	1 g	316.00	303.00	272.00	-
dmf-dG-5'-CE Phosphoramidite, Expedite	10-9201-05E	0.5 g	164.00	157.00	141.00	-
dmf-dG-5'-CPG	20-9202-10	1 g	633.00	608.00	544.00	-
dmf-dG-CE Phosphoramidite	10-1029-02	0.25 g	4.96	4.76	4.27	-
dmf-dG-CE Phosphoramidite	10-1029-05	0.5 g	9.29	8.92	7.99	-
dmf-dG-CE Phosphoramidite	10-1029-10	1 g	15.50	14.90	13.30	-
dmf-dG-CE Phosphoramidite	10-1029-1C	1 g	15.50	14.90	13.30	-
dmf-dG-CE Phosphoramidite	10-1029-1S	10 g	155.00	149.00	133.00	-
dmf-dG-CE Phosphoramidite	10-1029-20	2 g	31.00	29.80	26.70	-
dmf-dG-CE Phosphoramidite	10-1029-50	5 g	77.40	74.30	66.60	-
dmf-dG-CE Phosphoramidite	10-1029-5S	5 g	77.40	74.30	66.60	-
dmf-dG-CE Phosphoramidite	10-1029-C2	0.25 g	4.96	4.76	4.27	-
dmf-dG-CE Phosphoramidite	10-1029-C5	0.5 g	9.29	8.92	7.99	-
dmf-dG-CE Phosphoramidite, Expedite	10-1029-02E	0.25 g	4.96	4.76	4.27	-
dmf-dG-CE Phosphoramidite, Expedite	10-1029-C5E	0.5 g	9.29	8.92	7.99	-
dmf-dG-CPG	20-2029-01	0.1 g	13.30	12.80	11.40	-
dmf-dG-CPG	20-2029-02	0.25 g	17.70	17.00	15.20	-
dmf-dG-CPG	20-2029-10	1 g	66.40	63.70	57.10	-
dmf-dG-CPG	20-2129-13	1 x 10 µmol	88.50	85.00	76.10	-
dmf-dG-CPG	20-2129-41	4 x 1 µmol	31.00	29.80	26.70	-
dmf-dG-CPG	20-2129-42	4 x 0.2 µmol	31.00	29.80	26.70	-
dmf-dG-CPG	20-2129-45	4 x 40 nmol	31.00	29.80	26.70	-
dmf-dG-CPG	20-2229-14	1 x 15 µmol	133.00	128.00	114.00	-
dmf-dG-CPG	20-2229-41	4 x 1 µmol	22.10	21.20	19.00	-
dmf-dG-CPG	20-2229-42	4 x 0.2 µmol	15.50	14.90	13.30	-
dmf-dG-CPG	20-2229-45	4 x 40 nmol	15.50	14.90	13.30	-
dmf-dG-Q-CPG 500	21-2029-01	0.1 g	33.20	31.90	28.60	-
dmf-dG-Q-CPG 500	21-2029-10	1 g	277.00	266.00	238.00	-
dmf-dG-Q-CPG 500	21-2129-41	4 x 1 µmol	66.40	63.70	57.10	-
dmf-dG-Q-CPG 500	21-2129-42	4 x 0.2 µmol	44.30	42.50	38.10	-
dmf-dG-Q-CPG 500	21-2229-41	4 x 1 µmol	66.40	63.70	57.10	-
dmf-dG-Q-CPG 500	21-2229-42	4 x 0.2 µmol	44.30	42.50	38.10	-
dmf-dG-Q-CPG 500,	21-2029-02	0.25 g	66.40	63.70	57.10	-

Description	Reference	Scale/Quantity	Price List 2025			
			EUR	CHF	GBP	USD
dmf-G-LA Phosphoramidite	10-2029-05	0.5 g	108.00	104.00	92.90	-
dmf-G-LA-CE Phosphoramidite	10-2029-10	1g	216.00	207.00	186.00	-
dmf-G-LA-CE Phosphoramidite - Mermade	10-2029-05M	0.5 g	108.00	104.00	92.90	-
dmf-G-LA-CE Phosphoramidite - Mermade	10-2029-10M	1g	216.00	207.00	186.00	-
dmf-G-LA-CE Phosphoramidite, Expedite	10-2029-10E	1g	216.00	207.00	186.00	-
dmf-G-LA-CPG	20-2529-41E	4 x 1.0 µmol	93.50	89.80	80.40	-
dmf-isodG-CE Phosphoramidite	10-1078-02	0.25 g	432.00	415.00	372.00	-
dmf-isodG-CE Phosphoramidite	10-1078-02E	0.25 g	432.00	415.00	372.00	-
dmf-isodG-CE Phosphoramidite	10-1078-90	100 µmol	201.00	193.00	173.00	-
DNP-TEG Phosphoramidite	10-1985-02	0.25 g	971.00	932.00	835.00	-
DNP-TEG Phosphoramidite	10-1985-90	100 µmol	424.00	407.00	365.00	-
DNP-TEG Phosphoramidite	10-1985-95	50 µmol	237.00	228.00	204.00	-
DNP-TEG Phosphoramidite, Mermade	10-1985-90M	100 µmol	424.00	407.00	365.00	-
dP+dK-CE Phosphoramidite	10-1049-02	0.25 g	985.00	946.00	847.00	-
dP+dK-CE Phosphoramidite	10-1049-90	100 µmol	338.00	324.00	291.00	-
dP-CE Phosphoramidite	10-1047-02	0.25 g	999.00	959.00	859.00	-
dP-CE Phosphoramidite	10-1047-90	100 µmol	345.00	331.00	297.00	-
dSpacer CE Phosphoramidite	10-1914-02	0.25 g	415.00	398.00	357.00	-
dSpacer CE Phosphoramidite	10-1914-90	100 µmol	127.00	122.00	109.00	-
dSpacer CE Phosphoramidite	10-1914-90C	100 µmol	127.00	122.00	109.00	-
dSpacer CE Phosphoramidite, Expedite	10-1914-02E	0.25 g	415.00	398.00	357.00	-
dT-5'-CE Phosphoramidite	10-0301-02	0.25 g	85.20	81.80	73.30	-
dT-5'-CE Phosphoramidite	10-0301-02P	0.25 g	85.20	81.80	73.30	-
dT-5'-CE Phosphoramidite	10-0301-05	0.5 g	164.00	157.00	141.00	-
dT-5'-CE Phosphoramidite	10-0301-10	1g	316.00	303.00	272.00	-
dT-5'-CE Phosphoramidite, Expedite	10-0301-05E	0.5 g	164.00	157.00	141.00	-
dT-5'-CE Phosphoramidite, Expedite	10-0301-10E	1g	316.00	303.00	272.00	-
dT-5'-CE Phosphoramidite, Expedite - 0.25 g	10-0301-02E	0.25 g	85.20	81.80	73.30	-
dT-5'-CE Phosphoramidite, Mermade	10-0301-02M	0.25 g	85.20	81.80	73.30	-
dT-5'-CPG	20-0302-01	0.1 g	71.90	69.00	61.80	-
dT-5'-CPG	20-0302-10	1g	633.00	608.00	544.00	-
dT-5'-CPG	20-0312-13	1 x 10 µmol	324.00	311.00	279.00	-
dT-5'-CPG	20-0312-14	1 x 15 µmol	431.00	414.00	371.00	-
dT-5'-CPG	20-0312-41	4 x 1 µmol	144.00	138.00	124.00	-
dT-5'-CPG	20-0312-42	4 x 0.2 µmol	108.00	104.00	92.90	-
dT-5'-lcaa-CPG 500	20-0312-42C	0.2 µmol	108.00	104.00	92.90	-
dT-5'-lcaa-CPG 500, column	20-0312-41E	4 x 1 µmol	144.00	138.00	124.00	-
dT-Me Phosphoramidite	10-1130-02	0.25 g	395.00	379.00	340.00	-
dT-Me Phosphoramidite	10-1130-05	0.5 g	784.00	753.00	674.00	-
dT-Me Phosphoramidite	10-1130-05E	0.5 g	784.00	753.00	674.00	-
dT-Q-CPG 500	21-2030-01	0.1 g	33.20	31.90	28.60	-
dT-Q-CPG 500	21-2030-10	1g	277.00	266.00	238.00	-
dT-Q-CPG 500	21-2130-41	4 x 1 µmol	66.40	63.70	57.10	-
dT-Q-CPG 500	21-2130-42	4 x 0.2 µmol	44.30	42.50	38.10	-
dT-Q-CPG 500	21-2230-41	4 x 1 µmol	66.40	63.70	57.10	-
dT-Q-CPG 500,	21-2030-02	0.25 g	66.40	63.70	57.10	-
dU-CE Phosphoramidite	10-1050-02	0.25 g	88.50	85.00	76.10	-
dU-CE Phosphoramidite	10-1050-02P	0.25 g	88.50	85.00	76.10	-
dU-CE Phosphoramidite	10-1050-90	100 µmol	31.00	29.80	26.70	-
dU-CE Phosphoramidite Exp	10-1050-90E	100 µmol	31.00	29.80	26.70	-
dU-CE Phosphoramidite, Expedite - 0.25 g	10-1050-02E	0.25 g	88.50	85.00	76.10	-
dU-CPG 1000	20-2051-01	0.1 g	71.90	69.00	61.80	-
dU-CPG 1000	20-2151-41	4 x 1 µmol	288.00	276.00	248.00	-
dU-CPG 1000	20-2151-42	4 x 0.2 µmol	173.00	166.00	149.00	-
dU-CPG 500	20-2050-01	0.1 g	57.50	55.20	49.50	-
dU-CPG 500	20-2150-41	4 x 1 µmol	173.00	166.00	149.00	-
dU-CPG 500	20-2150-42	4 x 0.2 µmol	104.00	99.80	89.40	-
dX-CE Phosphoramidite	10-1537-02	0.25 g	1,424.00	1,367.00	1,225.00	-
dX-CE Phosphoramidite	10-1537-90	100 µmol	618.00	593.00	531.00	-
dX-CE Phosphoramidite	10-1537-95	50 µmol	324.00	311.00	279.00	-
Eclipse® Quencher Phosphoramidite, Expedite	10-5925-02E	0.25 g	2,876.00	2,761.00	2,473.00	-
EDTA-C2-dT-CE Phosphoramidite	10-1059-02	0.25 g	1,402.00	1,346.00	1,206.00	-
EDTA-C2-dT-CE Phosphoramidite	10-1059-90	100 µmol	712.00	684.00	612.00	-
EDTA-C2-dT-CE Phosphoramidite	10-1059-95	50 µmol	360.00	346.00	310.00	-
EDTA-C2-dT-CE Phosphoramidite, Expedite - 50 µmol	10-1059-95E	50 µmol	360.00	346.00	310.00	-
Empty MerMade Columns (200 nmol and 1 µmol)	20-0050-02	pk / 48	177.00	170.00	152.00	-
Epoch Eclipse QuencherCPG0.2µm(1pk/4)Mer	20-5925-42M	4 x 0.2 µmol	232.00	223.00	200.00	-
Epoch Eclipse QuencherCPG1µm(1pk/4)Mer	20-5925-41M	4 x 1 µmol	465.00	446.00	400.00	-
Epoch Eclipse™ Quencher CPG	20-5925-01	0.1 g	431.00	414.00	371.00	-
Epoch Eclipse™ Quencher CPG	20-5925-10	1g	3,962.00	3,804.00	3,407.00	-
Epoch Eclipse™ Quencher CPG	20-5925-13	1 x 10 µmol	1,431.00	1,374.00	1,231.00	-
Epoch Eclipse™ Quencher CPG	20-5925-14	1 x 15 µmol	2,150.00	2,064.00	1,849.00	-
Epoch Eclipse™ Quencher CPG	20-5925-41	4 x 1 µmol	465.00	446.00	400.00	-
Epoch Eclipse™ Quencher CPG	20-5925-42	4 x 0.2 µmol	232.00	223.00	200.00	-
Epoch Eclipse™ Quencher Phosphoramidite	10-5925-02	0.25 g	2,876.00	2,761.00	2,473.00	-
Epoch Eclipse™ Quencher Phosphoramidite	10-5925-90	100 µmol	1,122.00	1,077.00	965.00	-
Epoch Eclipse™ Quencher Phosphoramidite	10-5925-95	50 µmol	561.00	539.00	482.00	-
Epoch Redmond Red™ CPG	20-5920-01	0.1 g	324.00	311.00	279.00	-
Epoch Redmond Red™ CPG	20-5920-10	1g	3,063.00	2,940.00	2,634.00	-
Epoch Redmond Red™ CPG	20-5920-13	1 x 10 µmol	1,265.00	1,214.00	1,088.00	-
Epoch Redmond Red™ CPG	20-5920-14	1 x 15 µmol	1,898.00	1,822.00	1,632.00	-
Epoch Redmond Red™ CPG	20-5920-41	4 x 1 µmol	575.00	552.00	495.00	-
Epoch Redmond Red™ CPG	20-5920-42	4 x 0.2 µmol	252.00	242.00	217.00	-
Epoch Redmond Red™ Phosphoramidite	10-5920-02	0.25 g	2,696.00	2,588.00	2,319.00	-

