



# 2021 Price List

Products for DNA research exclusively  
distributed by Eurogentec



Description	Reference	Scale/Quantity	Price List 2021				Ice Shipping
			EUR	CHF	GBP	USD	
<b>Glen Research products for oligonucleotide synthesis</b>							<b>EA1</b>
<b>GlenDNA</b>							<b>EA1A12</b>
0.2 M Saccharin 1-Methylimidazole (SMI) in Acetonitrile	30-3182-52	200 ml	154.00	169.00	142.00	-	
0.2M Saccharin 1-Methylimidazole (SMI) in Acetonitrile	30-3180-52	200 mL	154.00	169.00	142.00	-	
1c-dC-CE Phosphoramidite	10-1015-40	4 g	56.50	62.20	52.00	-	
1-Me-A-CE Phosphoramidite	10-3501-02	0.25 g	1,310.00	1,441.00	1,205.00	-	
1-Me-A-CE Phosphoramidite	10-3501-95	50 µmol	275.00	303.00	253.00	-	
3'-Phosphate CPG (High Load)	25-2900-46	4 x 2.5 µmol	157.00	173.00	144.00	-	
5-aza-5,6-dihydro-dC-CE Phosphoramidite	10-1511-02	0.25 g	1,468.00	1,615.00	1,351.00	-	
5-aza-5,6-dihydro-dC-CE Phosphoramidite	10-1511-90	100 µmol	472.00	519.00	434.00	-	
5-aza-5,6-dihydro-dC-CE Phosphoramidite	10-1511-95	50 µmol	236.00	260.00	217.00	-	
5-I-U-CE Phosphoramidite	10-3091-95	50 µmol	128.00	141.00	118.00	-	
5-Me-dC Brancher Phosphoramidite	10-1018-02	0.25 g	662.00	728.00	609.00	-	
5-Me-dC Brancher Phosphoramidite	10-1018-90	100 µmol	269.00	296.00	247.00	-	
5-Me-dC Brancher Phosphoramidite, Expedite	10-1018-02E	0.25 g	662.00	728.00	609.00	-	
5-Me-U-CE Phosphoramidite	10-3050-95	50 µmol	124.00	136.00	114.00	-	
8-D-dmf-CE Phosphoramidite	10-1520-90	100 µmol	129.00	142.00	119.00	-	
8-D-dmf-dG-CE Phosphoramidite	10-1520-02	0.25 g	344.00	378.00	316.00	-	
AAA Trimer Phosphoramidite	13-1000-90	100 µmol	1,114.00	1,225.00	1,025.00	-	
AAA Trimer Phosphoramidite	13-1000-95	50 µmol	557.00	613.00	512.00	-	
AAC Trimer Phosphoramidite	13-1001-90	100 µmol	1,114.00	1,225.00	1,025.00	-	
AAC Trimer Phosphoramidite	13-1001-95	50 µmol	557.00	613.00	512.00	-	
AB3900 Ac-dC-CPG 1000A,1.0 µmol	20-2115-61	200 pce	706.00	777.00	650.00	-	
AB3900 dmf-dG-CPG 1000A,1.0 µmol	20-2129-61	200 pce	706.00	777.00	650.00	-	
AB3900 dt-CPG 1000A,1.0 µmol	20-2131-61	200 pce	706.00	777.00	650.00	-	
Ac-dC-CE Phosphoramidite	10-1015-02	0.25 g	3.53	3.88	3.25	-	
Ac-dC-CE Phosphoramidite	10-1015-05	0.5 g	7.06	7.77	6.50	-	
Ac-dC-CE Phosphoramidite	10-1015-05M	0.5 g	7.06	7.77	6.50	-	
Ac-dC-CE Phosphoramidite	10-1015-10	1 g	14.10	15.50	13.00	-	
Ac-dC-CE Phosphoramidite	10-1015-10M	1 g	14.10	15.50	13.00	-	
Ac-dC-CE Phosphoramidite	10-1015-1B	1 g	16.50	18.20	15.20	-	
Ac-dC-CE Phosphoramidite	10-1015-1C	1 g	14.10	15.50	13.00	-	
Ac-dC-CE Phosphoramidite	10-1015-1P	1 g	16.50	18.20	15.20	-	
Ac-dC-CE Phosphoramidite	10-1015-15	10 g	141.00	155.00	130.00	-	
Ac-dC-CE Phosphoramidite	10-1015-20	2 g	28.20	31.00	25.90	-	
Ac-dC-CE Phosphoramidite	10-1015-2C	2 g	28.20	31.00	25.90	-	
Ac-dC-CE Phosphoramidite	10-1015-50	5 g	70.60	77.70	65.00	-	
Ac-dC-CE Phosphoramidite	10-1015-B2	0.25 g	4.10	4.51	3.77	-	
Ac-dC-CE Phosphoramidite	10-1015-B5	0.5 g	8.30	9.13	7.64	-	
Ac-dC-CE Phosphoramidite	10-1015-C2	0.25 g	3.53	3.88	3.25	-	
Ac-dC-CE Phosphoramidite	10-1015-C5	0.5 g	7.06	7.77	6.50	-	
Ac-dC-CE Phosphoramidite	10-1015-P4	0.4 g	6.60	7.26	6.07	-	
Ac-dC-CE Phosphoramidite	10-1015-02M	0.25 g	3.53	3.88	3.25	-	
Ac-dC-CE Phosphoramidite, Expedite	10-1015-02E	0.25 g	3.53	3.88	3.25	-	
Ac-dC-CE Phosphoramidite, Expedite	10-1015-05E	0.5 g	7.06	7.77	6.50	-	
Ac-dC-CE Phosphoramidite, Pharmacia	10-1015-02P	0.25 g	3.53	3.88	3.25	-	
Ac-dC-CE Phosphoramidite, Pharmacia	10-1015-05P	0.5 g	7.06	7.77	6.50	-	
Ac-dC-CPG 1000	20-2015-01	0.1 g	5.81	6.39	5.35	-	
Ac-dC-CPG 1000	20-2015-02	0.25 g	12.90	14.20	11.90	-	
Ac-dC-CPG 1000	20-2015-10	1 g	48.40	53.20	44.50	-	
Ac-dC-CPG 1000	20-2015-61	48 x 1 µmol	242.00	266.00	223.00	-	
Ac-dC-CPG 1000	20-2015-62	200 x 0.2 µmol	605.00	666.00	557.00	-	
Ac-dC-CPG 1000	20-2115-13	1 x 10 µmol	80.60	88.70	74.20	-	
Ac-dC-CPG 1000	20-2115-41	4 x 1 µmol	28.20	31.00	25.90	-	
Ac-dC-CPG 1000	20-2115-42	4 x 0.2 µmol	28.20	31.00	25.90	-	
Ac-dC-CPG 1000	20-2115-45	4 x 40 nmol	28.20	31.00	25.90	-	
Ac-dC-CPG 1000	20-2115-62	200 x 200 nmol	524.00	576.00	482.00	-	
Ac-dC-CPG 1000	20-2115-65	200 x 40 nmol	484.00	532.00	445.00	-	
Ac-dC-CPG 1000	20-2215-14	1 x 15 µmol	121.00	133.00	111.00	-	
Ac-dC-CPG 1000	20-2215-41	4 x 1 µmol	20.20	22.20	18.60	-	
Ac-dC-CPG 1000	20-2215-42	4 x 0.2 µmol	14.10	15.50	13.00	-	
Ac-dC-CPG 1000	20-2215-45	4 x 40 nmol	14.10	15.50	13.00	-	
Ac-dC-CPG 1000	20-2515-P1	10 x 1.3 µmol	106.00	117.00	98.00	-	
Ac-dC-CPG 1000	20-2515-P2	10 x 0.2 µmol	106.00	117.00	98.00	-	
Ac-dC-CPG 500	20-2013-01	0.1 g	5.81	6.39	5.35	-	
Ac-dC-CPG 500	20-2013-02	0.25 g	12.90	14.20	11.90	-	
Ac-dC-CPG 500	20-2113-13	1 x 10 µmol	80.60	88.70	74.20	-	
Ac-dC-CPG 500	20-2113-41	4 x 1 µmol	28.20	31.00	25.90	-	
Ac-dC-CPG 500	20-2113-42	4 x 0.2 µmol	28.20	31.00	25.90	-	
Ac-dC-CPG 500	20-2213-14	1 x 15 µmol	121.00	133.00	111.00	-	
Ac-dC-CPG 500	20-2213-41	4 x 1 µmol	20.20	22.20	18.60	-	
Ac-dC-CPG 500	20-2213-42	4 x 0.2 µmol	14.10	15.50	13.00	-	
Ac-dC-PACE Phosphoramidite	10-1150-02	0.25 g	178.00	196.00	164.00	-	
Ac-dC-PACE Phosphoramidite	10-1150-05	0.5 g	357.00	393.00	328.00	-	
Ac-dC-PACE Phosphoramidite	10-1150-10	1 g	713.00	784.00	656.00	-	
Ac-G-CE Phosphoramidite	10-3025-10	1 g	90.70	99.80	83.40	-	
Ac-G-CE Phosphoramidite	10-3025-02	0.25 g	32.30	35.50	29.70	-	
ACT Trimer Phosphoramidite	13-1013-90	100 µmol	1,114.00	1,225.00	1,025.00	-	
ACT Trimer Phosphoramidite	13-1013-95	50 µmol	557.00	613.00	512.00	-	
ATC Trimer Phosphoramidite	13-1031-90	100 µmol	1,114.00	1,225.00	1,025.00	-	
ATC Trimer Phosphoramidite	13-1031-95	50 µmol	557.00	613.00	512.00	-	
ATG Trimer Phosphoramidite	13-1032-90	100 µmol	1,114.00	1,225.00	1,025.00	-	
ATG Trimer Phosphoramidite	13-1032-95	50 µmol	557.00	613.00	512.00	-	
Br-U-CE Phosphoramidite	10-3090-95	50 µmol	128.00	141.00	118.00	-	
CAG Trimer Phosphoramidite	13-1102-90	100 µmol	1,114.00	1,225.00	1,025.00	-	
CAG Trimer Phosphoramidite	13-1102-95	50 µmol	557.00	613.00	512.00	-	
CAT Trimer Phosphoramidite	13-1103-90	100 µmol	1,114.00	1,225.00	1,025.00	-	
CAT Trimer Phosphoramidite	13-1103-95	50 µmol	557.00	613.00	512.00	-	
CCG Trimer Phosphoramidite	13-1112-90	100 µmol	1,114.00	1,225.00	1,025.00	-	
CCG Trimer Phosphoramidite	13-1112-95	50 µmol	557.00	613.00	512.00	-	
CGT Trimer Phosphoramidite	13-1123-90	100 µmol	1,114.00	1,225.00	1,025.00	-	

Description	Reference	Scale/Quantity	Price List 2021				Ice Shipping
			EUR	CHF	GBP	USD	
CGT Trimer Phosphoramidite	13-1123-95	50 µmol	557.00	613.00	512.00	-	
CleanAmp™-Ac-dC-CE Phosphoramidite	10-1450-05	0.5 g	650.00	715.00	598.00	-	
CleanAmp™-Pac-dA-CE Phosphoramidite	10-1440-05	0.5 g	650.00	715.00	598.00	-	
CleanAmp™-Pac-dG-CE Phosphoramidite	10-1460-05	0.5 g	650.00	715.00	598.00	-	
CTG Trimer Phosphoramidite	13-1132-90	100 µmol	1,114.00	1,225.00	1,025.00	-	
CTG Trimer Phosphoramidite	13-1132-95	50 µmol	557.00	613.00	512.00	-	
dA LV-PS	26-2100-42	4 x 0.2 µmol	38.70	42.60	35.60	-	
dA LV-PS	26-2100-45	4 x 40 nmol	38.70	42.60	35.60	-	
dA PS	26-2600-62	200 x 200 nmol	650.00	715.00	598.00	-	
dA PS	26-2600-65	200 x 40 nmol	650.00	715.00	598.00	-	
dA, dC, dG, dT-CPG 1000	20-2241-41	4 x 1 µmol	20.20	22.20	18.60	-	
dA, dC, dG, dT-CPG 1000	20-2241-42	4 x 0.2 µmol	14.10	15.50	13.00	-	
dA, dC, dG, dT-CPG 1000	20-2241-45	4 x 40 nmol	14.10	15.50	13.00	-	
dA,dC,dG,dT CPG 1000	20-2141-41	4 x 1 µmol	28.20	31.00	25.90	-	
dA,dC,dG,dT CPG 1000	20-2141-42	4 x 0.2 µmol	28.20	31.00	25.90	-	
dA,dC,dG,dT CPG 1000	20-2141-45	4 x 40 nmol	28.20	31.00	25.90	-	
dA,dC,dG,dT- LV-PS 0.2µm Columns	26-2140-42	4 x 0.2 µmol	38.70	42.60	35.60	-	