Description	Reference	Scale/Quantity	Price List 2025			
			EUR	CHF	GBP	USD
Epoch Redmond Red™ Phosphoramidite	10-5920-90	100 µmol	1,050.00	1,008.00	903.00	-
Epoch Redmond Red™ Phosphoramidite	10-5920-95	50 µmol	539.00	517.00	464.00	-
Epoch Yakima Yellow™ CPG	20-5921-01	0.1 g	324.00	311.00	279.00	-
Epoch Yakima Yellow™ CPG	20-5921-10	1 g	3,063.00	2,940.00	2,634.00	-
Epoch Yakima Yellow™ CPG	20-5921-13	1 x 10 µmol	1,208.00	1,160.00	1,039.00	-
Epoch Yakima Yellow™ CPG	20-5921-14	1 x 15 µmol	1,798.00	1,726.00	1,546.00	-
Epoch Yakima Yellow™ CPG	20-5921-41	4 x 1 µmol	575.00	552.00	495.00	-
Epoch Yakima Yellow™ CPG	20-5921-42	4 x 0.2 µmol	252.00	242.00	217.00	-
Epoch Yakima Yellow™ Phosphoramidite	10-5921-02	0.25 g	1,007.00	967.00	866.00	-
Epoch Yakima Yellow™ Phosphoramidite	10-5921-90	100 µmol	575.00	552.00	495.00	-
Epoch Yakima Yellow™ Phosphoramidite	10-5921-95	50 µmol	288.00	276.00	248.00	-
Etheno-dA-CE Phosphoramidite	10-1006-02	0.25 g	475.00	456.00	409.00	-
Etheno-dA-CE Phosphoramidite	10-1006-90	100 µmol	180.00	173.00	155.00	-
Etheno-dA-CE Phosphoramidite, Expedite	10-1006-02E	0.25 g	475.00	456.00	409.00	-
Ferrocene-dT-CE Phosphoramidite	10-1576-02	0.25g	1,416.00	1,359.00	1,218.00	-
Ferrocene-dT-CE Phosphoramidite	10-1576-90	100 µmol	647.00	621.00	556.00	-
Fluorescein CPG	20-2963-42E	0.2 µmol	173.00	166.00	149.00	-
Fluorescein CPG, Expedite - 4 x 1 µmol	20-2964-41E	4 x 1 µmol	288.00	276.00	248.00	-
Fluorescein DT Phosphoramidite	10-1056-02	0.25 g	725.00	696.00	624.00	-
Fluorescein NHS ester	50-5901-22	2.4 mg	115.00	110.00	98.90	-
Fluorescein Phosphoramidite	10-1963-02	0.25 g	856.00	822.00	736.00	-
Fluorescein Phosphoramidite	10-1963-90	100 µmol	424.00	407.00	365.00	-
Fluorescein Phosphoramidite	10-1963-95	50 µmol	237.00	228.00	204.00	-
Fluorescein Phosphoramidite, Exp.	10-1963-02E	0.25 g	856.00	822.00	736.00	-
Fluorescein Phosphoramidite, Mermade	10-1963-02M	0.25 g	856.00	822.00	736.00	-
Fluorescein-dT Phosphoramidite	10-1056-90	100 µmol	442.00	424.00	380.00	-
Fluorescein-dT Phosphoramidite	10-1056-95	50 µmol	232.00	223.00	200.00	-
Fluorescein-dT Phosphoramidite, Expedite	10-1056-02E	0.25 g	725.00	696.00	624.00	-
Fluorescein-dT Phosphoramidite, Expedite	10-1056-95E	50 µmol	232.00	223.00	200.00	-
Fluorescein-dT Phosphoramidite, Mermade	10-1056-02M	0.25 g	725.00	696.00	624.00	-
Fmoc Amino-Modifier C6 dT	10-1536-02	0.25 g	590.00	566.00	507.00	-
Fmoc Amino-Modifier C6 dT	10-1536-90	100 µmol	288.00	276.00	248.00	-
Formyl dC	10-1564-02		5,076.00	4,873.00	4,365.00	-
G-2'-MOE-Phosphoramidite	10-3220-10	1.0 g	108.00	104.00	92.90	-
G-2'-MOE-Phosphoramidite - Mermade	10-3220-10M	1.0 g	108.00	104.00	92.90	-
GalNAc C3 CPG	20-2974-01	0.1 g	57.50	55.20	49.50	-
GalNAc C3 CPG	20-2974-10	1 g	460.00	442.00	396.00	-
GalNAc C3 CPG	20-2974-41E	4 x 1 µmol	144.00	138.00	124.00	-
GAT Trimer Phosphoramidite	13-1203-90	100 µmol	1,294.00	1,242.00	1,113.00	-
GCA Trimer Phosphoramidite	13-1210-90	100 µmol	1,294.00	1,242.00	1,113.00	-
GCG Trimer Phosphoramidite	13-1212-90	100 µmol	1,294.00	1,242.00	1,113.00	-
Glen Gel-Pak™ 0.2 Desalting Column	61-5002-05	Pk / 5	71.90	69.00	61.80	-
Glen Gel-Pak™ 1.0 Desalting Column	61-5010-05	Pk / 5	71.90	69.00	61.80	-
Glen UnySupport™ PS	26-5040-10	1 g	288.00	276.00	248.00	-
Glen-Pak™ DNA 3mg 384-Well Plate	60-5500-01	Pk / 1	597.00	573.00	513.00	-
Glen-Pak™ DNA purification cartridge	60-5100-10	Pk / 10	70.80	68.00	60.90	-
Glen-Pak™ DNA purification cartridge	60-5100-96	Pk / 96	448.00	430.00	385.00	-
Glen-Pak™ DNA purification cartridge (for use with disposable syringes)	60-5200-10	Pk / 10	122.00	117.00	105.00	-
Glycerol CPG column (1 µmole)Expedite	20-2902-41E		180.00	173.00	155.00	-
GTA Trimer Phosphoramidite	13-1230-90	100 µmol	1,294.00	1,242.00	1,113.00	-
High Load Glen UnySupport	25-5040-10	1 g	127.00	122.00	109.00	-
I-CE Phosphoramidite	10-3040-02	0.25 g	942.00	904.00	810.00	-
iPr-Pac-dG-CE Phosphoramidite	10-1621-02	0.25 g	26.60	25.50	22.90	-
iPr-Pac-dG-CE Phosphoramidite	10-1621-05	0.5 g	53.10	51.00	45.70	-
iPr-Pac-dG-CE Phosphoramidite	10-1621-10	1 g	106.00	102.00	91.20	-
iPr-Pac-dG-CE Phosphoramidite - Mermade	10-1621-05M	0.5 g	53.10	51.00	45.70	-
iPr-Pac-dG-CE Phosphoramidite, Expedite	10-1621-02E	0.25 g	26.60	25.50	22.90	-
iPr-Pac-dG-CE Phosphoramidite, Expedite	10-1621-05E	0.5 g	53.10	51.00	45.70	-
iPr-Pac-dG-CPG	20-2621-01	0.1 g	36.00	34.60	31.00	-
iPr-Pac-dG-CPG	20-2621-02	0.25 g	71.90	69.00	61.80	-
iPr-Pac-dG-CPG	20-2621-10	1 g	288.00	276.00	248.00	-
iPr-Pac-dG-CPG	20-2721-13	1 x 10 µmol	158.00	152.00	136.00	-
iPr-Pac-dG-CPG	20-2721-41	4 x 1 µmol	86.30	82.80	74.20	-
iPr-Pac-dG-CPG	20-2721-42	4 x 0.2 µmol	57.50	55.20	49.50	-
iPr-Pac-dG-CPG	20-2721-45	4 x 40 nmol	57.50	55.20	49.50	-
iPr-Pac-dG-CPG	20-2821-14	1 x 15 µmol	237.00	228.00	204.00	-
iPr-Pac-dG-CPG	20-2821-41	4 x 1 µmol	86.30	82.80	74.20	-
iPr-Pac-dG-CPG	20-2821-42	4 x 0.2 µmol	57.50	55.20	49.50	-
iPr-Pac-dG-CPG	20-2821-45	4 x 40 nmol	57.50	55.20	49.50	-
iPr-Pac-G-RNA-CPG	20-3321-10	1 g	511.00	491.00	439.00	-
Long Trebler Phosphoramidite	10-1925-02	0.25 g	431.00	414.00	371.00	-
Long Trebler Phosphoramidite	10-1925-90	100 µmol	288.00	276.00	248.00	-
Long Trebler Phosphoramidite, Expedite	10-1925-02E	0.25 g	431.00	414.00	371.00	-
Maleimide NHS Ester (SMCC)	50-1938-23	3.3 mg	86.30	82.80	74.20	-
Methacrylate C6 Phosphoramidite	10-1891-02	0.25 g	935.00	898.00	804.00	-
Methacrylate C6 Phosphoramidite	10-1891-90	100 µmol	158.00	152.00	136.00	-
Methylene Blue II Phosphoramidite	10-5961-02	0.25grams	2,157.00	2,071.00	1,855.00	-
Methylene Blue II Phosphoramidite	10-5961-90	100 µmol	856.00	822.00	736.00	-
Methylene Blue II Phosphoramidite	10-5961-95	50 µmol	446.00	428.00	384.00	-
Methylene Blue NHS Ester	50-1960-23	5.4 mg	971.00	932.00	835.00	-
MGB Eclipse CPG	20-5927-01	0.1 g	388.00	372.00	334.00	-
MGB Eclipse CPG - 1.0gram	20-5927-10	1.0 gram	3,725.00	3,576.00	3,204.00	-
MGB Eclipse CPG, Mermade	20-5927-42M	4 x 0.2 µmol	259.00	249.00	223.00	-
N2-Amino-Modifier C6 dG	10-1529-02	0.25 g	1,834.00	1,761.00	1,577.00	-

Description	Reference	Scale/Quantity	Price List 2025			
			EUR	CHF	GBP	USD
N2-Amino-Modifier C6 dG	10-1529-90	100 µmol	733.00	704.00	630.00	-
N2-Amino-Modifier C6 dG – 50 µmol	10-1529-95	50 µmol	381.00	366.00	328.00	-
N3-Cyanoethyl-dT	10-1531-02	0.25 g	863.00	828.00	742.00	-
N3-Cyanoethyl-dT	10-1531-90	100 µmol	288.00	276.00	248.00	-
N4-Ac-N4-Et-dC-CE Phosphoramidite	10-1513-02	0.25 g	1,021.00	980.00	878.00	-
N4-Et-dC-CE Phosphoramidite	10-1068-02	0.25 g	1,136.00	1,091.00	977.00	-
N4-Et-dC-CE Phosphoramidite	10-1068-90	100 µmol	374.00	359.00	322.00	-
N4-Et-dC-CE Phosphoramidite	10-1068-95	50 µmol	201.00	193.00	173.00	-
N6-Ac-N6-Me-dA-CE Phosphoramidite	10-1503-02	0.25 g	697.00	669.00	599.00	-
N6-Ac-N6-Me-dA-CE Phosphoramidite	10-1503-90	100 µmol	230.00	221.00	198.00	-
N6-Me-dA-CE Phosphoramidite	10-1003-02	0.25 g	834.00	801.00	717.00	-
N6-Me-dA-CE Phosphoramidite	10-1003-90	100 µmol	266.00	255.00	229.00	-
N6-Me-dA-CE Phosphoramidite, Expedite	10-1003-02E	0.25 g	834.00	801.00	717.00	-
N6-Methyl-A-CE Phosphoramidite	10-3005-90	0.25 g	920.00	883.00	791.00	-
N6-Methyl-A-CE Phosphoramidite	10-3005-95	50 µmol	475.00	456.00	409.00	-
N6-Methyl-A-CE Phosphoramidite Ex	10-3005-02E	0.25 g	2,193.00	2,105.00	1,886.00	-
NHS-Carboxy-dT	10-1535-02	0.25 g	841.00	807.00	723.00	-
NHS-Carboxy-dT	10-1535-90	100 µmol	338.00	324.00	291.00	-
NHS-Carboxy-dT - Expedite	10-1535-90E	100 µmol	338.00	324.00	291.00	-
N-POM Caged-dT-CE Phosphoramidite	10-1534-02	0.25 g	1,575.00	1,512.00	1,355.00	-
N-POM Caged-dT-CE Phosphoramidite	10-1534-90	100 µmol	618.00	593.00	531.00	-
O4-Triazolyl-dU-CE Phosphoramidite	10-1051-02	0.25 g	719.00	690.00	618.00	-
O4-Triazolyl-dU-CE Phosphoramidite	10-1051-90	100 µmol	230.00	221.00	198.00	-
O6-Me-dG-CE Phosphoramidite	10-1070-02	0.25 g	482.00	463.00	415.00	-
O6-Me-dG-CE Phosphoramidite	10-1070-90	100 µmol	173.00	166.00	149.00	-
O6-Me-dG-CE Phosphoramidite	10-1070-90C	100 µmol	173.00	166.00	149.00	-
O6-Phenyl-dI-CE Phosphoramidite	10-1042-02	0.25 g	654.00	628.00	562.00	-
O6-Phenyl-dI-CE Phosphoramidite	10-1042-90	100 µmol	230.00	221.00	198.00	-
O6-Phenyl-dI-CE Phosphoramidite, Pharmacia - 100 µmol	10-1042-90P	100 µmol	230.00	221.00	198.00	-
Oligo-Affinity Support (PS)	26-4001-01	0.1 g	302.00	290.00	260.00	-
Oligo-Affinity Support (PS)	26-4001-02	0.25 g	755.00	725.00	649.00	-
Oligo-Affinity Support (PS)	26-4001-10	1 g	3,020.00	2,899.00	2,597.00	-
Oligo-Affinity Support (PS)	26-4101-41	4 x 1 µmol	503.00	483.00	433.00	-
Oligo-Affinity Support (PS), Expedite	26-4101-41E	4 x 1 µmol	503.00	483.00	433.00	-
Pac-2-Amino-dA-CE Phosphoramidite	10-1585-90	100 µmol	180.00	173.00	155.00	-
Pac-2-Amino-dA-CE Phosphoramidite	10-1585-95	50 µmol	101.00	97.00	86.90	-
Pac-A-CE Phosphoramidite - Expedite	10-3000-10E	1 g	395.00	379.00	340.00	-
Pac-A-CE Phosphoramidite Expedite	10-3000-02E	0.25 g	108.00	104.00	92.90	-
Pac-A-RNA-CPG	20-3300-10	1 g	511.00	491.00	439.00	-
Pac-dA-CE Phosphoramidite	10-1601-02	0.25 g	26.60	25.50	22.90	-
Pac-dA-CE Phosphoramidite	10-1601-05	0.5 g	53.10	51.00	45.70	-
Pac-dA-CE Phosphoramidite	10-1601-10	1 g	106.00	102.00	91.20	-
Pac-dA-CE Phosphoramidite - Mermade	10-1601-05M	0.5 g	53.10	51.00	45.70	-
Pac-dA-CE Phosphoramidite - Mermade	10-1601-10M	1 g	106.00	102.00	91.20	-
Pac-dA-CE Phosphoramidite, Expedite	10-1601-02E	0.25 g	26.60	25.50	22.90	-
Pac-dA-CE Phosphoramidite, Expedite	10-1601-05E	0.5 g	53.10	51.00	45.70	-
Pac-dA-CPG	20-2601-01	0.1 g	36.00	34.60	31.00	-
Pac-dA-CPG	20-2601-02	0.25 g	71.90	69.00	61.80	-
Pac-dA-CPG	20-2601-10	1 g	288.00	276.00	248.00	-
Pac-dA-CPG	20-2701-13	1 x 10 µmol	158.00	152.00	136.00	-
Pac-dA-CPG	20-2701-41	4 x 1 µmol	86.30	82.80	74.20	-
Pac-dA-CPG	20-2701-42	4 x 0.2 µmol	57.50	55.20	49.50	-
Pac-dA-CPG	20-2701-45	4 x 40 nmol	57.50	55.20	49.50	-
Pac-dA-CPG	20-2801-14	1 x 15 µmol	237.00	228.00	204.00	-
Pac-dA-CPG	20-2801-41	4 x 1 µmol	86.30	82.80	74.20	-
Pac-dA-CPG	20-2801-42	4 x 0.2 µmol	57.50	55.20	49.50	-
Pac-dA-CPG	20-2801-45	4 x 40 nmol	57.50	55.20	49.50	-
Palmitate Phosphoramidite	10-1978-02	0.25 g	971.00	932.00	835.00	-
Palmitate Phosphoramidite	10-1978-95E		129.00	124.00	111.00	-
Palmitate Phosphoramidite - Mermade	10-1978-02M	0.25 g	971.00	932.00	835.00	-
PC Amino-Modifier Phosphoramidite	10-4906-02	0.25 g	1,136.00	1,091.00	977.00	-
PC Amino-Modifier Phosphoramidite	10-4906-90	100 µmol	280.00	269.00	241.00	-
PC Biotin Phosphoramidite	10-4950-02	0.25 g	899.00	863.00	773.00	-
PC Biotin Phosphoramidite	10-4950-90	100 µmol	374.00	359.00	322.00	-
PC Biotin Phosphoramidite	10-4950-95	50 µmol	201.00	193.00	173.00	-
PC Linker Phosphoramidite	10-4920-02	0.25 g	1,143.00	1,097.00	983.00	-
PC Linker Phosphoramidite	10-4920-90	100 µmol	367.00	352.00	316.00	-
PC Linker Phosphoramidite, Expedite	10-4920-90E	100 µmol	367.00	352.00	316.00	-
PC Spacer Phosphoramidite	10-4913-02	0.25 g	654.00	628.00	562.00	-
PC Spacer Phosphoramidite	10-4913-90	100 µmol	216.00	207.00	186.00	-
PC Spacer Phosphoramidite - Expedite	10-4913-90E	100 µmol	216.00	207.00	186.00	-
pdC-CE Phosphoramidite	10-1014-02	0.25 g	338.00	324.00	291.00	-
pdC-CE Phosphoramidite	10-1014-05	0.5 g	676.00	649.00	581.00	-
pdC-CE Phosphoramidite	10-1014-90	100 µmol	158.00	152.00	136.00	-
pdC-CE Phosphoramidite - Mermade	10-1014-05M	0.5 g	676.00	649.00	581.00	-
pdC-CE Phosphoramidite, Expedite	10-1014-05E	0.5 g	676.00	649.00	581.00	-
pdC-CE Phosphoramidite, Mermade	10-1014-02M	0.25 g	338.00	324.00	291.00	-
pdU-CE Phosphoramidite	10-1054-02	0.25 g	288.00	276.00	248.00	-
pdU-CE Phosphoramidite	10-1054-05	0.5 g	575.00	552.00	495.00	-
pdU-CE Phosphoramidite - Mermade	10-1054-05M	0.5 g	575.00	552.00	495.00	-
pdU-CE Phosphoramidite, Expedite	10-1054-05E	0.5 g	575.00	552.00	495.00	-
pdU-CE Phosphoramidite, Mermade	10-1054-02M	0.25 g	288.00	276.00	248.00	-
Perylene-dU	10-1591-02	0.25 g	1,186.00	1,139.00	1,020.00	-
Perylene-dU	10-1591-90	100 µmol	489.00	469.00	421.00	-