dA,dC,dG,dT- LV-PS 40nm Columns	26-2140-45	4 x 40 nmol	38.70	42.60	35.60	-	
dA,dC,dG,dT-CPG 2000	20-2142-42	4 x 0.2 µmol	30.80	33.90	28.30	-	
dA,dC,dG,dT-CPG 2000	20-2242-42	4 x 0.2 µmol	14.10	15.50	13.00	-	
dA,dC,dG,dT-CPG 500	20-2140-41	4 x 1 µmol	28.20	31.00	25.90	-	
dA,dC,dG,dT-CPG 500	20-2140-42	4 x 0.2 µmol	28.20	31.00	25.90	-	
dA,dC,dG,dT-CPG 500	20-2240-41	4 x 1 µmol	20.20	22.20	18.60	-	
dA,dC,dG,dT-CPG 500	20-2240-42	4 x 0.2 µmol	14.10	15.50	13.00	-	
dA-CE Phosphoramidite	10-1000-02	0.25 g	3.53	3.88	3.25	-	
dA-CE Phosphoramidite	10-1000-05	0.5 g	7.06	7.77	6.50	-	
dA-CE Phosphoramidite	10-1000-10	1 g	14.10	15.50	13.00	-	
dA-CE Phosphoramidite	10-1000-1B	1 g	15.40	16.90	14.20	-	
dA-CE Phosphoramidite	10-1000-1C	1 g	14.10	15.50	13.00	-	
dA-CE Phosphoramidite	10-1000-1P	1 g	15.40	16.90	14.20	-	
dA-CE Phosphoramidite	10-1000-1S	10 g	141.00	155.00	130.00	-	
dA-CE Phosphoramidite	10-1000-20	2 g	28.20	31.00	25.90	-	
dA-CE Phosphoramidite	10-1000-2C	2 g	28.20	31.00	25.90	-	
dA-CE Phosphoramidite	10-1000-40	4 g	56.50	62.20	52.00	-	
dA-CE Phosphoramidite	10-1000-50	5 g	70.60	77.70	65.00	-	
dA-CE Phosphoramidite	10-1000-5S	5 g	70.60	77.70	65.00	-	
dA-CE Phosphoramidite	10-1000-B2	0.25 g	3.90	4.29	3.59	-	
dA-CE Phosphoramidite	10-1000-B5	0.5 g	7.70	8.47	7.08	-	
dA-CE Phosphoramidite	10-1000-C2	0.25 g	3.53	3.88	3.25	-	
dA-CE Phosphoramidite	10-1000-C5	0.5 g	7.06	7.77	6.50	-	
dA-CE Phosphoramidite, Expedite	10-1000-05E	0.5 g	7.06	7.77	6.50	-	
dA-CE Phosphoramidite, Mermade	10-1000-02M	0.25 g	3.53	3.88	3.25	-	
dA-CE Phosphoramidite, Mermade	10-1000-05M	0.5 g	7.06	7.77	6.50	-	
dA-CE Phosphoramidite, Mermade	10-1000-10M	1 g	14.10	15.50	13.00	-	
dA-CE Phosphoramidite, Pharmacia	10-1000-P4	0.4 g	6.20	6.82	5.70	-	
dA-CPG 1 µmol columns	20-2101-61	200 pce	706.00	777.00	650.00	-	
dA-CPG 1000	20-2001-01	0.1 g	5.81	6.39	5.35	-	
dA-CPG 1000	20-2001-02	0.25 g	12.90	14.20	11.90	-	
dA-CPG 1000	20-2001-10	1 g	48.40	53.20	44.50	-	
dA-CPG 1000	20-2001-61	48 x 1 µmol	242.00	266.00	223.00	-	
dA-CPG 1000	20-2001-62	200 x 0.2 µmol	605.00	666.00	557.00	-	
dA-CPG 1000	20-2101-13	1 x 10 µmol	80.60	88.70	74.20	-	
dA-CPG 1000	20-2101-41	4 x 1 µmol	28.20	31.00	25.90	-	
dA-CPG 1000	20-2101-42	4 x 0.2 µmol	28.20	31.00	25.90	-	
dA-CPG 1000	20-2101-45	4 x 40 nmol	28.20	31.00	25.90	-	
dA-CPG 1000	20-2101-62	200 x 200 nm	524.00	576.00	482.00	-	
dA-CPG 1000	20-2101-65	200 x 40 nmol	484.00	532.00	445.00	-	
dA-CPG 1000	20-2201-14	1 x 15 µmol	121.00	133.00	111.00	-	
dA-CPG 1000	20-2201-41	4 x 1 µmol	20.20	22.20	18.60	-	
dA-CPG 1000	20-2201-42	4 x 0.2 µmol	14.10	15.50	13.00	-	
dA-CPG 1000	20-2201-45	4 x 40 nmol	14.10	15.50	13.00	-	
dA-CPG 1000, Expedite	20-2201-42E	4 x 0.2 µmol	14.10	15.50	13.00	-	
dA-CPG 2000	20-2002-01	0.1 g	15.10	16.60	13.90	-	
dA-CPG 2000	20-2002-02	0.25 g	30.20	33.20	27.80	-	
dA-CPG 2000	20-2002-10	1 g	106.00	117.00	98.00	-	
dA-CPG 2000	20-2102-42	4 x 0.2 µmol	28.20	31.00	25.90	-	
dA-CPG 2000	20-2202-42	4 x 0.2 µmol	14.10	15.50	13.00	-	
dA-CPG 500	20-2000-01	0.1 g	5.81	6.39	5.35	-	
dA-CPG 500	20-2000-02	0.25 g	12.90	14.20	11.90	-	
dA-CPG 500	20-2000-10	1 g	48.40	53.20	44.50	-	
dA-CPG 500	20-2100-13	1 x 10 µmol	80.60	88.70	74.20	-	
dA-CPG 500	20-2100-41	4 x 1 µmol	28.20	31.00	25.90	-	
dA-CPG 500	20-2100-42	4 x 0.2 µmol	28.20	31.00	25.90	-	
dA-CPG 500	20-2200-14	1 x 15 µmol	121.00	133.00	111.00	-	
dA-CPG 500	20-2200-41	4 x 1 µmol	20.20	22.20	18.60	-	
dA-CPG 500	20-2200-42	4 x 0.2 µmol	14.10	15.50	13.00	-	
dA-CPG 500	20-2500-P2	10 x 0.2 µmol	106.00	117.00	98.00	-	
dA-High Load-CPG	25-2000-02	0.25 g	16.10	17.70	14.80	-	
dA-High Load-CPG	25-2000-10	1 g	58.10	63.90	53.50	-	
dA-High Load-CPG	25-2100-17	1 x 25 µmol	101.00	111.00	93.00	-	
dA-High Load-CPG	25-2100-46	4 x 2.5 µmol	60.50	66.60	55.70	-	
dA-High Load-CPG	25-2200-18	1 x 35 µmol	149.00	164.00	137.00	-	
dA-High Load-CPG	25-2200-46	4 x 2.5 µmol	60.50	66.60	55.70	-	
dA-PACE Phosphoramidite	10-1140-02	0.25 g	178.00	196.00	164.00	-	
dA-PACE Phosphoramidite	10-1140-05	0.5 g	357.00	393.00	328.00	-	
dA-PACE Phosphoramidite	10-1140-10	1 g	713.00	784.00	656.00	-	
dA-Thiophosphoramidite	10-1700-90	100 µmol	229.00	252.00	211.00	-	
dA-Thiophosphoramidite,	10-1700-02	0.25 g	550.00	605.00	506.00	-	
dC LV-PS	26-2110-42	4 x 0.2 µmol	38.70	42.60	35.60	-	
dC LV-PS	26-2110-45	4 x 40 nmol	38.70	42.60	35.60	-	
dC PS	26-2610-62	200 x 200 nmol	650.00	715.00	598.00	-	

Description	Reference	Scale/Quantity	Price List 2021				Ice Shipping
			EUR	CHF	GBP	USD	
dC PS	26-2610-65	200 x 40 nmol	650.00	715.00	598.00	-	
dC Thiophosphoramidite,	10-1710-02	0.25 g	550.00	605.00	506.00	-	✓
dC-CE Phosphoramidite	10-1010-02	0.25 g	3.53	3.88	3.25	-	
dC-CE Phosphoramidite	10-1010-02M	0.25 g	3.53	3.88	3.25	-	
dC-CE Phosphoramidite	10-1010-05	0.5 g	7.06	7.77	6.50	-	
dC-CE Phosphoramidite	10-1010-05M	0.5 g	7.06	7.77	6.50	-	
dC-CE Phosphoramidite	10-1010-10	1 g	14.10	15.50	13.00	-	
dC-CE Phosphoramidite	10-1010-10M	1 g	14.10	15.50	13.00	-	
dC-CE Phosphoramidite	10-1010-1B	1 g	15.40	16.90	14.20	-	
dC-CE Phosphoramidite	10-1010-1C	1 g	14.10	15.50	13.00	-	
dC-CE Phosphoramidite	10-1010-1P	1 g	15.40	16.90	14.20	-	
dC-CE Phosphoramidite	10-1010-15	10 g	141.00	155.00	130.00	-	
dC-CE Phosphoramidite	10-1010-20	2 g	28.20	31.00	25.90	-	
dC-CE Phosphoramidite	10-1010-2C	2 g	28.20	31.00	25.90	-	
dC-CE Phosphoramidite	10-1010-40	4 g	56.50	62.20	52.00	-	
dC-CE Phosphoramidite	10-1010-50	5 g	70.60	77.70	65.00	-	
dC-CE Phosphoramidite	10-1010-55	5 g	70.60	77.70	65.00	-	
dC-CE Phosphoramidite	10-1010-B2	0.25 g	3.90	4.29	3.59	-	
dC-CE Phosphoramidite	10-1010-B5	0.5 g	7.70	8.47	7.08	-	
dC-CE Phosphoramidite	10-1010-C2	0.25 g	3.53	3.88	3.25	-	
dC-CE Phosphoramidite	10-1010-C5	0.5 g	7.06	7.77	6.50	-	
dC-CE Phosphoramidite	10-1010-P4	0.4 g	6.20	6.82	5.70	-	
dC-CPG 1000	20-2011-01	0.1 g	5.81	6.39	5.35	-	
dC-CPG 1000	20-2011-02	0.25 g	12.90	14.20	11.90	-	
dC-CPG 1000	20-2011-10	1 g	48.40	53.20	44.50	-	
dC-CPG 1000	20-2111-13	1 x 10 µmol	80.60	88.70	74.20	-	
dC-CPG 1000	20-2111-41	4 x 1 µmol	28.20	31.00	25.90	-	
dC-CPG 1000	20-2111-42	4 x 0.2 µmol	28.20	31.00	25.90	-	
dC-CPG 1000	20-2111-45	4 x 40 nmol	28.20	31.00	25.90	-	
dC-CPG 1000	20-2211-14	1 x 15 µmol	121.00	133.00	111.00	-	
dC-CPG 1000	20-2211-41	4 x 1 µmol	20.20	22.20	18.60	-	
dC-CPG 1000	20-2211-42	4 x 0.2 µmol	14.10	15.50	13.00	-	
dC-CPG 1000	20-2211-45	4 x 40 nmol	14.10	15.50	13.00	-	
dC-CPG 1000, Expedite	20-2211-42E	4 x 0.2 µmol	14.10	15.50	13.00	-	
dC-CPG 2000	20-2012-01	0.1 g	15.10	16.60	13.90	-	
dC-CPG 2000	20-2012-02	0.25 g	30.20	33.20	27.80	-	
dC-CPG 2000	20-2012-10	1 g	106.00	117.00	98.00	-	
dC-CPG 2000	20-2112-42	4 x 0.2 µmol	28.20	31.00	25.90	-	
dC-CPG 2000	20-2212-42	4 x 0.2 µmol	14.10	15.50	13.00	-	
dC-CPG 500	20-2010-01	0.1 g	5.81	6.39	5.35	-	
dC-CPG 500	20-2010-02	0.25 g	12.90	14.20	11.90	-	
dC-CPG 500	20-2010-10	1 g	48.40	53.20	44.50	-	
dC-CPG 500	20-2110-13	1 x 10 µmol	80.60	88.70	74.20	-	
dC-CPG 500	20-2110-41	4 x 1 µmol	28.20	31.00	25.90	-	
dC-CPG 500	20-2110-42	4 x 0.2 µmol	28.20	31.00	25.90	-	
dC-CPG 500	20-2210-14	1 x 15 µmol	121.00	133.00	111.00	-	
dC-CPG 500	20-2210-41	4 x 1 µmol	20.20	22.20	18.60	-	
dC-CPG 500	20-2210-42	4 x 0.2 µmol	14.10	15.50	13.00	-	
dC-CPG 500	20-2510-P2	10 x 0.2 µmol	106.00	117.00	98.00	-	
dC-High Load-CPG	25-2010-02	0.25 g	16.10	17.70	14.80	-	
dC-High Load-CPG	25-2010-10	1 g	58.10	63.90	53.50	-	
dC-High Load-CPG	25-2110-17	1 x 25 µmol	101.00	111.00	93.00	-	
dC-High Load-CPG	25-2110-46	4 x 2.5 µmol	60.50	66.60	55.70	-	
dC-High Load-CPG	25-2210-18	1 x 35 µmol	149.00	164.00	137.00	-	