Description	Reference	Scale/Quantity	Price List 2025			
			EUR	CHF	GBP	USD
Perylene-dU	10-1591-95	50 µmol	244.00	234.00	210.00	-
Perylene-dU-CE Phosphoramidite, Expedite	10-1591-90E	100 µmol	489.00	469.00	421.00	-
Protected Biotin Serinol Phosphoramidite	10-1993-02	0.25 g	971.00	932.00	835.00	-
Protected Biotin Serinol Phosphoramidite	10-1993-90	100 µmol	424.00	407.00	365.00	-
Protected Biotin Serinol Phosphoramidite	10-1993-95	50 µmol	237.00	228.00	204.00	-
Protected BiotinLC Serinol Phosphoramidite	10-1995-02	0.25 g	971.00	932.00	835.00	-
Protected BiotinLC Serinol Phosphoramidite	10-1995-90	100 µmol	525.00	504.00	452.00	-
Protected BiotinLC Serinol Phosphoramidite	10-1995-95	50 µmol	295.00	283.00	254.00	-
Protected BiotinLC Serinol Phosphoramidite, Mermade	10-1995-02M	0.25 g	971.00	932.00	835.00	-
PseudoUridine-CE Phosphoramidite	10-3055-02	0.25 g	1,790.00	1,718.00	1,539.00	-
PseudoUridine-CE Phosphoramidite	10-3055-02M	0.25 g	1,790.00	1,718.00	1,539.00	-
PseudoUridine-CE Phosphoramidite	10-3055-90	100 µmol	633.00	608.00	544.00	-
PseudoUridine-CE Phosphoramidite	10-3055-95	50 µmol	331.00	318.00	285.00	-
Psoralen Azide	50-2009-92	25 µmol	165.00	158.00	142.00	-
Psoralen C2 Phosphoramidite	10-1982-02	0.25 g	791.00	759.00	680.00	-
Psoralen C2 Phosphoramidite	10-1982-90	100 µmol	280.00	269.00	241.00	-
Psoralen C6 Phosphoramidite	10-1983-02	0.25 g	784.00	753.00	674.00	-
Psoralen C6 Phosphoramidite	10-1983-90	100 µmol	295.00	283.00	254.00	-
Puromycin, column, Expedite - 4 x 1 µmol	20-4140-41E	4 x 1 µmol	302.00	290.00	260.00	-
Puromycin-CPG	20-4040-01	0.1 g	187.00	180.00	161.00	-
Puromycin-CPG	20-4040-10	1 g	1,611.00	1,547.00	1,385.00	-
Puromycin-CPG	20-4140-13	1 x 10 µmol	539.00	517.00	464.00	-
Puromycin-CPG	20-4140-14	1 x 15 µmol	805.00	773.00	692.00	-
Puromycin-CPG	20-4140-41	4 x 1 µmol	302.00	290.00	260.00	-
Puromycin-CPG	20-4140-42	4 x 0.2 µmol	180.00	173.00	155.00	-
Pyrene-dU	10-1590-02	0.25 g	971.00	932.00	835.00	-
Pyrene-dU	10-1590-90	100 µmol	374.00	359.00	322.00	-
Pyrene-dU	10-1590-95	50 µmol	187.00	180.00	161.00	-
Pyrene-dU-CE Phosphoramidite, Expedite	10-1590-90E	100 µmol	374.00	359.00	322.00	-
Pyrrolo-C-TOM-CE Phosphoramidite, Pharmacia - 100 µmol	10-3017-90P	100 µmol	712.00	684.00	612.00	-
Pyrrolo-dC-CE Phosphoramidite	10-1017-02	0.25 g	1,136.00	1,091.00	977.00	-
Pyrrolo-dC-CE Phosphoramidite	10-1017-90	100 µmol	360.00	346.00	310.00	-
Pyrrolo-dC-CE Phosphoramidite	10-1017-95	50 µmol	194.00	186.00	167.00	-
Quasar® 570-dT Phosphoramidite	10-5953-90	100 µmol	899.00	863.00	773.00	-
Quasar® 670-dT Phosphoramidite	10-5955-02	0.25 g	1,539.00	1,477.00	1,324.00	-
Quasar® 670-dT Phosphoramidite	10-5955-90	100 µmol	949.00	911.00	816.00	-
Quasar® 670-dT Phosphoramidite	10-5955-95	50 µmol	496.00	476.00	427.00	-
Quenching buffer for Glen-Pak RNA purification	60-4120-80	1 l	177.00	170.00	152.00	-
Quenching buffer for Glen-Pak RNA purification	60-4120-82	250 mL	70.80	68.00	60.90	-
Replacement Filters-Expedite	20-0021-0F	Pk / 20	22.10	21.20	19.00	-
Replacement Frits - TWIST 10 µmol / 15 µmol	20-0040-0F	Pk / 20	26.60	25.50	22.90	-
Ribo-tC® Phosphoramidite - 0.25 g	10-3517-02	0.25 g	2,136.00	2,051.00	1,837.00	-
ROX NHS Ester - 3.2mg	50-5911-22	3.2mg	194.00	186.00	167.00	-
rSpacer TBDMS CE Phosphoramidite	10-3915-02	0.25 g	597.00	573.00	513.00	-
S-Bz-Thiol-Modifier C6-dT	10-1538-02	0.25 g	978.00	939.00	841.00	-
S-Bz-Thiol-Modifier C6-dT	10-1538-90	100 µmol	431.00	414.00	371.00	-
S-Bz-Thiol-Modifier C6-dT	10-1538-95	50 µmol	230.00	221.00	198.00	-
SIMA (HEX) Phosphoramidite	10-5905-90	100 µmol	273.00	262.00	235.00	-
SIMA (HEX) Phosphoramidite,	10-5905-02	0.25 g	618.00	593.00	531.00	-
SIMA (HEX) Phosphoramidite, Expedite	10-5905-90E	100 µmol	273.00	262.00	235.00	-
SIMA (HEX)-dT Phosphoramidite	10-5945-90	100 µmol	971.00	932.00	835.00	-
Solid Chemical Phosphorylation Reagent II - Expedite	10-1902-02E	0.25 g	288.00	276.00	248.00	-
Solid Chemical Phosphor Reagent II,	10-1902-02	0.25 g	288.00	276.00	248.00	-
SolidChemicalPhosphor Reagent II,	10-1902-90	100 µmol	86.30	82.80	74.20	-
Spacer C12 CE Phosphoramidite	10-1928-02	0.25 g	345.00	331.00	297.00	-
Spacer C12 CE Phosphoramidite	10-1928-90	100 µmol	137.00	132.00	118.00	-
Spacer Phosphoramidite 18	10-1918-02	0.25 g	265.00	254.00	228.00	-
Spacer Phosphoramidite 18	10-1918-02M	0.25 g	273.00	273.00	246.00	-
Spacer Phosphoramidite 18	10-1918-90	100 µmol	105.00	101.00	90.30	-
Spacer Phosphoramidite 18, Expedite	10-1918-02E	0.25 g	265.00	254.00	228.00	-
Spacer Phosphoramidite 9	10-1909-02	0.25 g	345.00	331.00	297.00	-
Spacer Phosphoramidite 9	10-1909-90	100 µmol	108.00	104.00	92.90	-
Spacer Phosphoramidite 9, Expedite - 0.25 g	10-1909-02E	0.25 g	345.00	331.00	297.00	-
Spacer Phosphoramidite 9, Merm.	10-1909-02M	0.25 g	345.00	331.00	297.00	-
Spacer Phosphoramidite C3	10-1913-02	0.25 g	345.00	331.00	297.00	-
Spacer Phosphoramidite C3	10-1913-90	100 µmol	108.00	104.00	92.90	-
Spacer Phosphoramidite C3, Expedite	10-1913-90E	100 µmol	108.00	104.00	92.90	-
Spacer Phosphoramidite C3, Expedite - 0.25 g	10-1913-02E	0.25 g	345.00	331.00	297.00	-
Spacer Phosphoramidite C3, Mermade	10-1913-02M	0.25 g	345.00	331.00	297.00	-
Spermine Phosphoramidite	10-1939-02	0.25 g	848.00	814.00	729.00	-
Spermine Phosphoramidite	10-1939-90	100 µmol	446.00	428.00	384.00	-
Spermine Phosphoramidite	10-1939-95	50 µmol	237.00	228.00	204.00	-
Stearyl Phosphoramidite	10-1979-02	0.25 g	259.00	249.00	223.00	-
Stearyl Phosphoramidite, Expedite	10-1979-02E	0.25 g	259.00	249.00	223.00	-
SulfoCyanine 3 NHS Ester	50-5913-22	3.8 mg	230.00	221.00	198.00	-
SulfoCyanine 5 NHS Ester	50-5915-22	3.9 mg	237.00	228.00	204.00	-
Symmetric Doubler Phosphoramidite	10-1920-02	0.25 g	360.00	346.00	310.00	-
Symmetric Doubler Phosphoramidite	10-1920-90	100 µmol	223.00	214.00	192.00	-
Symmetric Doubler Phosphoramidite, Expedite	10-1920-90E	100 µmol	223.00	214.00	192.00	-
TAMRA NHS Ester	50-5910-66	60 µl	446.00	428.00	384.00	-
TAMRA-dT	10-1057-02	0.25 g	1,402.00	1,346.00	1,206.00	-
TAMRA-dT	10-1057-90	100 µmol	712.00	684.00	612.00	-
TAMRA-dT	10-1057-95	50 µmol	360.00	346.00	310.00	-
tC®-CE Phosphoramidite	10-1517-02	250 mg	2,301.00	2,209.00	1,979.00	-