dC-High Load-CPG	25-2210-46	4 x 2.5 µmol	60.50	66.60	55.70	-	
dC-Thiophosphoramidite	10-1710-90	100 µmol	229.00	252.00	211.00	-	✓
dDs-CE Phosphoramidite	10-1521-02	0.25 g	601.00	661.00	553.00	-	
dDs-CE Phosphoramidite	10-1521-90	100 µmol	208.00	229.00	191.00	-	
dG LV-PS	26-2120-42	4 x 0.2 µmol	38.70	42.60	35.60	-	
dG LV-PS	26-2120-45	4 x 40 nmol	38.70	42.60	35.60	-	
dG-CE Phosphoramidite	10-1020-02	0.25 g	3.53	3.88	3.25	-	
dG-CE Phosphoramidite	10-1020-05	0.5 g	7.06	7.77	6.50	-	
dG-CE Phosphoramidite	10-1020-05M	0.5 g	7.06	7.77	6.50	-	
dG-CE Phosphoramidite	10-1020-10	1 g	14.10	15.50	13.00	-	
dG-CE Phosphoramidite	10-1020-10M	1 g	14.10	15.50	13.00	-	
dG-CE Phosphoramidite	10-1020-1B	1 g	15.40	16.90	14.20	-	
dG-CE Phosphoramidite	10-1020-1C	1 g	14.10	15.50	13.00	-	
dG-CE Phosphoramidite	10-1020-1P	1 g	15.40	16.90	14.20	-	
dG-CE Phosphoramidite	10-1020-15	10 g	141.00	155.00	130.00	-	
dG-CE Phosphoramidite	10-1020-20	2 g	28.20	31.00	25.90	-	
dG-CE Phosphoramidite	10-1020-2C	2 g	28.20	31.00	25.90	-	
dG-CE Phosphoramidite	10-1020-40	4 g	56.50	62.20	52.00	-	
dG-CE Phosphoramidite	10-1020-50	5 g	70.60	77.70	65.00	-	
dG-CE Phosphoramidite	10-1020-55	5 g	70.60	77.70	65.00	-	
dG-CE Phosphoramidite	10-1020-B2	0.25 g	3.90	4.29	3.59	-	
dG-CE Phosphoramidite	10-1020-B5	0.5 g	7.70	8.47	7.08	-	
dG-CE Phosphoramidite	10-1020-C2	0.25 g	3.53	3.88	3.25	-	
dG-CE Phosphoramidite	10-1020-P4	0.4 g	6.20	6.82	5.70	-	
dG-CE Phosphoramidite,	10-1020-02M	0.25 g	3.53	3.88	3.25	-	
dG-CE Phosphoramidite, Expedite	10-1020-C5	0.5 g	7.06	7.77	6.50	-	
dG-CPG 1000	20-2021-01	0.1 g	5.81	6.39	5.35	-	
dG-CPG 1000	20-2021-02	0.25 g	12.90	14.20	11.90	-	
dG-CPG 1000	20-2021-10	1 g	48.40	53.20	44.50	-	
dG-CPG 1000	20-2021-61	48 x 1 µmol	242.00	266.00	223.00	-	
dG-CPG 1000	20-2021-62	200 x 0.2 µmol	605.00	666.00	557.00	-	
dG-CPG 1000	20-2121-13	1 x 10 µmol	80.60	88.70	74.20	-	
dG-CPG 1000	20-2121-41	4 x 1 µmol	28.20	31.00	25.90	-	
dG-CPG 1000	20-2121-42	4 x 0.2 µmol	28.20	31.00	25.90	-	
dG-CPG 1000	20-2121-45	4 x 40 nmol	28.20	31.00	25.90	-	
dG-CPG 1000	20-2221-14	1 x 15 µmol	121.00	133.00	111.00	-	
dG-CPG 1000	20-2221-41	4 x 1 µmol	20.20	22.20	18.60	-	

Description	Reference	Scale/Quantity	Price List 2021				Ice Shipping
			EUR	CHF	GBP	USD	
dG-CPG 1000	20-2221-42	4 x 0.2 µmol	14.10	15.50	13.00	-	
dG-CPG 1000	20-2221-45	4 x 40 nmol	14.10	15.50	13.00	-	
dG-CPG 1000, Expedite	20-2221-42E	4 x 0.2 µmol	14.10	15.50	13.00	-	
dG-CPG 2000	20-2022-01	0.1 g	15.10	16.60	13.90	-	
dG-CPG 2000	20-2022-02	0.25 g	30.20	33.20	27.80	-	
dG-CPG 2000	20-2022-10	1 g	106.00	117.00	98.00	-	
dG-CPG 2000	20-2122-42	4 x 0.2 µmol	28.20	31.00	25.90	-	
dG-CPG 2000	20-2222-42	4 x 0.2 µmol	14.10	15.50	13.00	-	
dG-CPG 500	20-2020-01	0.1 g	5.81	6.39	5.35	-	
dG-CPG 500	20-2020-02	0.25 g	12.90	14.20	11.90	-	
dG-CPG 500	20-2020-10	1 g	48.40	53.20	44.50	-	
dG-CPG 500	20-2120-13	1 x 10 µmol	80.60	88.70	74.20	-	
dG-CPG 500	20-2120-41	4 x 1 µmol	28.20	31.00	25.90	-	
dG-CPG 500	20-2120-42	4 x 0.2 µmol	28.20	31.00	25.90	-	
dG-CPG 500	20-2220-14	1 x 15 µmol	121.00	133.00	111.00	-	
dG-CPG 500	20-2220-41	4 x 1 µmol	20.20	22.20	18.60	-	
dG-CPG 500	20-2220-42	4 x 0.2 µmol	14.10	15.50	13.00	-	
dG-CPG 500	20-2520-P2	10 x 0.2 µmol	106.00	117.00	98.00	-	
dG-High Load CPG	25-2020-02	0.25 g	16.10	17.70	14.80	-	
dG-High Load CPG	25-2020-10	1 g	58.10	63.90	53.50	-	
dG-High Load CPG	25-2120-17	1 x 25 µmol	101.00	111.00	93.00	-	
dG-High Load CPG	25-2120-46	4 x 2.5 µmol	60.50	66.60	55.70	-	
dG-High Load CPG	25-2120-46M	4 x 2.5 µmol	60.50	66.60	55.70	-	
dG-High Load CPG	25-2220-18	1 x 35 µmol	149.00	164.00	137.00	-	
dG-High Load CPG	25-2220-46	4 x 2.5 µmol	60.50	66.60	55.70	-	
dG-PACE Phosphoramidite	10-1160-02	0.25 g	178.00	196.00	164.00	-	
dG-PACE Phosphoramidite	10-1160-05	0.5 g	357.00	393.00	328.00	-	
dG-PACE Phosphoramidite	10-1160-10	1 g	713.00	784.00	656.00	-	
dG-Thiophosphoramidite	10-1720-90	100 µmol	229.00	252.00	211.00	-	
dG-Thiophosphoramidite,	10-1720-02	0.25 g	550.00	605.00	506.00	-	
dmf-dG-CE Phosphoramidite	10-1029-05M	0.5 g	7.06	7.77	6.50	-	
dmf-dG-CE Phosphoramidite	10-1029-10M	1 g	14.10	15.50	13.00	-	
dmf-dG-CE Phosphoramidite	10-1029-2C	2 g	28.20	31.00	25.90	-	
dmf-dG-CE Phosphoramidite	10-1029-40	4 g	56.50	62.20	52.00	-	
dmf-dG-CE Phosphoramidite,	10-1029-02M	0.25 g	3.53	3.88	3.25	-	
dmf-dG-CE Phosphoramidite, Expedite	10-1029-05E	0.5 g	7.06	7.77	6.50	-	
dmf-dG-CPG	20-2129-62	200 x 200 nm	524.00	576.00	482.00	-	
dmf-dG-CPG	20-2129-65	200 x 40 nmol	484.00	532.00	445.00	-	
dPa-CE Phosphoramidite	10-1523-90	100 µmol	186.00	205.00	171.00	-	
dPa-CE Phosphoramidite,	10-1523-02	0.25 g	601.00	661.00	553.00	-	
dT LV-PS	26-2130-42	4 x 0.2 µmol	38.70	42.60	35.60	-	
dT LV-PS	26-2130-45	4 x 40 nmol	38.70	42.60	35.60	-	
dT-CE Phosphoramidite	10-1030-02	0.25 g	3.53	3.88	3.25	-	
dT-CE Phosphoramidite	10-1030-05	0.5 g	7.06	7.77	6.50	-	
dT-CE Phosphoramidite	10-1030-05M	0.5 g	7.06	7.77	6.50	-	
dT-CE Phosphoramidite	10-1030-10	1 g	14.10	15.50	13.00	-	
dT-CE Phosphoramidite	10-1030-10M	1 g	14.10	15.50	13.00	-	
dT-CE Phosphoramidite	10-1030-1B	1 g	15.40	16.90	14.20	-	
dT-CE Phosphoramidite	10-1030-1C	1 g	14.10	15.50	13.00	-	
dT-CE Phosphoramidite	10-1030-1P	1 g	15.40	16.90	14.20	-	
dT-CE Phosphoramidite	10-1030-1S	10 g	141.00	155.00	130.00	-	
dT-CE Phosphoramidite	10-1030-20	2 g	28.20	31.00	25.90	-	
dT-CE Phosphoramidite	10-1030-2C	2 g	28.20	31.00	25.90	-	
dT-CE Phosphoramidite	10-1030-40	4 g	56.50	62.20	52.00	-	
dT-CE Phosphoramidite	10-1030-50	5 g	70.60	77.70	65.00	-	
dT-CE Phosphoramidite	10-1030-5S	5 g	70.60	77.70	65.00	-	
dT-CE Phosphoramidite	10-1030-B2	0.25 g	3.90	4.29	3.59	-	
dT-CE Phosphoramidite	10-1030-B5	0.5 g	7.70	8.47	7.08	-	
dT-CE Phosphoramidite	10-1030-C2	0.25 g	3.53	3.88	3.25	-	
dT-CE Phosphoramidite	10-1030-C5	0.5 g	7.06	7.77	6.50	-	
dT-CE Phosphoramidite	10-1030-P4	0.4 g	6.20	6.82	5.70	-	
dT-CE Phosphoramidite,	10-1030-02M	0.25 g	3.53	3.88	3.25	-	
dT-CE Phosphoramidite, Expedite	10-1030-C7	0.75 g	22.50	24.80	20.70	-	
dT-CPG 1000	20-2031-01	0.1 g	5.81	6.39	5.35	-	
dT-CPG 1000	20-2031-02	0.25 g	12.90	14.20	11.90	-	
dT-CPG 1000	20-2031-10	1 g	48.40	53.20	44.50	-	
dT-CPG 1000	20-2031-61	48 x 1 µmol	242.00	266.00	223.00	-	
dT-CPG 1000	20-2031-62	200 x 0.2 µmol	605.00	666.00	557.00	-	
dT-CPG 1000	20-2131-13	1 x 10 µmol	80.60	88.70	74.20	-	
dT-CPG 1000	20-2131-41	4 x 1 µmol	28.20	31.00	25.90	-	
dT-CPG 1000	20-2131-42	4 x 0.2 µmol	28.20	31.00	25.90	-	
dT-CPG 1000	20-2131-45	4 x 40 nmol	28.20	31.00	25.90	-	
dT-CPG 1000	20-2131-62	200 x 200 nm	524.00	576.00	482.00	-	
dT-CPG 1000	20-2131-65	200 x 40 nmol	484.00	532.00	445.00	-	
dT-CPG 1000	20-2231-14	1 x 15 µmol	121.00	133.00	111.00	-	
dT-CPG 1000	20-2231-41	4 x 1 µmol	20.20	22.20	18.60	-	
dT-CPG 1000	20-2231-42	4 x 0.2 µmol	14.10	15.50	13.00	-	
dT-CPG 1000	20-2231-45	4 x 40 nmol	14.10	15.50	13.00	-	
dT-CPG 1000, Expedite	20-2231-42E	4 x 0.2 µmol	14.10	15.50	13.00	-	
dT-CPG 2000	20-2032-01	0.1 g	15.10	16.60	13.90	-	
dT-CPG 2000	20-2032-02	0.25 g	30.20	33.20	27.80	-	
dT-CPG 2000	20-2032-10	1 g	106.00	117.00	98.00	-	
dT-CPG 2000	20-2132-42	4 x 0.2 µmol	28.20	31.00	25.90	-	
dT-CPG 2000	20-2232-42	4 x 0.2 µmol	14.10	15.50	13.00	-	
dT-CPG 500	20-2030-01	0.1 g	5.81	6.39	5.35	-	
dT-CPG 500	20-2030-02	0.25 g	12.90	14.20	11.90	-	
dT-CPG 500	20-2030-10	1 g	48.40	53.20	44.50	-	
dT-CPG 500	20-2130-13	1 x 10 µmol	80.60	88.70	74.20	-	
dT-CPG 500	20-2130-41	4 x 1 µmol	28.20	31.00	25.90	-	
dT-CPG 500	20-2130-42	4 x 0.2 µmol	28.20	31.00	25.90	-	
dT-CPG 500	20-2230-14	1 x 15 µmol	121.00	133.00	111.00	-	

Description	Reference	Scale/Quantity	Price List 2021				Ice Shipping