Description	Reference	Scale/Quantity	Price List 2025			
			EUR	CHF	GBP	USD
tC ^o -CE Phosphoramidite	10-1517-90	100 µmol	755.00	725.00	649.00	-
tC ^o -CE Phosphoramidite	10-1517-95	50 µmol	395.00	379.00	340.00	-
tC-CE Phosphoramidite	10-1516-90	100 µmol	705.00	677.00	606.00	-
tC-CE Phosphoramidite	10-1516-95	50 µmol	360.00	346.00	310.00	-
tC-CE Phosphoramidite - 0.25 g	10-1516-02	0.25 g	2,100.00	2,016.00	1,806.00	-
tC-CE Phosphoramidite, Expedite	10-1516-02E	0.25 g	2,100.00	2,016.00	1,806.00	-
tC-CE Phosphoramidite, Expedite	10-1516-90E	100 µmol	705.00	677.00	606.00	-
tNitro-CE Phosphoramidite	10-1518-95	50 µmol	467.00	448.00	402.00	-
TEMPO Azide	50-2007-90	100 µmoles	503.00	483.00	433.00	-
TEMPO Azide, 25 µmol	50-2007-92	Pk / 1	165.00	158.00	142.00	-
Thiazole Orange NHS Ester	50-1970-23	5.4 mg	86.30	82.80	74.20	-
Thiol-Modifier C6 S-S	10-1936-02	0.25 g	319.00	306.00	274.00	-
Thiol-Modifier C6 S-S	10-1936-90	100 µmol	133.00	128.00	114.00	-
Thiol-Modifier C6 S-S - Expedite	10-1936-90E	100 µmol	133.00	128.00	114.00	-
Thiol-Modifier C6 S-S, Cyclone	10-1936-90C	100 µmol	133.00	128.00	114.00	-
Thiol-Modifier C6 S-S, Expedite	10-1936-02E	0.25 g	319.00	306.00	274.00	-
Thiol-Modifier C6 S-S, Mermade	10-1936-02M	0.25 g	319.00	306.00	274.00	-
THPTA	50-1004-90	100 µmol	259.00	249.00	223.00	-
Thymidine Glycol CE Phosphoramidite	10-1096-02	0.25 g	1,632.00	1,567.00	1,404.00	-
Thymidine Glycol CE Phosphoramidite	10-1096-90	100 µmol	690.00	662.00	593.00	-
Thymidine Glycol CE Phosphoramidite	10-1096-95	50 µmol	360.00	346.00	310.00	-
TIPS-5-Ethynyl-dU-CE Phosphoramidite	10-1555-02	0.25 g	1,402.00	1,346.00	1,206.00	-
TIPS-5-Ethynyl-dU-CE Phosphoramidite	10-1555-90	100 µmol	532.00	511.00	458.00	-
T-LA-CE Phosphoramidite	10-2030-10	1 g	216.00	207.00	186.00	-
T-LA-CE Phosphoramidite - Mermade	10-2030-05M	0.5 g	108.00	104.00	92.90	-
T-LA-CE Phosphoramidite - Mermade	10-2030-10M		216.00	207.00	186.00	-
T-LA-CE Phosphoramidite, Expedite	10-2030-10E	1 g	216.00	207.00	186.00	-
T-LA-CPG	20-2531-41E	4 x 1.0 µmol	93.50	89.80	80.40	-
TMP-F-dU-CE Phosphoramidite	10-1016-02	0.25 g	834.00	801.00	717.00	-
TMP-F-dU-CE Phosphoramidite	10-1016-90	100 µmol	309.00	297.00	266.00	-
Treble Phosphoramidite	10-1922-02	0.25 g	662.00	636.00	569.00	-
Treble Phosphoramidite	10-1922-90	100 µmol	374.00	359.00	322.00	-
Treble Phosphoramidite, Expedite	10-1922-90E	100 µmol	374.00	359.00	322.00	-
Trimer Phosphoramidite Mix 2	13-1992-90	100 µmol	1,869.00	1,794.00	1,607.00	-
Trimer Phosphoramidite Mix 2	13-1992-95	50 µmol	956.00	918.00	822.00	-
Trimer Phosphoramidite Mix 2 - Mix of 19 Trimers	13-1992-90E	100 µmol	1,869.00	1,794.00	1,607.00	-
TTT Trimer Phosphoramidite	13-1333-90	100 µmol	1,294.00	1,242.00	1,113.00	-
U-CE Phosphoramidite - Expedite	10-3030-10E	1 g	99.60	95.60	85.70	-
UniCap Phosphoramidite	10-4410-02	0.25 g	93.50	89.80	80.40	-
UniCap Phosphoramidite	10-4410-05	0.5 g	173.00	166.00	149.00	-
UniCap Phosphoramidite	10-4410-10	1 g	345.00	331.00	297.00	-
UniCap Phosphoramidite	10-4410-20	2 g	690.00	662.00	593.00	-
UniCap Phosphoramidite, Mermade	10-4410-10M	1 g	345.00	331.00	297.00	-
UniCap Phosphoramidite, Mermade	10-4410-20M	2 g	690.00	662.00	593.00	-
Universal Support III PS	26-5010-01	0.1 g	36.00	34.60	31.00	-
Universal Support III PS	26-5010-02	0.25 g	71.90	69.00	61.80	-
Universal Support III PS	26-5110-41	4 x 1 µmol	53.10	51.00	45.70	-
Universal Support III PS	26-5110-42	4 x 0.2 µmol	35.40	34.00	30.40	-
Universal Support III PS	26-5110-52	10 x 200 nmol	97.40	93.50	83.80	-
Universal Support III PS	26-5110-55	10 x 40 nmol	97.40	93.50	83.80	-
Universal Support III PS	26-5210-41	4 x 1 µmol	53.10	51.00	45.70	-
Universal Support III PS, Expedite	26-5210-42	4 x 0.2 µmol	35.40	34.00	30.40	-
Universal-CE Phosphoramidite	10-5000-90	100 µmol	64.70	62.10	55.60	-
U-TOM-CE Phosphoramidite - Expedite	10-3034-02E	0.25 g	83.00	79.70	71.40	-
U-TOM-CE Phosphoramidite, Expedite	10-3034-10E	1 g	304.00	292.00	261.00	-
Yakima Yellow™ Phosphoramidite - Expedite	10-5921-90E	100 µmol	575.00	552.00	495.00	-
Zebularine-CE Phosphoramidite	10-3011-02	0.25 g	1,201.00	1,153.00	1,033.00	-
Zebularine-CE Phosphoramidite	10-3011-90	100 µmol	403.00	387.00	347.00	-
Zebularine-CE Phosphoramidite	10-3011-95	50 µmol	216.00	207.00	186.00	-
α-Tocopherol-TEG Phosphoramidite, Mermade	10-1977-90M	100 µmol	475.00	456.00	409.00	-
β-L-Ac-dC-CE Phosphoramidite	10-2115-02	0.25 g	108.00	104.00	92.90	-
β-L-Ac-dC-CE Phosphoramidite	10-2115-05	0.5 g	216.00	207.00	186.00	-
β-L-Ac-dC-CE Phosphoramidite	10-2115-10	1.0 g	431.00	414.00	371.00	-
β-L-dT-CE Phosphoramidite	10-2130-02	0.25 g	108.00	104.00	92.90	-
β-L-dT-CE Phosphoramidite	10-2130-05	0.5 g	216.00	207.00	186.00	-
β-L-dT-CE Phosphoramidite	10-2130-10	1.0 g	431.00	414.00	371.00	-
β-L-iPr-Pac-dG-CE Phosphoramidite	10-2121-02	0.25 g	216.00	207.00	186.00	-
β-L-iPr-Pac-dG-CE Phosphoramidite	10-2121-05	0.5 g	431.00	414.00	371.00	-
β-L-iPr-Pac-dG-CE Phosphoramidite	10-2121-10	1.0 g	863.00	828.00	742.00	-
β-L-Pac-dA-CE Phosphoramidite	10-2101-05	0.5 g	431.00	414.00	371.00	-
β-L-Pac-dA-CE Phosphoramidite	10-2101-10	1.0 g	863.00	828.00	742.00	-
Glen RNA Synthesis						
1-Methyl-PseudoUridine Phosphoram. 0.25g	10-3056-02E	0.25 g	4,034.00	3,873.00	3,469.00	-
1-Methyl-PseudoUridine Phosphoram. 0.25g	10-3056-02M	0.25 g	4,034.00	3,873.00	3,469.00	-
2'-F-A-CE Phosphoramidite, Expedite	10-3400-02E	0.25 g	108.00	104.00	92.90	-
2'-F-U-CE Phosphoramidite, Expedite	10-3430-02E	0.25 g	71.90	69.00	61.80	-
2'-OME-ibu-G-CE Phosphoramidite, Expedite	10-3120-02E	0.25 g	30.40	29.20	26.10	-
2,6-Diaminopurine-TOM-CE Phosphoramidite	10-3085-02	0.25 g	1,402.00	1,346.00	1,206.00	-
2-Aminopurine-TBDMs-CE Phosphoramidite	10-3070-02	0.25 g	1,632.00	1,567.00	1,404.00	-
2-Aminopurine-TBDMs-CE Phosphoramidite	10-3070-90	100 µmol	633.00	608.00	544.00	-
2-Aminopurine-TBDMs-CE Phosphoramidite	10-3070-95	50 µmol	331.00	318.00	285.00	-
2'-F-Ac-C-CE Phosphoramidite	10-3415-02	0.25 g	71.90	69.00	61.80	-
2'-F-Ac-C-CE Phosphoramidite	10-3415-05	0.5 g	144.00	138.00	124.00	-
2'-F-Ac-C-CE Phosphoramidite, Expedite	10-3415-05E	0.5 g	144.00	138.00	124.00	-

EA1E1Z

Description	Reference	Scale/Quantity	Price List 2025			
			EUR	CHF	GBP	USD
2'-F-Ac-C-CE Phosphoramidite, Expedite	10-3415-02E	0.25 g	71.90	69.00	61.80	-
2'-F-I-CE Phosphoramidite - 100 µmol	10-3440-90	100 µmol	137.00	132.00	118.00	-
2'-F-U-ANA-CE Phosphoramidite - 0.25 g	10-3830-02	0.25 g	424.00	407.00	365.00	-
2'-F-U-CE Phosphoramidite	10-3430-02	0.25 g	71.90	69.00	61.80	-
2'-F-U-CE Phosphoramidite	10-3430-05	0.5 g	144.00	138.00	124.00	-
2'ome a thiophosphoramidite	10-3170-90		431.00	414.00	371.00	-
2'-OMe-2,6-Diaminopurine-CE Phosphoramidite	10-3124-02	0.25 g	1,769.00	1,698.00	1,521.00	-
2'-OMe-2,6-Diaminopurine-CE Phosphoramidite	10-3124-90	100 µmol	662.00	636.00	569.00	-
2'-OMe-2,6-Diaminopurine-CE Phosphoramidite	10-3124-95	50 µmol	345.00	331.00	297.00	-
2'-OMe-5-Br-U-CE Phosphoramidite	10-3190-02	0.25 g	1,301.00	1,249.00	1,119.00	-
2'-OMe-5-Br-U-CE Phosphoramidite	10-3190-90	100 µmol	475.00	456.00	409.00	-
2'-OMe-5-Me-C-CE Phosphoramidite	10-3160-02	0.25 g	971.00	932.00	835.00	-
2'-OMe-5-Me-C-CE Phosphoramidite	10-3160-90	100 µmol	345.00	331.00	297.00	-
2'-OMe-5-Me-U-CE Phosphoramidite	10-3131-02	0.25 g	518.00	497.00	445.00	-
2'-OMe-5-Me-U-CE Phosphoramidite	10-3131-90	100 µmol	216.00	207.00	186.00	-
2'-OMe-5-Me-U-CE Phosphoramidite, Mermade	10-3131-02M	0.25 g	518.00	497.00	445.00	-
2'-OMe-Ac-C-CE Phosphoramidite	10-3115-02	0.25 g	30.40	29.20	26.10	-
2'-OMe-Ac-C-CE Phosphoramidite	10-3115-05	0.5 g	60.80	58.40	52.30	-
2'-OMe-Ac-C-CE Phosphoramidite	10-3115-10	1 g	122.00	117.00	105.00	-
2'-OMe-Ac-C-CE Phosphoramidite	10-3115-90	100 µmol	19.90	19.10	17.10	-
2'-OMe-Ac-C-CE Phosphoramidite, Expedite	10-3115-02E	0.25 g	30.40	29.20	26.10	-
2'-OMe-Ac-C-CE Phosphoramidite, Expedite	10-3115-05E	0.5 g	60.80	58.40	52.30	-
2'-OMe-Ac-C-RNA-CPG	20-3615-01	0.1 g	57.50	55.20	49.50	-
2'-OMe-Ac-C-RNA-CPG	20-3615-02	0.25 g	137.00	132.00	118.00	-
2'-OMe-Ac-C-RNA-CPG	20-3615-10	1 g	511.00	491.00	439.00	-
2'-OMe-Ac-C-RNA-CPG	20-3715-13	1 x 10 µmol	324.00	311.00	279.00	-
2'-OMe-Ac-C-RNA-CPG	20-3715-14	1 x 15 µmol	431.00	414.00	371.00	-
2'-OMe-Ac-C-RNA-CPG	20-3715-41	4 x 1 µmol	144.00	138.00	124.00	-
2'-OMe-Ac-C-RNA-CPG	20-3715-42	4 x 0.2 µmol	108.00	104.00	92.90	-
2'-OMe-A-CE Phosphoramidite	10-3100-02E	0.25 g	30.40	29.20	26.10	-
2'-OMe-A-CE Phosphoramidite	10-3100-05	0.5 g	60.80	58.40	52.30	-
2'-OMe-A-CE Phosphoramidite	10-3100-10	1 g	122.00	117.00	105.00	-
2'-OMe-A-CE Phosphoramidite	10-3100-90	100 µmol	19.90	19.10	17.10	-
2'-OMe-A-CE Phosphoramidite Ex	10-3100-90C		19.90	19.10	17.10	-
2'-OMe-A-CE Phosphoramidite, Expedite	10-3100-02	0.25 g	30.40	29.20	26.10	-
2'-OMe-A-CE Phosphoramidite, Expedite	10-3100-05E	0.5 g	60.80	58.40	52.30	-
2'-OMe-A-RNA-CPG	20-3600-01	0.1 g	57.50	55.20	49.50	-
2'-OMe-A-RNA-CPG	20-3600-02	0.25 g	137.00	132.00	118.00	-
2'-OMe-A-RNA-CPG	20-3600-10	1 g	511.00	491.00	439.00	-
2'-OMe-A-RNA-CPG	20-3700-13	1 x 10 µmol	324.00	311.00	279.00	-
2'-OMe-A-RNA-CPG	20-3700-14	1 x 15 µmol	431.00	414.00	371.00	-
2'-OMe-A-RNA-CPG	20-3700-41	4 x 1 µmol	144.00	138.00	124.00	-
2'-OMe-A-RNA-CPG	20-3700-42	4 x 0.2 µmol	108.00	104.00	92.90	-
2'-OMe-A-RNA-CPG, Expedite	20-3700-42E	4 x 0.2 µmol	108.00	104.00	92.90	-
2'OMe-A-Thiophosphoramidite	10-3170-02	0.25 g	1,035.00	994.00	890.00	-
2'-OMe-C-RNA 500 Å - 0.2 µmol	20-3710-42E	0.2 µmol	108.00	104.00	92.90	-
2'-OMe-C-RNA-CPG	20-3610-01	0.1 g	57.50	55.20	49.50	-
2'-OMe-C-RNA-CPG	20-3610-02	0.25 g	137.00	132.00	118.00	-
2'-OMe-C-RNA-CPG	20-3610-10	1 g	511.00	491.00	439.00	-
2'-OMe-C-RNA-CPG	20-3710-13	1 x 10 µmol	324.00	311.00	279.00	-
2'-OMe-C-RNA-CPG	20-3710-14	1 x 15 µmol	431.00	414.00	371.00	-
2'-OMe-C-RNA-CPG	20-3710-41	4 x 1 µmol	144.00	138.00	124.00	-
2'-OMe-C-RNA-CPG	20-3710-42	4 x 0.2 µmol	108.00	104.00	92.90	-
2'-OMe-G-CE Phosphoramidite	10-3121-02	0.25 g	49.80	47.80	42.80	-
2'-OMe-G-CE Phosphoramidite	10-3121-05	0.5 g	99.60	95.60	85.70	-
2'-OMe-G-CE Phosphoramidite	10-3121-10	1 g	199.00	191.00	171.00	-
2'-OMe-G-CE Phosphoramidite	10-3121-90	100 µmol	24.30	23.30	20.90	-
2'-OMe-G-CE Phosphoramidite Ex	10-3121-90C		24.30	23.30	20.90	-
2'-OMe-G-CE Phosphoramidite, Expedite	10-3121-02E	0.25 g	49.80	47.80	42.80	-
2'-OMe-G-CE Phosphoramidite, Expedite	10-3121-05E	0.5 g	99.60	95.60	85.70	-
2'-OMe-G-RNA 500 Å - 0.2 µmol	20-3721-42E	0.2 µmol	108.00	104.00	92.90	-
2'-OMe-G-RNA-CPG	20-3621-01	0.1 g	57.50	55.20	49.50	-
2'-OMe-G-RNA-CPG	20-3621-02	0.25 g	137.00	132.00	118.00	-
2'-OMe-G-RNA-CPG	20-3621-10	1 g	511.00	491.00	439.00	-
2'-OMe-G-RNA-CPG	20-3721-13	1 x 10 µmol	324.00	311.00	279.00	-
2'-OMe-G-RNA-CPG	20-3721-14	1 x 15 µmol	431.00	414.00	371.00	-
2'-OMe-G-RNA-CPG	20-3721-41	4 x 1 µmol	144.00	138.00	124.00	-
2'-OMe-G-RNA-CPG	20-3721-42	4 x 0.2 µmol	108.00	104.00	92.90	-
2'-OMe-G-Thiophosphoramidite	10-3172-02	0.25 g	1,035.00	994.00	890.00	-
2'-oMe-ibu-G-CE Phosphoramidite	10-3120-02	0.25 g	30.40	29.20	26.10	-
2'-OMe-I-CE Phosphoramidite	10-3140-02	0.25 g	604.00	580.00	519.00	-
2'-OMe-I-CE Phosphoramidite	10-3140-90	100 µmol	237.00	228.00	204.00	-
2'-OMe-iPr-Pac-G-CE Phosphoramidite	10-3621-02	0.25 g	89.90	86.30	77.30	-
2'-OMe-iPr-Pac-G-CE Phosphoramidite	10-3621-05	0.5 g	180.00	173.00	155.00	-
2'-OMe-iPr-Pac-G-CE Phosphoramidite	10-3621-10	1 g	360.00	346.00	310.00	-
2'-OMe-Pac-A-CE Phosphoramidite	10-3601-02	0.25 g	89.90	86.30	77.30	-
2'-OMe-Pac-A-CE Phosphoramidite	10-3601-05	0.5 g	180.00	173.00	155.00	-
2'-OMe-U-CE Phosphoramidite	10-3130-02	0.25 g	30.40	29.20	26.10	-
2'-OMe-U-CE Phosphoramidite	10-3130-05	0.5 g	60.80	58.40	52.30	-
2'-OMe-U-CE Phosphoramidite	10-3130-10	1 g	122.00	117.00	105.00	-
2'-OMe-U-CE Phosphoramidite	10-3130-90	100 µmol	19.90	19.10	17.10	-
2'-OMe-U-CE Phosphoramidite, Expedite	10-3130-02E	0.25 g	30.40	29.20	26.10	-
2'-OMe-U-CE Phosphoramidite, Expedite - 0.5 g	10-3130-05E	0.5 g	60.80	58.40	52.30	-
2'-OMe-U-RNA 500 Å - 0.2 µmol	20-3730-42E	0.2 µmol	108.00	104.00	92.90	-