			EUR	CHF	GBP	USD	
dT-CPG 500	20-2230-41	4 x 1 µmol	20.20	22.20	18.60	-	
dT-CPG 500	20-2230-42	4 x 0.2 µmol	14.10	15.50	13.00	-	
dT-CPG 500	20-2530-P2	10 x 0.2 µmol	106.00	117.00	98.00	-	
dT-High Load-CPG	25-2030-02	0.25 g	16.10	17.70	14.80	-	
dT-High Load-CPG	25-2030-10	1 g	58.10	63.90	53.50	-	
dT-High Load-CPG	25-2130-17	1 x 25 µmol	101.00	111.00	93.00	-	
dT-High Load-CPG	25-2130-46	4 x 2.5 µmol	60.50	66.60	55.70	-	
dT-High Load-CPG	25-2230-18	1 x 35 µmol	149.00	164.00	137.00	-	
dT-High Load-CPG	25-2230-46	4 x 2.5 µmol	60.50	66.60	55.70	-	
dT-PACE Phosphoramidite	10-1170-02	0.25 g	178.00	196.00	164.00	-	
dT-PACE Phosphoramidite	10-1170-05	0.5 g	357.00	393.00	328.00	-	
dT-PACE Phosphoramidite	10-1170-10	1 g	713.00	784.00	656.00	-	
dT-PS	26-2630-62	200 x 200 nmol	650.00	715.00	598.00	-	
dT-PS	26-2630-65	200 x 40 nmol	650.00	715.00	598.00	-	
dT-Q-CPG 500	21-2230-42	4 x 0.2 µmol	32.30	35.50	29.70	-	
dT-Thiophosphoramidite	10-1730-90	100 µmol	229.00	252.00	211.00	-	
dT-Thiophosphoramidite,	10-1730-02	0.25 g	550.00	605.00	506.00	-	
Empty MerMade Columns - 50 nmol	20-0050-05	Pk / 48	161.00	177.00	148.00	-	
Empty Synthesis Columns - TWIST 10um/15um	20-0040-00	Pk / 10	242.00	266.00	223.00	-	
Empty Synthesis Columns, 1µmol Expedite Style	20-0021-01	Pk / 10	38.70	42.60	35.60	-	
Empty Synthesis Columns, 40 nmol, 0.2 µmol Expedite Style	20-0021-02	Pk / 10	38.70	42.60	35.60	-	
Empty Synthesis Columns-TWIST 40nm, 0.2um or 1um	20-0030-00	Pk / 10	48.40	53.20	44.50	-	
F-DMT-dA-CE Phosphoramidite	10-1400-02	0.25 g	200.00	220.00	184.00	-	
F-DMT-dA-CE Phosphoramidite	10-1400-90	100 µmol	107.00	118.00	98.00	-	
F-DMT-dC-CE Phosphoramidite	10-1410-02	0.25 g	200.00	220.00	184.00	-	
F-DMT-dC-CE Phosphoramidite	10-1410-90	100 µmol	107.00	118.00	98.00	-	
F-DMT-dG-CE Phosphoramidite	10-1420-02	0.25 g	200.00	220.00	184.00	-	
F-DMT-dG-CE Phosphoramidite	10-1420-90	100 µmol	107.00	118.00	98.00	-	
F-DMT-dT-CE Phosphoramidite	10-1430-02	0.25 g	200.00	220.00	184.00	-	
F-DMT-dT-CE Phosphoramidite	10-1430-90	100 µmol	107.00	118.00	98.00	-	
GAA Trimer Phosphoramidite	13-1200-90	100 µmol	1,114.00	1,225.00	1,025.00	-	
GAA Trimer Phosphoramidite	13-1200-95	50 µmol	557.00	613.00	512.00	-	
GAC Trimer Phosphoramidite	13-1201-90	100 µmol	1,114.00	1,225.00	1,025.00	-	
GAC Trimer Phosphoramidite	13-1201-95	50 µmol	557.00	613.00	512.00	-	
GCT Trimer Phosphoramidite	13-1213-90	100 µmol	1,114.00	1,225.00	1,025.00	-	
GCT Trimer Phosphoramidite	13-1213-95	50 µmol	557.00	613.00	512.00	-	
GGT Trimer Phosphoramidite	13-1223-90	100 µmol	1,114.00	1,225.00	1,025.00	-	
GGT Trimer Phosphoramidite	13-1223-95	50 µmol	557.00	613.00	512.00	-	
Glen Gel-Pak™ 2.5 Desalting Column	61-5025-05	Pk / 5	78.60	86.50	72.30	-	
Glen Gel-Pak™ 2.5 Desalting Column	61-5025-25	Pk / 25	393.00	432.00	362.00	-	
Glen UnySupport PS	26-5140-52	10 x 200 nmol	80.60	88.70	74.20	-	
Glen UnySupport PS	26-5140-55	10 x 40 nmol	80.60	88.70	74.20	-	
Glen UnySupport™ 1000	20-5041-01	0.1 g	11.10	12.20	10.20	-	
Glen UnySupport™ 1000	20-5041-02	0.25 g	25.20	27.70	23.20	-	
Glen UnySupport™ 1000	20-5041-10	1 g	95.80	105.40	88.10	-	
Glen UnySupport™ 1000	20-5141-41	4 x 1 µmol	48.40	53.20	44.50	-	
Glen UnySupport™ 1000	20-5241-42	4 x 0.2 µmol	32.30	35.50	29.70	-	
Glen UnySupport™ 500	20-5040-01	0.1 g	11.10	12.20	10.20	-	
Glen UnySupport™ 500	20-5040-02	0.25 g	25.20	27.70	23.20	-	
Glen UnySupport™ 500	20-5040-10	1 g	95.80	105.40	88.10	-	
Glen UnySupport™ 500	20-5140-13	1 x 10 µmol	88.00	96.80	81.00	-	
Glen UnySupport™ 500	20-5140-41	4 x 1 µmol	53.00	58.30	48.80	-	
Glen UnySupport™ 500	20-5140-42	4 x 0.2 µmol	35.20	38.70	32.40	-	
Glen UnySupport™ 500	20-5140-45	4 x 40 nmol	35.20	38.70	32.40	-	
Glen UnySupport™ 500	20-5140-91	96 x 1 µmol	578.00	636.00	532.00	-	
Glen UnySupport™ 500	20-5140-92	96 x 0.2 µmol	385.00	424.00	354.00	-	
Glen UnySupport™ 500	20-5140-95	96 x 40 nmol	385.00	424.00	354.00	-	
Glen UnySupport™ 500	20-5240-14	1 x 15 µmol	132.00	145.00	121.00	-	
Glen UnySupport™ 500	20-5240-42	4 x 0.2 µmol	35.20	38.70	32.40	-	
Glen UnySupport™ 500	20-5240-45	4 x 40 nmol	35.20	38.70	32.40	-	
Glen UnySupport™ FC	22-5141-41	4 x 1.0 µmol	50.00	55.00	46.00	-	
Glen UnySupport™ FC	22-5141-55	10 x 40 nmol	86.00	94.60	79.10	-	
Glen UnySupport™ FC	22-5141-95	96 x 40 nmol	376.00	414.00	346.00	-	
GTT Trimer Phosphoramidite	13-1233-90	100 µmol	1,114.00	1,225.00	1,025.00	-	
GTT Trimer Phosphoramidite	13-1233-95	50 µmol	557.00	613.00	512.00	-	
I-CE Phosphoramidite	10-3040-90	100 µmol	249.00	274.00	229.00	-	
I-CE Phosphoramidite	10-3040-95	50 µmol	124.00	136.00	114.00	-	
N4-Ac-N4-Et-dC-CE Phosphoramidite	10-1513-90	100 µmol	295.00	325.00	271.00	-	
Pac-ds-CE Phosphoramidite	10-1522-90	100 µmol	243.00	267.00	224.00	-	
Pac-ds-CE Phosphoramidite,	10-1522-02	0.25 g	601.00	661.00	553.00	-	
pdU-CE Phosphoramidite	10-1054-90	100 µmol	85.20	93.70	78.40	-	
TAC Trimer Phosphoramidite	13-1301-90	100 µmol	1,114.00	1,225.00	1,025.00	-	
TAC Trimer Phosphoramidite	13-1301-95	50 µmol	557.00	613.00	512.00	-	
TCT Trimer Phosphoramidite	13-1313-90	100 µmol	1,114.00	1,225.00	1,025.00	-	
TCT Trimer Phosphoramidite	13-1313-95	50 µmol	557.00	613.00	512.00	-	
TGC Trimer Phosphoramidite	13-1321-90	100 µmol	1,114.00	1,225.00	1,025.00	-	
TGC Trimer Phosphoramidite	13-1321-95	50 µmol	557.00	613.00	512.00	-	
TGG Trimer Phosphoramidite	13-1322-90	100 µmol	1,114.00	1,225.00	1,025.00	-	
TGG Trimer Phosphoramidite	13-1322-95	50 µmol	557.00	613.00	512.00	-	
Trimer Phosphoramidite Mix 1	13-1991-90	100 µmol	1,638.00	1,802.00	1,507.00	-	
Trimer Phosphoramidite Mix 1	13-1991-95	50 µmol	819.00	901.00	753.00	-	
TTC Trimer Phosphoramidite	13-1331-90	100 µmol	1,114.00	1,225.00	1,025.00	-	
TTC Trimer Phosphoramidite	13-1331-95	50 µmol	557.00	613.00	512.00	-	
Universal Support II	20-5110-91	96 x 1 µmol	578.00	636.00	532.00	-	
Universal Support III	26-5010-10	1 g	164.00	180.00	151.00	-	
UnyLinker™ CPG 500	20-5030-01	0.1 g	12.10	13.30	11.10	-	
UnyLinker™ CPG 500	20-5030-02	0.25 g	27.50	30.30	25.30	-	
UnyLinker™ CPG 500	20-5030-10	1 g	105.00	116.00	97.00	-	
UnyLinker™ CPG 500	20-5130-13	1 x 10 µmol	88.00	96.80	81.00	-	
UnyLinker™ CPG 500	20-5130-41	4 x 1 µmol	53.00	58.30	48.80	-	
UnyLinker™ CPG 500	20-5130-42	4 x 0.2 µmol	35.20	38.70	32.40	-	

Description	Reference	Scale/Quantity	Price List 2021				Ice Shipping
			EUR	CHF	GBP	USD	
UnyLinker™ CPG 500	20-5130-45	4 x 40 nmol	35.20	38.70	32.40	-	
UnyLinker™ CPG 500	20-5230-14	1 x 15 μmol	132.00	145.00	121.00	-	
UnyLinker™ CPG 500	20-5230-41	4 x 1 μmol	53.00	58.30	48.80	-	
UnyLinker™ CPG 500	20-5230-42	4 x 0.2 μmol	35.20	38.70	32.40	-	
UnyLinker™ CPG 500	20-5230-45	4 x 40 nmol	35.20	38.70	32.40	-	
<b>Glen Solvant</b>							<b>EA1B1Z</b>
0.02M Iodine in Tetrahydrofuran/ Water/ Pyridine	40-4132-52	200 mL	28.20	31.00	25.90	-	
0.02M Iodine in Tetrahydrofuran/ Water/ Pyridine	40-4132-55	240 mL	37.00	40.70	34.00	-	
0.02M Iodine in Tetrahydrofuran/ Water/ Pyridine	40-4132-57	450 mL	67.70	74.50	62.30	-	
0.02M Iodine in Tetrahydrofuran/ Water/ Pyridine	40-4132-62	2000 mL	229.00	252.00	211.00	-	
0.02M Iodine in Tetrahydrofuran/ Water/ Pyridine	40-4132-66	60 mL	14.10	15.50	13.00	-	
0.02M Iodine in Tetrahydrofuran/Pyridine/Water	40-4330-52	200 mL	21.20	23.30	19.50	-	
0.02M Iodine in Tetrahydrofuran/Pyridine/Water	40-4330-57	450 mL	50.80	55.90	46.70	-	
0.02M Iodine in Tetrahydrofuran/Pyridine/Water	40-4330-61	960 mL	109.00	120.00	100.00	-	
0.02M Iodine in Tetrahydrofuran/Pyridine/Water	40-4330-62	2000 mL	229.00	252.00	211.00	-	
0.05 M Potassium Carbonate in Methanol	60-4600-30	30 mL	21.20	23.30	19.50	-	