Description	Reference	Scale/Quantity	Price List 2025			
			EUR	CHF	GBP	USD
2'-OMe-U-RNA-CPG	20-3630-02	0.25 g	137.00	132.00	118.00	-
2'-OMe-U-RNA-CPG	20-3630-10	1g	511.00	491.00	439.00	-
2'-OMe-U-RNA-CPG	20-3730-13	1 x 10 µmol	324.00	311.00	279.00	-
2'-OMe-U-RNA-CPG	20-3730-14	1 x 15 µmol	431.00	414.00	371.00	-
2'-OMe-U-RNA-CPG	20-3730-41	4 x 1 µmol	144.00	138.00	124.00	-
2'-OMe-U-RNA-CPG	20-3730-42	4 x 0.2 µmol	108.00	104.00	92.90	-
2'OMe-U-Thiophosphoramidite	10-3173-02	0.25 g	1,035.00	994.00	890.00	-
3'-Phosphate PS 0.1g	26-2900-01	0.1 g	108.00	104.00	92.90	-
3'-Propargyl-5-Me-dC CPG - 1 g	20-2982-10	1g	2,682.00	2,575.00	2,307.00	-
3'-PT-Amino-Modifier C6 PS 0.1 g	26-2956-01	0.1 g	180.00	173.00	155.00	-
4-Thio-U-TOM-CE Phosphoramidite	10-3052-02	0.25 g	1,402.00	1,346.00	1,206.00	-
4-Thio-U-TOM-CE Phosphoramidite	10-3052-90	100 µmol	611.00	587.00	525.00	-
4-Thio-U-TOM-CE Phosphoramidite	10-3052-95	50 µmol	306.00	294.00	263.00	-
5-Br-U-CE Phosphoramidite	10-3090-02	0.25 g	748.00	718.00	643.00	-
5-Br-U-CE Phosphoramidite	10-3090-90	100 µmol	309.00	297.00	266.00	-
5-I-U-CE Phosphoramidite	10-3091-02	0.25 g	848.00	814.00	729.00	-
5-I-U-CE Phosphoramidite	10-3091-90	100 µmol	345.00	331.00	297.00	-
5-Me-C-2'-MOE-Phosphoramidite - 0.5 g	10-3211-05	0.5 g	53.90	51.70	46.40	-
5-Me-C-2'-MOE-Phosphoramidite - 0.5 g	10-3211-05E	0.5 g	53.90	51.70	46.40	-
5-Me-C-2'-MOE-Phosphoramidite - 2.0 g	10-3211-20	2.0 g	216.00	207.00	186.00	-
5-Me-C-TOM-CE Phosphoramidite	10-3064-02	0.25 g	683.00	656.00	587.00	-
5-Me-C-TOM-CE Phosphoramidite	10-3064-90	100 µmol	273.00	262.00	235.00	-
5-Me-C-TOM-CE Phosphoramidite	10-3064-95	50 µmol	137.00	132.00	118.00	-
5-Me-U-2'-MOE-Phosphoramidite - 0.5 g	10-3231-05	0.5 g	53.90	51.70	46.40	-
5-Me-U-2'-MOE-Phosphoramidite - 0.5 g	10-3231-05E	0.5 g	53.90	51.70	46.40	-
5-Me-U-2'-MOE-Phosphoramidite - 2.0 g	10-3231-20	2.0 g	216.00	207.00	186.00	-
5-Me-U-CE Phosphoramidite	10-3050-02	0.25 g	683.00	656.00	587.00	-
5-Me-U-CE Phosphoramidite	10-3050-90	100 µmol	273.00	262.00	235.00	-
6-Fluorescein Phosphoramidite - 0.25 g	10-1964-02E	0.25 g	856.00	822.00	736.00	-
6-Fluorescein Phosphoramidite, Expedite	10-1964-95E	50 µmol	237.00	228.00	204.00	-
A-2'-MOE-Phosphoramidite - 0.5 g	10-3200-05	0.5 g	53.90	51.70	46.40	-
A-2'-MOE-Phosphoramidite - 0.5 g	10-3200-05E	0.5 g	53.90	51.70	46.40	-
A-2'-MOE-Phosphoramidite - 2.0 g	10-3200-20	2.0 g	216.00	207.00	186.00	-
Ac-A-RNA-CPG	20-3304-01	0.1 g	57.50	55.20	49.50	-
Ac-A-RNA-CPG	20-3304-10	1g	511.00	491.00	439.00	-
Ac-C-CE Phosphoramidite	10-3015-02	0.25 g	35.40	34.00	30.40	-
Ac-C-CE Phosphoramidite	10-3015-05	0.5 g	53.10	51.00	45.70	-
Ac-C-CE Phosphoramidite	10-3015-10	1g	99.60	95.60	85.70	-
Ac-C-CE Phosphoramidite - Expedite - 0.25 g	10-3015-02E	0.25 g	35.40	34.00	30.40	-
Ac-C-CE Phosphoramidite MerMade	10-3015-10M	1g	99.60	95.60	85.70	-
Ac-C-CE Phosphoramidite, Expedite	10-3015-05E	0.5 g	53.10	51.00	45.70	-
Ac-C-RNA 500 Å	20-3415-41A	4x1 µmol	144.00	138.00	124.00	-
Ac-C-RNA 500 Å - 1 µmol column	20-3415-41E	144.00	138.00	124.00	-	
Ac-C-RNA 500 Å - 0.2µmol col (4/pack)	20-3415-42A	108.00	104.00	92.90	-	
Ac-C-RNA 500Å-0.2µmol col.Exp.	20-3415-42E	108.00	104.00	92.90	-	
Ac-C-RNA-CPG	20-3315-01	0.1 g	57.50	55.20	49.50	-
Ac-C-RNA-CPG	20-3315-02	0.25 g	137.00	132.00	118.00	-
Ac-C-RNA-CPG	20-3315-10	1g	511.00	491.00	439.00	-
Ac-C-RNA-CPG	20-3415-13	1 x 10 µmol	324.00	311.00	279.00	-
Ac-C-RNA-CPG	20-3415-14	1 x 15 µmol	431.00	414.00	371.00	-
Ac-C-RNA-CPG	20-3415-41	4 x 1 µmol	144.00	138.00	124.00	-
Ac-C-RNA-CPG	20-3415-42	4 x 0.2 µmol	108.00	104.00	92.90	-
Ac-G-CE Phosphoramidite MerMade	10-3025-10M	1g	99.60	95.60	85.70	-
Ac-G-RNA-CPG	20-3324-02	0.25 g	137.00	132.00	118.00	-
Ac-G-RNA-CPG	20-3324-10	1g	511.00	491.00	439.00	-
Amino-Modifier Serinol Phosphoramidite	10-1997-02M	0.25 g	856.00	822.00	736.00	-
A-RNA-TOM-CPG	20-3304-02	0.25 g	137.00	132.00	118.00	-
A-RNA-TOM-CPG	20-3404-13	1 x 10 µmol	324.00	311.00	279.00	-
A-RNA-TOM-CPG	20-3404-14	1 x 15 µmol	431.00	414.00	371.00	-
A-RNA-TOM-CPG	20-3404-41	4 x 1 µmol	144.00	138.00	124.00	-
A-RNA-TOM-CPG	20-3404-42	4 x 0.2 µmol	108.00	104.00	92.90	-
a-Tocopherol-TEG Phosphoramidite 0.25g	10-1977-02M	1g	949.00	911.00	816.00	-
A-TOM-CE Phosphoramidite	10-3004-02	0.25 g	83.00	79.70	71.40	-
A-TOM-CE Phosphoramidite	10-3004-05	0.5 g	166.00	159.00	143.00	-
A-TOM-CE Phosphoramidite	10-3004-10	1g	304.00	292.00	261.00	-
A-TOM-CE Phosphoramidite MerMade	10-3004-10M	1g	304.00	292.00	261.00	-
A-TOM-CE Phosphoramidite, Expedite	10-3004-10E	1g	304.00	292.00	261.00	-
A-TOM-CE Phosphoramidite, Expedite - 0.5 g	10-3004-05E	0.5 g	166.00	159.00	143.00	-
A-TOM-RNA 1000 (0.2µm col.) 4p	20-3404-42E	108.00	104.00	92.90	-	
beta-L-Pac-dA-CE Phosphoramidite, 0.25 g	10-2101-02	0.25 g	216.00	207.00	186.00	-
BHQ-2-dT, Expedite - 50 µmol	10-5942-95E	50 µmol	532.00	511.00	458.00	-
Bz-A-CE Phosphoramidite	10-3003-02	0.25 g	35.40	34.00	30.40	-
Bz-A-CE Phosphoramidite	10-3003-05	0.5 g	53.10	51.00	45.70	-
Bz-A-CE Phosphoramidite	10-3003-10	1g	99.60	95.60	85.70	-
Bz-A-CE Phosphoramidite MerMade	10-3003-10M	1g	99.60	95.60	85.70	-
Bz-A-RNA 500 Å - 0.2 mol col.	20-3403-42E	108.00	104.00	92.90	-	
Bz-A-RNA-CPG	20-3303-01	0.1 g	57.50	55.20	49.50	-
Bz-A-RNA-CPG	20-3303-02	0.25 g	137.00	132.00	118.00	-
Bz-A-RNA-CPG	20-3303-10	1g	511.00	491.00	439.00	-
Bz-A-RNA-CPG	20-3403-13	1 x 10 µmol	324.00	311.00	279.00	-
Bz-A-RNA-CPG	20-3403-14	1 x 15 µmol	431.00	414.00	371.00	-
Bz-A-RNA-CPG	20-3403-41	4 x 1 µmol	144.00	138.00	124.00	-
Bz-A-RNA-CPG	20-3403-42	4 x 0.2 µmol	108.00	104.00	92.90	-
Bz-A-RNA-CPG, Expedite - 4 x 1 µmol	20-3403-41E	4 x 1 µmol	144.00	138.00	124.00	-