0.05 M Sulfurizing Reagent II in pyridine/acetonitrile, Expedite	40-4137-52E	200 ml	141.00	155.00	130.00	-	
0.05M I2 in Pyridine/H2O	40-4035-71	1000 mL	159.00	175.00	146.00	-	
0.10 M CSO in Anhydrous Acetonitrile, Expedite	40-4631-52E	200 ml	60.00	66.00	55.20	-	
0.25 M 5-Benzylthio-1H-Tetrazole in Acetonitrile	30-3172-52	200 mL	70.60	77.70	65.00	-	
0.25 M 5-Benzylthio-1H-tetrazole in Acetonitrile	30-3170-45	45 mL	28.20	31.00	25.90	-	
0.25 M 5-Benzylthio-1H-tetrazole in Acetonitrile	30-3170-52	200 mL	70.60	77.70	65.00	-	
0.25 M 5-Benzylthio-1H-tetrazole in Acetonitrile	30-3170-57	450 mL	141.00	155.00	130.00	-	
0.25 M 5-Benzylthio-1H-tetrazole in Acetonitrile	30-3170-62	2000 mL	536.00	590.00	493.00	-	
0.25 M 5-Ethylthio-1H-Tetrazole in Anhydrous Acetonitrile	30-3140-45	45 mL	28.20	31.00	25.90	-	
0.25 M 5-Ethylthio-1H-Tetrazole in Anhydrous Acetonitrile	30-3140-52	200 mL	70.60	77.70	65.00	-	
0.25 M 5-Ethylthio-1H-Tetrazole in Anhydrous Acetonitrile	30-3140-57	450 mL	141.00	155.00	130.00	-	
0.25 M 5-Ethylthio-1H-Tetrazole in Anhydrous Acetonitrile	30-3140-61	960 mL	258.00	284.00	237.00	-	
0.25 M 5-Ethylthio-1H-Tetrazole in Anhydrous Acetonitrile	30-3140-62	2000 mL	536.00	590.00	493.00	-	
0.25 M 5-Ethylthio-1H-Tetrazole in Anhydrous Acetonitrile	30-3142-52	200 mL	70.60	77.70	65.00	-	
0.25 M DCI in Anhydrous Acetonitrile	30-3150-45	45 mL	28.20	31.00	25.90	-	
0.25 M DCI in Anhydrous Acetonitrile	30-3150-52	200 mL	70.60	77.70	65.00	-	
0.25 M DCI in Anhydrous Acetonitrile	30-3150-57	450 mL	141.00	155.00	130.00	-	
0.25 M DCI in Anhydrous Acetonitrile	30-3150-61	960 mL	281.00	309.00	259.00	-	
0.25 M DCI in Anhydrous Acetonitrile	30-3150-62	2000 mL	536.00	590.00	493.00	-	
0.25 M DCI in Anhydrous Acetonitrile	30-3152-52	200 mL	70.60	77.70	65.00	-	
0.25M 5-Benzylthio-1H-Tetrazole in Acetonitrile, Expedite	30-3172-52E	200 mL	70.60	77.70	65.00	-	
0.2M Saccharin 1-Methylimidazole (SMI) in Acetonitrile	30-3180-45	45 mL	61.50	67.70	56.60	-	
0.40M Tetrazole in Acetonitrile	30-3105-71	1000 mL	268.00	295.00	247.00	-	
0.5 M CSO in Anhydrous Acetonitrile	40-4632-62	2 L	1,764.00	1,940.00	1,623.00	-	
0.5 M CSO in Anhydrous Acetonitrile (Expedite)	40-4632-52E	200 mL	176.00	194.00	162.00	-	
0.5M CSO in Anhydrous Acetonitrile	40-4632-52	200 mL	176.00	194.00	162.00	-	
10% 1-Methylimidazole in Tetrahydrofuran	40-4120-52	200 mL	21.20	23.30	19.50	-	
10% 1-Methylimidazole in Tetrahydrofuran	40-4120-57	450 mL	50.80	55.90	46.70	-	
10% 1-Methylimidazole in Tetrahydrofuran	40-4120-62	2000 mL	229.00	252.00	211.00	-	
10% 1-Methylimidazole in Tetrahydrofuran/Pyridine	40-4122-51	150 mL	23.10	25.40	21.30	-	
10% 1-Methylimidazole in Tetrahydrofuran/Pyridine	40-4122-52	200 mL	28.20	31.00	25.90	-	
10% 1-Methylimidazole in Tetrahydrofuran/Pyridine	40-4122-57	450 mL	67.70	74.50	62.30	-	
10% 1-Methylimidazole in Tetrahydrofuran/Pyridine	40-4122-62	2000 mL	229.00	252.00	211.00	-	
10% 1-Methylimidazole in Tetrahydrofuran/Pyridine	40-4122-66	60 mL	14.10	15.50	13.00	-	
16% 1-Methylimidazole in Tetrahydrofuran	40-4220-45	45 mL	14.10	15.50	13.00	-	
16% 1-Methylimidazole in Tetrahydrofuran	40-4220-52	200 mL	28.20	31.00	25.90	-	
16% 1-Methylimidazole in Tetrahydrofuran	40-4220-57	450 mL	67.70	74.50	62.30	-	
16% 1-Methylimidazole in Tetrahydrofuran	40-4220-61	960 mL	144.00	158.00	132.00	-	
16% 1-Methylimidazole in Tetrahydrofuran	40-4220-62	2000 mL	300.00	330.00	276.00	-	
2% Aqueous Trifluoroacetic Acid	60-4040-57	450 mL	29.00	31.90	26.70	-	
2.0M Triethylamine Acetate, HPLC grade, pH=7	60-4110-52	200 mL	48.40	53.20	44.50	-	
2.0M Triethylamine Acetate, HPLC grade, pH=7	60-4110-57	450 mL	96.80	106.50	89.10	-	
2.0M Triethylamine Acetate, HPLC grade, pH=7	60-4110-60	960 mL	161.00	177.00	148.00	-	
2.0M Triethylamine Acetate, HPLC grade, pH=7	60-4110-62	2000 mL	323.00	355.00	297.00	-	
2.5% Dichloroacetic acid in Dichloromethane	40-4042-57	450 mL	25.40	27.90	23.40	-	
2.5% Dichloroacetic acid in Dichloromethane	40-4042-62	2000 mL	102.00	112.00	94.00	-	
3% TCA/DCM	40-4140-68	180 mL	12.70	14.00	11.70	-	
3% TCA/DCM, Expedite	40-4140-61E	960 mL	52.90	58.20	48.70	-	
3% Aqueous Trifluoroacetic Acid	60-4041-57	450 mL	31.70	34.90	29.20	-	
3% DCA in Toluene	40-4240-71	1000 mL	102.00	112.00	94.00	-	
3% Dichloroacetic acid in Dichloromethane	40-4040-57	450 mL	25.40	27.90	23.40	-	
3% Dichloroacetic acid in Dichloromethane	40-4040-61	960 mL	52.90	58.20	48.70	-	
3% Dichloroacetic acid in Dichloromethane	40-4040-62	2000 mL	102.00	112.00	94.00	-	
3% Dichloroacetic acid in Dichloromethane	40-4040-71	1000 mL	56.50	62.20	52.00	-	
3% TCA/DCM	40-4140-57	450 mL	25.40	27.90	23.40	-	
3% TCA/DCM	40-4140-61	960 mL	52.90	58.20	48.70	-	
3% TCA/DCM	40-4140-62	2000 mL	102.00	112.00	94.00	-	
3% TCA/DCM	40-4140-71	1000 mL	56.50	62.20	52.00	-	
4% Aqueous Trifluoroacetic Acid	60-4042-57	450 mL	31.70	34.90	29.20	-	
4,5-Dicyanoimidazole (DCI), Crystalline	30-3050-10	1 g	24.70	27.20	22.70	-	
4,5-Dicyanoimidazole (DCI), Crystalline	30-3050-25	25 g	353.00	388.00	325.00	-	
5% Phenoxyacetic anhydride in THF	40-4212-52	200 mL	141.00	155.00	130.00	-	
5% Phenoxyacetic anhydride in THF	40-4212-57	450 mL	318.00	350.00	293.00	-	
5% Phenoxyacetic anhydride in THF, Expedite	40-4212-52E	200 mL	141.00	155.00	130.00	-	
5% Phenoxyacetic anhydride in THF/Pyridine	40-4210-52	200 mL	141.00	155.00	130.00	-	
5% Phenoxyacetic anhydride in THF/Pyridine	40-4210-57	450 mL	318.00	350.00	293.00	-	
5-Benzylthio-1H-tetrazole (BTT)	30-3070-10	1 g	24.70	27.20	22.70	-	
5-Benzylthio-1H-tetrazole (BTT)	30-3070-20	2 g	42.30	46.50	38.90	-	
5-Benzylthio-1H-tetrazole (BTT)	30-3070-25	25 g	353.00	388.00	325.00	-	
5-Ethylthio-1H-Tetrazole	30-3040-10	1 g	24.70	27.20	22.70	-	
5-Ethylthio-1H-Tetrazole	30-3040-20	2 g	42.30	46.50	38.90	-	
5-Ethylthio-1H-Tetrazole	30-3040-25	25 g	353.00	388.00	325.00	-	
6.5% Dimethylaminopyridine in Tetrahydrofuran	40-4020-52	200 mL	29.60	32.60	27.20	-	
Acetonitrile, anhydrous	40-4050-45	60 mL	8.47	9.32	7.79	-	
Acetonitrile, anhydrous	40-4050-50	100 mL	11.30	12.40	10.40	-	

Description	Reference	Scale/Quantity	Price List 2021				Ice Shipping
			EUR	CHF	GBP	USD	
Acetonitrile, anhydrous	40-4050-53	300 mL	28.20	31.00	25.90	-	
Acetonitrile, anhydrous	40-4050-57	450 mL	35.30	38.80	32.50	-	
Acetonitrile/Melm	40-4015-71	1000 mL	112.00	123.00	103.00	-	
Acetonitrile/Pyridine (1:1)	40-4500-51	150 mL	21.60	23.80	19.90	-	
Acetonitrile/Pyridine (1:1)	40-4500-60	960 mL	93.00	102.30	85.60	-	
Acetonitrile/Pyridine, anhydrous (1:1)	40-4550-50	100 mL	18.50	20.40	17.00	-	
Acetonitrile/Pyridine, anhydrous (95:5)	40-4551-50	100 mL	18.50	20.40	17.00	-	
Cap Mix (40-4026-75+40-4027-75ref)	40-4028-71	500 mL	134.00	147.00	123.00	-	
Conjugation Reagent (TCEP)	50-1003-01	1 g	107.00	118.00	98.00	-	
DCM	40-4100-62	2000 mL	93.00	102.30	85.60	-	
Dichloroacetic acid (DCA)	40-4044-50	100 mL	38.50	42.40	35.40	-	
Dichloroacetic acid (DCA)	40-4044-54	120 mL	46.20	50.80	42.50	-	
Fluorous Chemical Phosphorylation Reagent II	10-1904-02	0.25 g	279.00	307.00	257.00	-	
Fluorous Chemical Phosphorylation Reagent II	10-1904-90	100 µmol	136.00	150.00	125.00	-	
Isopropyl Phosphite/Acetonitrile/Pyridine	40-4510-45	45 mL	27.70	30.50	25.50	-	
Loading Buffer,	61-4120-52		50.00	55.00	46.00	-	
Oligo Modifying Reagent	50-1001-01	150 µmol	415.00	457.00	382.00	-	
Peptide Modifying Reagent	50-1002-01	1 g	358.00	394.00	329.00	-	
Saccharin 1-Methylimidazole (SMI)	30-3080-10	1 g	23.00	25.30	21.20	-	
Sublimed 1H-Tetrazole in Anhydrous Acetonitrile	30-3100-45	45 mL	28.20	31.00	25.90	-	