Description	Reference	Scale/Quantity	Price List 2025			
			EUR	CHF	GBP	USD
Chemical Phosphorylation Reage	10-1900-90E		49.80	47.80	42.80	-
C-TOM-CE Phosphoramidite	10-3014-02	0.25 g	83.00	79.70	71.40	-
C-TOM-CE Phosphoramidite	10-3014-05	0.5 g	166.00	159.00	143.00	-
C-TOM-CE Phosphoramidite	10-3014-10	1 g	304.00	292.00	261.00	-
C-TOM-CE Phosphoramidite MerMade	10-3014-10M	1 g	304.00	292.00	261.00	-
C-TOM-CE Phosphoramidite, Expedite	10-3014-10E	1 g	304.00	292.00	261.00	-
C-TOM-CE Phosphoramidite, Expedite - 0.5 g	10-3014-05E	0.5 g	166.00	159.00	143.00	-
Cyanine 3 CPG - 4 x 0.2 µmol	20-5913-42	4 x 0.2 µmol	129.00	124.00	111.00	-
DBCO-Serinol Phosphoramidite 100 µmol	10-1998-90	100 µmol	518.00	497.00	445.00	-
dmf-dG-5'-CE Phosphoramidite, Exp. 0.5 g	10-9201-05E	0.5 g	164.00	157.00	141.00	-
dT-5'-CE Phosphoramidite, Expedite 0.5 g	10-0301-05E	0.5 g	164.00	157.00	141.00	-
dX-CE Phosphoramidite - 0.25 g	10-1537-02	0.25 g	1,424.00	1,367.00	1,225.00	-
dX-CE Phosphoramidite - 100 µmol	10-1537-90	100 µmol	618.00	593.00	531.00	-
Fluorescein Phosphoramidite 0.25 g	10-1963-02M		856.00	822.00	736.00	-
Fluorescein-dT Phosphoramidite, Expedite	10-1056-95E	50 µmol	232.00	223.00	200.00	-
G-2'-MOE-Phosphoramidite - 0.5 g	10-3220-05	0.5 g	53.90	51.70	46.40	-
G-2'-MOE-Phosphoramidite - 0.5 g	10-3220-05E		53.90	51.70	46.40	-
G-2'-MOE-Phosphoramidite - 2.0 g	10-3220-20	2.0 g	216.00	207.00	186.00	-
Glen Gel-Pak™ 0.2 Desalting Column	61-5002-50	Pk / 50	647.00	621.00	556.00	-
Glen reagent Bz-A-CE Phosphoramidite, 0.5grams, Expedite	10-3003-05E		53.10	51.00	45.70	-
G-RNA-TOM-CPG	20-3324-01	0.1 g	57.50	55.20	49.50	-
G-RNA-TOM-CPG	20-3424-13	1 x 10 µmol	324.00	311.00	279.00	-
G-RNA-TOM-CPG	20-3424-14	1 x 15 µmol	431.00	414.00	371.00	-
G-RNA-TOM-CPG	20-3424-41	4 x 1 µmol	144.00	138.00	124.00	-
G-RNA-TOM-CPG	20-3424-42	4 x 0.2 µmol	108.00	104.00	92.90	-
G-TOM-CE Phosphoramidite	10-3024-02	0.25 g	83.00	79.70	71.40	-
G-TOM-CE Phosphoramidite	10-3024-05	0.5 g	166.00	159.00	143.00	-
G-TOM-CE Phosphoramidite	10-3024-10	1 g	304.00	292.00	261.00	-
G-TOM-CE Phosphoramidite MerMade	10-3024-10M	1 g	304.00	292.00	261.00	-
G-TOM-CE Phosphoramidite, Expedite	10-3024-10E	1 g	304.00	292.00	261.00	-
G-TOM-CE Phosphoramidite, Expedite - 0.5 g	10-3024-05E	0.5 g	166.00	159.00	143.00	-
G-TOM-RNA 1000 (0.2µm col.) 4p	20-3424-42E		108.00	104.00	92.90	-
G-TOM-RNA 1000 (1µm col.) 4pcs	20-3424-41E		144.00	138.00	124.00	-
iPr-Pac-G-CE Phosphoramidite	10-3021-02	0.25 g	115.00	110.00	98.90	-
iPr-Pac-G-CE Phosphoramidite	10-3021-05	0.5 g	230.00	221.00	198.00	-
iPr-Pac-G-CE Phosphoramidite	10-3021-10	1 g	460.00	442.00	396.00	-
iPr-Pac-G-CE Phosphoramidite, Expedite	10-3021-05E	0.5 g	230.00	221.00	198.00	-
iPr-Pac-G-CE Phosphoramidite, Expedite - 1 g	10-3021-10E	1 g	460.00	442.00	396.00	-
iPr-Pac-G-CE Phosphoramidite, Pharmacia - 0.5 g	10-3021-05P	0.5 g	230.00	221.00	198.00	-
iPr-Pac-G-RNA 500 Å	20-3421-41A	4x1 µmol	144.00	138.00	124.00	-
iPr-Pac-G-RNA 500 Å - 1 µmol T	20-3421-41P		144.00	138.00	124.00	-
iPr-Pac-G-RNA 500 Å -0.2µmol col (4/pack)	20-3421-42A		108.00	104.00	92.90	-
iPr-Pac-G-RNA-CPG	20-3321-01	0.1 g	57.50	55.20	49.50	-
iPr-Pac-G-RNA-CPG	20-3321-02	0.25 g	137.00	132.00	118.00	-
iPr-Pac-G-RNA-CPG	20-3421-13	1 x 10 µmol	324.00	311.00	279.00	-
iPr-Pac-G-RNA-CPG	20-3421-14	1 x 15 µmol	431.00	414.00	371.00	-
iPr-Pac-G-RNA-CPG	20-3421-41	4 x 1 µmol	144.00	138.00	124.00	-
iPr-Pac-G-RNA-CPG	20-3421-42	4 x 0.2 µmol	108.00	104.00	92.90	-
N6-Methyl-A-CE Phosphoramidite	10-3005-02	0.25 g	2,193.00	2,105.00	1,886.00	-
Pac-A-CE Phosphoramidite	10-3000-02	0.25 g	108.00	104.00	92.90	-
Pac-A-CE Phosphoramidite	10-3000-05	0.5 g	216.00	207.00	186.00	-
Pac-A-CE Phosphoramidite	10-3000-10	1 g	395.00	379.00	340.00	-
Pac-A-CE Phosphoramidite, Expedite	10-3000-05E	0.5 g	216.00	207.00	186.00	-
Pac-A-RNA 500 Å	20-3400-41A	4x1 µmol	144.00	138.00	124.00	-
Pac-A-RNA 500 Å - 0.2 µmol col(4/pack)	20-3400-42A		108.00	104.00	92.90	-
Pac-A-RNA-CPG	20-3300-01	0.1 g	57.50	55.20	49.50	-
Pac-A-RNA-CPG	20-3300-02	0.25 g	137.00	132.00	118.00	-
Pac-A-RNA-CPG	20-3400-13	1 x 10 µmol	324.00	311.00	279.00	-
Pac-A-RNA-CPG	20-3400-14	1 x 15 µmol	431.00	414.00	371.00	-
Pac-A-RNA-CPG	20-3400-41	4 x 1 µmol	144.00	138.00	124.00	-
Pac-A-RNA-CPG	20-3400-42	4 x 0.2 µmol	108.00	104.00	92.90	-
PseudoUridine-CE Phosphoramidite 0.25 g	10-3055-02E		1,790.00	1,718.00	1,539.00	-
Pyrrolo-C-TOM-CE Phosphoramidite	10-3017-02	0.25 g	1,798.00	1,726.00	1,546.00	-
Pyrrolo-C-TOM-CE Phosphoramidite	10-3017-90	100 µmol	712.00	684.00	612.00	-
Pyrrolo-C-TOM-CE Phosphoramidite	10-3017-95	50 µmol	370.00	355.00	318.00	-
SIMA (HEX) Phosphoramidite, 100 µmol	10-5905-90E	100 µmol	273.00	262.00	235.00	-
SIMA (HEX) Phosphoramidite, 50 µmol	10-5905-95	50 µmol	137.00	132.00	118.00	-
U-CE Phosphoramidite	10-3030-02	0.25 g	35.40	34.00	30.40	-
U-CE Phosphoramidite	10-3030-05	0.5 g	53.10	51.00	45.70	-
U-CE Phosphoramidite	10-3030-10	1 g	99.60	95.60	85.70	-
U-CE Phosphoramidite MerMade	10-3030-10M	1 g	99.60	95.60	85.70	-
U-CE Phosphoramidite, Expedite	10-3030-05E	0.5 g	53.10	51.00	45.70	-
U-CE Phosphoramidite, Expedite - 0.25 g	10-3030-02E	0.25 g	35.40	34.00	30.40	-
U-RNA 500 Å - 0.2 µmol col (4/pack)	20-3430-42A		108.00	104.00	92.90	-
U-RNA 500 Å - 1.0 µmol columns	20-3430-41E	0.2 µmol	144.00	138.00	124.00	-
U-RNA-CPG	20-3330-01	0.1 g	57.50	55.20	49.50	-
U-RNA-CPG	20-3330-02	0.25 g	137.00	132.00	118.00	-
U-RNA-CPG	20-3330-10	1 g	511.00	491.00	439.00	-
U-RNA-CPG	20-3430-13	1 x 10 µmol	324.00	311.00	279.00	-
U-RNA-CPG	20-3430-14	1 x 15 µmol	431.00	414.00	371.00	-
U-RNA-CPG	20-3430-41	4 x 1 µmol	144.00	138.00	124.00	-
U-RNA-CPG	20-3430-42	4 x 0.2 µmol	108.00	104.00	92.90	-
U-RNA-CPG, Expedite - 4 x 0.2 µmol	20-3430-42E	4 x 0.2 µmol	108.00	104.00	92.90	-
U-TOM-CE Phosphoramidite	10-3034-02	0.25 g	83.00	79.70	71.40	-

Description	Reference	Scale/Quantity	Price List 2025			
			EUR	CHF	GBP	USD
U-TOM-CE Phosphoramidite	10-3034-05	0.5 g	166.00	159.00	143.00	-
U-TOM-CE Phosphoramidite	10-3034-10	1 g	304.00	292.00	261.00	-
U-TOM-CE Phosphoramidite MerMade	10-3034-10M	1 g	304.00	292.00	261.00	-
U-TOM-CE Phosphoramidite, Expedite - 0.5 g	10-3034-05E	0.5 g	166.00	159.00	143.00	-

All minor bases, RNA products and modifiers are packaged in septum capped vials suitable for ABI and other instruments. If you would like another type of vial/column

please choose the reference ending by Monomers

E = Expedite

B = Beckman Oligo 1000

P = Pharmacia Gene Assembler

M = MerMade Columns