Sublimed 1H-Tetrazole in Anhydrous Acetonitrile	30-3100-52	200 mL	70.60	77.70	65.00	-	
Sublimed 1H-Tetrazole in Anhydrous Acetonitrile	30-3100-57	450 mL	141.00	155.00	130.00	-	
Sublimed 1H-Tetrazole in Anhydrous Acetonitrile	30-3100-62	2000 mL	536.00	590.00	493.00	-	
Sublimed 1H-Tetrazole in Anhydrous Acetonitrile	30-3102-52	200 mL	77.00	84.70	70.80	-	
Sublimed 1H-Tetrazole in Anhydrous Acetonitrile	30-3102-66	60 mL	38.50	42.40	35.40	-	
Sublimed 1H-Tetrazole in Anhydrous Acetonitrile, Expedite	30-3102-52E	200 mL	77.00	84.70	70.80	-	
Sulfurizing Reagent II	40-4037-10	1 g	35.30	38.80	32.50	-	
Sulfurizing Reagent II	40-4037-20	2 g	70.60	77.70	65.00	-	
Sulfurizing Solution	40-4137-51	150 mL	70.60	77.70	65.00	-	
Sulfurizing Solution	40-4137-52	200 mL	141.00	155.00	130.00	-	
Sulfurizing Solution	40-4137-57	450 mL	691.00	760.00	636.00	-	
Tetrahydrofuran/ 2,6-Lutidine/ Acetic Anhydride	40-4010-52	200 mL	21.20	23.30	19.50	-	
Tetrahydrofuran/ 2,6-Lutidine/ Acetic Anhydride	40-4010-57	450 mL	50.80	55.90	46.70	-	
Tetrahydrofuran/ 2,6-Lutidine/ Acetic Anhydride	40-4010-61	960 mL	109.00	120.00	100.00	-	
Tetrahydrofuran/ 2,6-Lutidine/ Acetic Anhydride	40-4010-62	2000 mL	229.00	252.00	211.00	-	
Tetrahydrofuran/ Acetic Anhydride	40-4012-50	100 mL	15.40	16.90	14.20	-	
Tetrahydrofuran/ Acetic Anhydride	40-4012-52	200 mL	21.20	23.30	19.50	-	
Tetrahydrofuran/ Acetic Anhydride	40-4012-57	450 mL	50.80	55.90	46.70	-	
Tetrahydrofuran/ Acetic Anhydride	40-4012-62	2000 mL	194.00	213.00	178.00	-	
Tetrahydrofuran/ Acetic Anhydride	40-4012-66	60 mL	10.60	11.70	9.80	-	
Tetrahydrofuran/ Pyridine/ Acetic Anhydride	40-4110-45	45 mL	11.30	12.40	10.40	-	
Tetrahydrofuran/ Pyridine/ Acetic Anhydride	40-4110-52	200 mL	21.20	23.30	19.50	-	
Tetrahydrofuran/ Pyridine/ Acetic Anhydride	40-4110-57	450 mL	50.80	55.90	46.70	-	
Tetrahydrofuran/ Pyridine/ Acetic Anhydride	40-4110-62	2000 mL	229.00	252.00	211.00	-	
<b>Glen Purifications</b>							<b>EA1C12</b>
Adapter Rack for 96 well manifolds, Pk/1	60-0010-01		16.10	17.70	14.80	-	
Fluoro-Pak™ columns (with Luer adaptor), Pk/10	61-2100-10A		143.00	157.00	132.00	-	
Fluoro-Pak™ columns, Pk/10	61-2100-10		143.00	157.00	132.00	-	
Fluoro-Pak™ II columns (with Luer adaptor), Pk/10	61-4100-10A		193.00	212.00	178.00	-	
Fluoro-Pak™ II columns, Pk/10	61-4100-10		193.00	212.00	178.00	-	
Glen-Pak 50 mg DNA Purification Cartridge	60-5000-96	1 pack of 96 pce	335.00	369.00	308.00	-	
Glen-Pak™ DNA 30 mg 96-Well Plate	60-5400-01	Pk / 1	479.00	527.00	441.00	-	
Glen-Pak™ DNA Cartridge 3 g	60-5300-01	Pk / 1	121.00	133.00	111.00	-	
Glen-Pak™ DNA purification cartridge	60-5100-30	Pk / 30	161.00	177.00	148.00	-	
Glen-Pak™ DNA purification cartridge (for use with disposable syringes)	60-5200-01	Pk / 1	10.10	11.10	9.30	-	
Glen-Pak™ RNA purification cartridge	60-6100-30	Pk / 30	181.00	199.00	167.00	-	
Glen-Pak™ RNA purification cartridge (for use with disposable syringes)	60-6200-01	Pk / 1	12.60	13.90	11.60	-	
Glen UnySupport™ 1000	20-5241-41	4 x 1 µmol	48.40	53.20	44.50	-	
Poly-Pak Cartridges	60-1100-01	Pk / 1	6.45	7.10	5.93	-	
Poly-Pak Cartridges	60-1100-10	Pk / 10	64.50	71.00	59.30	-	
Poly-Pak II Cartridge	60-3100-01	Pk / 1	9.68	10.65	8.91	-	
Poly-Pak II Cartridge	60-3100-10	Pk / 10	96.80	106.50	89.10	-	
Poly-Pak Packing	60-1000-05	5 g	212.00	233.00	195.00	-	
Poly-Pak Packing	60-1000-25	25 g	1,058.00	1,164.00	973.00	-	
Seal for unused wells on Adapter Rack, Pk/1	60-0020-01		24.20	26.60	22.30	-	
<b>Glen Research reagents for cartridge purification</b>							<b>EA1D12</b>
1-Ethynyl-dSpacer CE Phosphoramidite	10-1910-90	100 µmol	446.00	491.00	410.00	-	
1-Me-A-CE Phosphoramidite	10-3501-90	100 µmol	524.00	576.00	482.00	-	
1-Me-dA-CE Phosphoramidite	10-1501-02	0.25 g	983.00	1,081.00	904.00	-	
1-Me-dA-CE Phosphoramidite	10-1501-90	100 µmol	328.00	361.00	302.00	-	
1-Me-dA-CE Phosphoramidite	10-1501-95	50 µmol	164.00	180.00	151.00	-	
1-Methyl-PseudoUridine Phosphoramidite	10-3056-95	50 µmol	550.00	605.00	506.00	-	
2'-F-5-Me-U-ANA-CE Phosphoramidite	10-3850-02	0.25 g	328.00	361.00	302.00	-	
2'-F-5-Me-U-ANA-CE Phosphoramidite	10-3850-05	0.5 g	655.00	721.00	603.00	-	
2'-F-G-CE Phosphoramidite, Expedite	10-3420-02E	0.25 g	98.30	108.10	90.40	-	
2,2'-Dipicolylamin Phosphoramidite	10-5801-02	250 mg	895.00	985.00	823.00	-	
2,2'-Dipicolylamin Phosphoramidite	10-5801-90	100 µmol	286.00	315.00	263.00	-	
2,2'-Dipicolylamin Phosphoramidite	10-5801-95	50 µmol	150.00	165.00	138.00	-	
2',3'-ddA-CE Phosphoramidite	10-7001-02	0.25 g	714.00	785.00	657.00	-	
2',3'-ddA-CE Phosphoramidite	10-7001-90	100 µmol	170.00	187.00	156.00	-	
2',3'-ddC-CE Phosphoramidite	10-7101-02	0.25 g	714.00	785.00	657.00	-	
2',3'-ddC-CE Phosphoramidite	10-7101-90	100 µmol	170.00	187.00	156.00	-	
2',3'-ddC-CPG	20-2017-01	0.1 g	302.00	332.00	278.00	-	
2',3'-ddC-CPG	20-2117-41	4 x 1 µmol	605.00	666.00	557.00	-	
2',3'-ddC-CPG	20-2117-42	4 x 0.2 µmol	202.00	222.00	186.00	-	
2',3'-ddC-CPG, Mermade	20-2117-42M	4 x 0.2 µmol	202.00	222.00	186.00	-	
2',3'-ddG-CE Phosphoramidite	10-7201-02	0.25 g	885.00	974.00	814.00	-	
2',3'-ddG-CE Phosphoramidite	10-7201-90	100 µmol	190.00	209.00	175.00	-	
2',3'-ddT-CE Phosphoramidite	10-7301-02	0.25 g	714.00	785.00	657.00	-	
2',3'-ddT-CE Phosphoramidite	10-7301-90	100 µmol	170.00	187.00	156.00	-	
2,6-Diaminopurine riboside a-thiotriphosphate (0.25mM) (DTPaS)	80-3305-01	0.1 mL	143.00	157.00	132.00	-	



Description	Reference	Scale/Quantity	Price List 2021				Ice Shipping
			EUR	CHF	GBP	USD	
2,6-Diaminopurine-TOM-CE Phosphoramidite	10-3085-90	100 µmol	557.00	613.00	512.00	-	
2,6-Diaminopurine-TOM-CE Phosphoramidite	10-3085-95	50 µmol	278.00	306.00	256.00	-	
2-Amino-dA-CE Phosphoramidite	10-1085-02	0.25 g	328.00	361.00	302.00	-	
2-Amino-dA-CE Phosphoramidite	10-1085-90	100 µmol	164.00	180.00	151.00	-	
2-Amino-dA-CE Phosphoramidite	10-1085-95	50 µmol	91.70	100.90	84.40	-	
2-Aminopurine riboside a-thiotriphosphate (20mM) (2-AP-TTPaS)	80-3304-01	0.1 mL	143.00	157.00	132.00	-	
2-Aminopurine-CE Phosphoramidite	10-1046-02	0.25 g	465.00	512.00	428.00	-	
2-Aminopurine-CE Phosphoramidite	10-1046-90	100 µmol	177.00	195.00	163.00	-	
2'-deoxyAdenosine a-thiotriphosphate (15mM) (2'dATTPaS)	80-1000-01	0.1 mL	143.00	157.00	132.00	-	
2'-deoxyCytidine a-thiotriphosphate (15mM) (2'-dCTTPaS)	80-1010-01	0.1 mL	143.00	157.00	132.00	-	
2'-deoxyGuanosine a-thiotriphosphate (5mM) (2'dGTPPaS)	80-1020-01	0.1 mL	143.00	157.00	132.00	-	
2'-DeoxyNebularine-CE Phosphoramidite (Purine)	10-1041-02	0.25 g	334.00	367.00	307.00	-	
2'-DeoxyNebularine-CE Phosphoramidite (Purine)	10-1041-90	100 µmol	138.00	152.00	127.00	-	
2'-DeoxyNebularine-CE Phosphoramidite (Purine), Expedite	10-1041-02E	0.25 g	334.00	367.00	307.00	-	
2'-deoxypseudoU-CE Phosphoramidite	10-1055-02	0.25 g	1,389.00	1,528.00	1,278.00	-	
2'-deoxypseudoU-CE Phosphoramidite	10-1055-90	100 µmol	465.00	512.00	428.00	-	
2'-deoxypseudoU-CE Phosphoramidite	10-1055-95	50 µmol	246.00	271.00	226.00	-	
2'-deoxyUridine a-thiotriphosphate (5mM) (2'dUTTP)	80-1040-01	0.1 mL	143.00	157.00	132.00	-	
2'-F-A-ANA-CE Phosphoramidite	10-3800-02	0.25 g	655.00	721.00	603.00	-	
2'-F-Ac-C-ANA-CE Phosphoramidite	10-3815-02	0.25 g	328.00	361.00	302.00	-	
2'-F-A-CE Phosphoramidite	10-3400-02	0.25 g	98.30	108.10	90.40	-	
2'-F-A-CE Phosphoramidite	10-3400-90	100 µmol	72.00	79.20	66.20	-	
2'-F-dI-CE Phosphoramidite	10-1082-02	0.25 g	1,278.00	1,406.00	1,176.00	-	
2'-F-dI-CE Phosphoramidite	10-1082-90	100 µmol	472.00	519.00	434.00	-	
2'-F-dI-CE Phosphoramidite	10-1082-95	50 µmol	236.00	260.00	217.00	-	
2'-F-dI-CE Phosphoramidite - Expedite	10-1082-95E	50 µmol	236.00	260.00	217.00	-	✓
2'-F-G-ANA-CE Phosphoramidite	10-3820-02	0.25 g	655.00	721.00	603.00	-	
2'-F-G-ANA-CE Phosphoramidite	10-3820-90	100 µmol	229.00	252.00	211.00	-	
2'-F-G-CE Phosphoramidite	10-3420-02	0.25 g	98.30	108.10	90.40	-	
2'-F-G-CE Phosphoramidite	10-3420-05	0.5 g	197.00	217.00	181.00	-	
2'-F-G-CE Phosphoramidite	10-3420-90	100 µmol	72.00	79.20	66.20	-	
2'-OMe-A-CE Phosphoramidite	10-3100-SP	2 g / 60 ml bottle	400.00	440.00	368.00	-	
2'-OMe-G-CE Phosphoramidite	10-3121-SP	2 g / 60 ml bottle	400.00	440.00	368.00	-	
2'-OMe-Pac-A-CE Phosphoramidite	10-3601-10	1 g	328.00	361.00	302.00	-	
2'-O-Methyl-Uridine thiotriphosphate (20mM)	80-1141-01	0.1 mL	143.00	157.00	132.00	-	
2'-OMe-U-RNA-CPG	20-3630-01	0.1 g	52.40	57.60	48.20	-	
2-Thio-dT-CE Phosphoramidite	10-1036-02	0.25 g	885.00	974.00	814.00	-	
2-Thio-dT-CE Phosphoramidite	10-1036-90	100 µmol	387.00	426.00	356.00	-	
2-Thio-dT-CE Phosphoramidite	10-1036-95	50 µmol	216.00	238.00	199.00	-	
3' - (6 - FAM)CPG -0.2 µm(1pk/4)MerMade	20-2961-42M	4 x 0.2 µmol	121.00	133.00	111.00	-	
3' - Amino-ModifierC7 CPG1000-(1pk/4)Mer	20-2958-41M	4 x 1 µmol	155.00	171.00	143.00	-	
3' - CPR II CPG	20-2903-01	0.1 g	91.70	100.90	84.40	-	
3' -Amino-ModifierC7 CPG1000-(1pk/4)Mer.	20-2958-42M	4 x 0.2 µmol	85.70	94.30	78.80	-	
3' Thiol-Modifier 6 S-S CPG	20-2938-01	0.1 g	98.30	108.10	90.40	-	
3' Thiol-Modifier 6 S-S CPG	20-2938-10	1 g	983.00	1,081.00	904.00	-	✓
3'-(6-FAM) CPG	20-2961-01	0.1 g	157.00	173.00	144.00	-	
3'-(6-FAM) CPG	20-2961-10	1 g	1,304.00	1,434.00	1,200.00	-	
3'-(6-FAM) CPG	20-2961-13	1 x 10 µmol	393.00	432.00	362.00	-	
3'-(6-FAM) CPG	20-2961-14	1 x 15 µmol	590.00	649.00	543.00	-	
3'-(6-FAM) CPG	20-2961-41	4 x 1 µmol	222.00	244.00	204.00	-	
3'-(6-FAM) CPG	20-2961-42	4 x 0.2 µmol	121.00	133.00	111.00	-	
3'-(6-FAM) PS	26-2961-01	0.1 g	170.00	187.00	156.00	-	
3'-(6-FAM) PS	26-2961-10	1 g	1,369.00	1,506.00	1,259.00	-	
3'-(6-FAM) PS	26-2961-52	10 x 200 nmol	393.00	432.00	362.00	-	
3'-(6-FAM) PS	26-2961-55	10 x 40 nmol	393.00	432.00	362.00	-	
3'-(6-Fluorescein) CPG	20-2964-01	0.1 g	157.00	173.00	144.00	-	
3'-(6-Fluorescein) CPG	20-2964-10	1 g	1,304.00	1,434.00	1,200.00	-	
3'-(6-Fluorescein) CPG	20-2964-13	1 x 10 µmol	393.00	432.00	362.00	-	
3'-(6-Fluorescein) CPG	20-2964-14	1 x 15 µmol	590.00	649.00	543.00	-	
3'-(6-Fluorescein) CPG	20-2964-41	4 x 1 µmol	262.00	288.00	241.00	-	
3'-(6-Fluorescein) CPG	20-2964-42	4 x 0.2 µmol	157.00	173.00	144.00	-	
3'-6-Fluorescein Serinol CPG	20-2994-01	0.1 g	157.00	173.00	144.00	-	
3'-6-Fluorescein Serinol CPG	20-2994-10	1 g	1,304.00	1,434.00	1,200.00	-	
3'-6-Fluorescein Serinol CPG	20-2994-13	1 x 10 µmol	393.00	432.00	362.00	-	
3'-6-Fluorescein Serinol CPG	20-2994-14	1 x 15 µmol	590.00	649.00	543.00	-	
3'-6-Fluorescein Serinol CPG	20-2994-41	4 x 1 µmol	262.00	288.00	241.00	-	
3'-6-Fluorescein Serinol CPG	20-2994-42	4 x 0.2 µmol	157.00	173.00	144.00	-	
3'-Acridine CPG	20-2973-01	0.1 g	157.00	173.00	144.00	-	
3'-Acridine CPG	20-2973-10	1 g	1,304.00	1,434.00	1,200.00	-	
3'-Acridine CPG	20-2973-13	1 x 10 µmol	393.00	432.00	362.00	-	
3'-Acridine CPG	20-2973-14	1 x 15 µmol	590.00	649.00	543.00	-	
3'-Acridine CPG	20-2973-41	4 x 1 µmol	262.00	288.00	241.00	-	
3'-Acridine CPG	20-2973-42	4 x 0.2 µmol	157.00	173.00	144.00	-	
3'-Alkyne-Modifier Serinol CPG	20-2992-01	0.1 g	138.00	152.00	127.00	-	
3'-Alkyne-Modifier Serinol CPG	20-2992-10	1 g	1,048.00	1,153.00	964.00	-	
3'-Alkyne-Modifier Serinol CPG	20-2992-41	4 x 1 µmol	229.00	252.00	211.00	-	
3'-Alkyne-Modifier Serinol CPG	20-2992-42	4 x 0.2 µmol	131.00	144.00	121.00	-	
3'-Alkyne-Modifier Serinol CPG	20-2992-42	4 x 0.2 µmol	131.00	144.00	121.00	-	
3'-Amino-Modifier C6 dC CPG	20-2019-01	0.1 g	157.00	173.00	144.00	-	
3'-Amino-Modifier C6 dC CPG	20-2019-10	1 g	1,304.00	1,434.00	1,200.00	-	
3'-Amino-Modifier C6 dC CPG	20-2019-13	1 x 10 µmol	393.00	432.00	362.00	-	
3'-Amino-Modifier C6 dC CPG	20-2019-14	1 x 15 µmol	590.00	649.00	543.00	-	
3'-Amino-Modifier C6 dC CPG	20-2019-41	4 x 1 µmol	262.00	288.00	241.00	-	
3'-Amino-Modifier C6 dC CPG	20-2019-42	4 x 0.2 µmol	157.00	173.00	144.00	-	
3'-Amino-Modifier C6 dT CPG	20-2039-01	0.1 g	126.00	139.00	116.00	-	
3'-Amino-Modifier C6 dT CPG	20-2039-10	1 g	1,048.00	1,153.00	964.00	-	
3'-Amino-Modifier C6 dT CPG	20-2039-13	1 x 10 µmol	315.00	347.00	290.00	-	
3'-Amino-Modifier C6 dT CPG	20-2039-14	1 x 15 µmol	472.00	519.00	434.00	-	
3'-Amino-Modifier C6 dT CPG	20-2039-41	4 x 1 µmol	210.00	231.00	193.00	-	
3'-Amino-Modifier C6 dT CPG	20-2039-42	4 x 0.2 µmol	126.00	139.00	116.00	-	
3'-Amino-Modifier C6 dT CPG, Expedite	20-2039-41E	4 x 1 µmol	210.00	231.00	193.00	-	

Description	Reference	Scale/Quantity	Price List 2021				Ice Shipping
			EUR	CHF	GBP	USD	
3'-Amino-Modifier C7 CPG	20-2957-01	0.1 g	136.00	150.00	125.00	-	
3'-Amino-Modifier C7 CPG	20-2957-10	1 g	966.00	1,063.00	889.00	-	
3'-Amino-Modifier C7 CPG	20-2957-13	1 x 10 µmol	358.00	394.00	329.00	-	
3'-Amino-Modifier C7 CPG	20-2957-14	1 x 15 µmol	537.00	591.00	494.00	-	
3'-Amino-Modifier C7 CPG	20-2957-41	4 x 1 µmol	200.00	220.00	184.00	-	
3'-Amino-Modifier C7 CPG	20-2957-42	4 x 0.2 µmol	122.00	134.00	112.00	-	
3'-Amino-Modifier C7 CPG 1000	20-2958-01	0.1 g	124.00	136.00	114.00	-	
3'-Amino-Modifier C7 CPG 1000	20-2958-10	1 g	885.00	974.00	814.00	-	
3'-Amino-Modifier C7 CPG 1000	20-2958-13	1 x 10 µmol	328.00	361.00	302.00	-	
3'-Amino-Modifier C7 CPG 1000	20-2958-14	1 x 15 µmol	491.00	540.00	452.00	-	
3'-Amino-Modifier C7 CPG 1000	20-2958-41	4 x 1 µmol	155.00	171.00	143.00	-	
3'-Amino-Modifier C7 CPG 1000	20-2958-42	4 x 0.2 µmol	85.70	94.30	78.80	-	
3'-Amino-Modifier Serinol CPG	20-2997-01	0.1 g	124.00	136.00	114.00	-	
3'-Amino-Modifier Serinol CPG	20-2997-10	1 g	885.00	974.00	814.00	-	
3'-Amino-Modifier Serinol CPG	20-2997-13	1 x 10 µmol	328.00	361.00	302.00	-	
3'-Amino-Modifier Serinol CPG	20-2997-14	1 x 15 µmol	491.00	540.00	452.00	-	
3'-Amino-Modifier Serinol CPG	20-2997-41	4 x 1 µmol	183.00	201.00	168.00	-	
3'-Amino-Modifier Serinol CPG	20-2997-42	4 x 0.2 µmol	111.00	122.00	102.00	-	
3'-BBQ 650 CPG	20-5934-42	4 x 0.2 µmol	138.00	152.00	127.00	-	
3'-BHQ-1 CPG	20-5931-01	0.1 g	360.00	396.00	331.00	-	
3'-BHQ-1 CPG	20-5931-10	1 g	3,014.00	3,315.00	2,773.00	-	
3'-BHQ-1 CPG	20-5931-13	1 x 10 µmol	944.00	1,038.00	868.00	-	
3'-BHQ-1 CPG	20-5931-14	1 x 15 µmol	1,409.00	1,550.00	1,296.00	-	
3'-BHQ-1 CPG	20-5931-41	4 x 1 µmol	366.00	403.00	337.00	-	
3'-BHQ-1 CPG	20-5931-42	4 x 0.2 µmol	138.00	152.00	127.00	-	
3'-BHQ-1 CPG, Mermade	20-5931-42M	4 x 0.2 µmol	138.00	152.00	127.00	-	
3'-BHQ-1 CPG-1.0gram	20-5931-10D		3,014.00	3,315.00	2,773.00	-	
3'-BHQ-2 CPG	20-5932-01	0.1 g	360.00	396.00	331.00	-	
3'-BHQ-2 CPG	20-5932-10	1 g	3,014.00	3,315.00	2,773.00	-	
3'-BHQ-2 CPG	20-5932-10D	1 g	3,014.00	3,315.00	2,773.00	-	
3'-BHQ-2 CPG	20-5932-13	1 x 10 µmol	944.00	1,038.00	868.00	-	
3'-BHQ-2 CPG	20-5932-14	1 x 15 µmol	1,409.00	1,550.00	1,296.00	-	
3'-BHQ-2 CPG	20-5932-41	4 x 1 µmol	344.00	378.00	316.00	-	
3'-BHQ-2 CPG	20-5932-42	4 x 0.2 µmol	138.00	152.00	127.00	-	
3'-BHQ-2 CPG, Mermade	20-5932-42M	4 x 0.2 µmol	138.00	152.00	127.00	-	
3'-BHQ-3 CPG	20-5933-01	0.1 g	360.00	396.00	331.00	-	
3'-BHQ-3 CPG	20-5933-10	1 g	3,014.00	3,315.00	2,773.00	-	
3'-BHQ-3 CPG	20-5933-13	1 x 10 µmol	1,081.00	1,189.00	995.00	-	
3'-BHQ-3 CPG	20-5933-14	1 x 15 µmol	1,625.00	1,788.00	1,495.00	-	
3'-BHQ-3 CPG	20-5933-41	4 x 1 µmol	491.00	540.00	452.00	-	
3'-BHQ-3 CPG	20-5933-42	4 x 0.2 µmol	151.00	166.00	139.00	-	
3'-BHQ-3 CPG-4x1.0um	20-5933-41E		491.00	540.00	452.00	-	
3'-BiotinTEG CPG	20-2955-01	0.1 g	157.00	173.00	144.00	-	
3'-BiotinTEG CPG	20-2955-10	1 g	1,304.00	1,434.00	1,200.00	-	
3'-BiotinTEG CPG	20-2955-13	1 x 10 µmol	393.00	432.00	362.00	-	
3'-BiotinTEG CPG	20-2955-14	1 x 15 µmol	590.00	649.00	543.00	-	
3'-BiotinTEG CPG	20-2955-41	4 x 1 µmol	222.00	244.00	204.00	-	
3'-BiotinTEG CPG	20-2955-42	4 x 0.2 µmol	121.00	133.00	111.00	-	
3'-BiotinTEG PS	26-2955-52	10 x 200 nmol	393.00	432.00	362.00	-	
3'-BiotinTEG PS	26-2955-55	10 x 40 nmol	393.00	432.00	362.00	-	
3'-Cholesteryl-TEG CPG	20-2975-01	0.1 g	111.00	122.00	102.00	-	
3'-Cholesteryl-TEG CPG	20-2975-10	1 g	917.00	1,009.00	844.00	-	
3'-Cholesteryl-TEG CPG	20-2975-10	1 g	917.00	1,009.00	844.00	-	
3'-Cholesteryl-TEG CPG	20-2975-13	1 x 10 µmol	275.00	303.00	253.00	-	
3'-Cholesteryl-TEG CPG	20-2975-14	1 x 15 µmol	413.00	454.00	380.00	-	
3'-Cholesteryl-TEG CPG	20-2975-41	4 x 1 µmol	183.00	201.00	168.00	-	
3'-Cholesteryl-TEG CPG	20-2975-42	4 x 0.2 µmol	110.00	121.00	101.00	-	
3'-CPR II CPG	20-2903-10	1 g	629.00	692.00	579.00	-	
3'-CPR II CPG, Mermade	20-2903-41M	4 x 0.1 µmol	131.00	144.00	121.00	-	
3'-CPR II CPG, Mermade	20-2903-42M	4 x 0.2 µmol	78.60	86.50	72.30	-	
3-Cyanovinylcarbazole Phosphoramidite (CNVK), Expedite	10-4960-90E	100 µmol	511.00	562.00	470.00	-	
3'-DabcyI CPG	20-5912-01	0.1 g	145.00	160.00	133.00	-	
3'-DabcyI CPG	20-5912-10	1 g	1,179.00	1,297.00	1,085.00	-	
3'-DabcyI CPG	20-5912-13	1 x 10 µmol	423.00	465.00	389.00	-	
3'-DabcyI CPG	20-5912-14	1 x 15 µmol	605.00	666.00	557.00	-	
3'-DabcyI CPG	20-5912-41	4 x 1 µmol	222.00	244.00	204.00	-	
3'-DabcyI CPG	20-5912-42	4 x 0.2 µmol	121.00	133.00	111.00	-	
3'-DabcyI PS	26-5912-52	10 x 200 nmol	393.00	432.00	362.00	-	
3'-DabcyI PS	26-5912-55	10 x 40 nmol	393.00	432.00	362.00	-	
3'-DabcyI-CPG	20-5912-SP	4 x 1.3 µmol	333.00	366.00	306.00	-	
3'-DabcyI-CPG-0.2µmol col.(1pk/4)Mermade	20-5912-42M	4 x 0.2 µmol	121.00	133.00	111.00	-	
3'-DabcyI-CPG-1000 col.	20-5912-41E	4 x 1 µmol	222.00	244.00	204.00	-	
3'-DabcyI-CPG1000 col1µM(1pk/4)Mermade	20-5912-41M	4 x 1 µmol	222.00	244.00	204.00	-	
3'-Dabsyl CPG	20-5911-01	0.1 g	157.00	173.00	144.00	-	
3'-Dabsyl CPG	20-5911-10	1 g	1,278.00	1,406.00	1,176.00	-	
3'-Dabsyl CPG	20-5911-13	1 x 10 µmol	459.00	505.00	422.00	-	
3'-Dabsyl CPG	20-5911-14	1 x 15 µmol	655.00	721.00	603.00	-	
3'-Dabsyl CPG	20-5911-41	4 x 1 µmol	222.00	244.00	204.00	-	
3'-Dabsyl CPG	20-5911-42	4 x 0.2 µmol	121.00	133.00	111.00	-	
3'-Dabsyl-CPG Col 0.2 µmol pk/4 Mermade	20-5911-42M	0.2 µmol	121.00	133.00	111.00	-	
3'-Dabsyl-CPG col.	20-5911-41E	4 x 1 µmol	222.00	244.00	204.00	-	
3'-Dabsyl-CPG col. Expedite	20-5911-42E	0.2 µmol	121.00	133.00	111.00	-	
3'-Dabsyl-CPG-Col-1µM (4/pack)Mermade	20-5911-41M	4 x 1 µmol	222.00	244.00	204.00	-	
3'-dA-CE Phosphoramidite	10-1004-02	0.25 g	1,638.00	1,802.00	1,507.00	-	
3'-dA-CE Phosphoramidite	10-1004-90	100 µmol	590.00	649.00	543.00	-	
3'-dA-CE Phosphoramidite	10-1004-95	50 µmol	295.00	325.00	271.00	-	
3'-dA-CPG	20-2004-01	0.1 g	688.00	757.00	633.00	-	
3'-dA-CPG	20-2104-41	4 x 1 µmol	885.00	974.00	814.00	-	
3'-dA-CPG	20-2104-42	4 x 0.2 µmol	315.00	347.00	290.00	-	
3'-dC-CE Phosphoramidite	10-1064-02	0.25 g	2,490.00	2,739.00	2,291.00	-	
3'-dC-CE Phosphoramidite	10-1064-90	100 µmol	858.00	944.00	789.00	-	

Description	Reference	Scale/Quantity	Price List 2021				Ice Shipping
			EUR	CHF	GBP	USD	
3'-dC-CE Phosphoramidite	10-1064-95	50 µmol	426.00	469.00	392.00	-	
3'-dC-CPG	20-2064-01	0.1 g	393.00	432.00	362.00	-	
3'-dC-CPG	20-2164-41	4 x 1 µmol	786.00	865.00	723.00	-	
3'-dC-CPG	20-2164-42	4 x 0.2 µmol	262.00	288.00	241.00	-	
3-deaza-dA-CE Phosphoramidite	10-1088-02	0.25 g	1,396.00	1,536.00	1,284.00	-	
3-deaza-dA-CE Phosphoramidite	10-1088-90	100 µmol	508.00	559.00	467.00	-	
3-deaza-dA-CE Phosphoramidite	10-1088-95	50 µmol	254.00	279.00	234.00	-	
3'-dG-CE Phosphoramidite	10-1074-02	0.25 g	1,638.00	1,802.00	1,507.00	-	
3'-dG-CE Phosphoramidite	10-1074-90	100 µmol	590.00	649.00	543.00	-	
3'-dG-CE Phosphoramidite	10-1074-95	50 µmol	295.00	325.00	271.00	-	
3'-dG-CPG	20-2074-01	0.1 g	393.00	432.00	362.00	-	
3'-dG-CPG	20-2174-41	4 x 1 µmol	786.00	865.00	723.00	-	
3'-dG-CPG	20-2174-42	4 x 0.2 µmol	262.00	288.00	241.00	-	
3'-dT-CE Phosphoramidite	10-1084-02	0.25 g	2,064.00	2,270.00	1,899.00	-	
3'-dT-CE Phosphoramidite	10-1084-90	100 µmol	642.00	706.00	591.00	-	
3'-dT-CE Phosphoramidite	10-1084-95	50 µmol	334.00	367.00	307.00	-	
3'-dT-CPG	20-2084-01	0.1 g	393.00	432.00	362.00	-	
3'-dT-CPG	20-2184-41	4 x 1 µmol	786.00	865.00	723.00	-	
3'-dT-CPG	20-2184-42	4 x 0.2 µmol	262.00	288.00	241.00	-	
3'-DTPA CPG	20-2937-01	0.1 g	179.00	197.00	165.00	-	
3'-DTPA CPG	20-2937-10	1 g	1,281.00	1,409.00	1,179.00	-	
3'-DTPA CPG	20-2937-14	1 x 15 µmol	744.00	818.00	684.00	-	
3'-DTPA CPG	20-2937-41	4 x 1 µmol	265.00	292.00	244.00	-	
3'-DTPA CPG	20-2937-42	4 x 0.2 µmol	157.00	173.00	144.00	-	
3'-Fluorescein CPG	20-2963-01	0.1 g	157.00	173.00	144.00	-	
3'-Fluorescein CPG	20-2963-10	1 g	1,304.00	1,434.00	1,200.00	-	
3'-Fluorescein CPG	20-2963-13	1 x 10 µmol	393.00	432.00	362.00	-	
3'-Fluorescein CPG	20-2963-14	1 x 15 µmol	590.00	649.00	543.00	-	
3'-Fluorescein CPG	20-2963-41	4 x 1 µmol	262.00	288.00	241.00	-	
3'-Fluorescein CPG	20-2963-42	4 x 0.2 µmol	157.00	173.00	144.00	-	
3'-Fluorescein CPG, Expedite	20-2963-41E	4 x 1 µmol	262.00	288.00	241.00	-	
3'-Fluorescein-dT CPG	20-2056-01	0.1 g	157.00	173.00	144.00	-	
3'-Fluorescein-dT CPG	20-2056-10	1 g	1,304.00	1,434.00	1,200.00	-	
3'-Fluorescein-dT CPG	20-2056-13	1 x 10 µmol	393.00	432.00	362.00	-	
3'-Fluorescein-dT CPG	20-2056-14	1 x 15 µmol	590.00	649.00	543.00	-	
3'-Fluorescein-dT CPG	20-2056-41	4 x 1 µmol	262.00	288.00	241.00	-	
3'-Fluorescein-dT CPG	20-2056-42	4 x 0.2 µmol	157.00	173.00	144.00	-	
3'-Glyceryl CPG	20-2902-01	0.1 g	111.00	122.00	102.00	-	
3'-Glyceryl CPG	20-2902-10	1 g	786.00	865.00	723.00	-	
3'-Glyceryl CPG	20-2902-13	1 x 10 µmol	295.00	325.00	271.00	-	
3'-Glyceryl CPG	20-2902-14	1 x 15 µmol	459.00	505.00	422.00	-	
3'-Glyceryl CPG	20-2902-41	4 x 1 µmol	164.00	180.00	151.00	-	
3'-Glyceryl CPG	20-2902-42	4 x 0.2 µmol	98.30	108.10	90.40	-	
3'-Glyceryl CPG, column, Pharmacia	20-2902-41P	1 µmol (4 x 1 pce)	164.00	180.00	151.00	-	
3'-Phosphate CPG	20-2900-01	0.1 g	91.70	100.90	84.40	-	
3'-Phosphate CPG	20-2900-10	1 g	629.00	692.00	579.00	-	
3'-Phosphate CPG	20-2900-13	1 x 10 µmol	236.00	260.00	217.00	-	
3'-Phosphate CPG	20-2900-14	1 x 15 µmol	367.00	404.00	338.00	-	
3'-Phosphate CPG	20-2900-41	4 x 1 µmol	111.00	122.00	102.00	-	
3'-Phosphate CPG	20-2900-42	4 x 0.2 µmol	60.50	66.60	55.70	-	
3'-Phosphate CPG (High Load)	25-2900-01	0.1 g	111.00	122.00	102.00	-	
3'-Phosphate CPG (High Load)	25-2900-10	1 g	786.00	865.00	723.00	-	
3'-Phosphate CPG, column	20-2900-41P	4 x 1 µmol	111.00	122.00	102.00	-	
3'-Phosphate PS	26-2900-52	10 x 200 nmol	197.00	217.00	181.00	-	
3'-Phosphate PS	26-2900-55	10 x 40 nmol	197.00	217.00	181.00	-	
3'-PhosphateCPG-0.2µm.col(1pk/4)Mermade	20-2900-42M	0.2 µmol	60.50	66.60	55.70	-	
3'-PhosphateCPG-1 µm.col(1pk/4)Mermade	20-2900-41M	1 µmol	111.00	122.00	102.00	-	
3'-Propargyl-5-Me-dC CPG	20-2982-01	0.1 g	249.00	274.00	229.00	-	
3'-Propargyl-5-Me-dC CPG, column	20-2982-41	4 x 1 µmol	498.00	548.00	458.00	-	
3'-Propargyl-5-Me-dC CPG, column	20-2982-42	4 x 0.2 µmol	197.00	217.00	181.00	-	
3'-Protected Biotin Serinol CPG	20-2993-01	0.1 g	157.00	173.00	144.00	-	
3'-Protected Biotin Serinol CPG	20-2993-10	1 g	1,304.00	1,434.00	1,200.00	-	
3'-Protected Biotin Serinol CPG	20-2993-13	1 x 10 µmol	393.00	432.00	362.00	-	
3'-Protected Biotin Serinol CPG	20-2993-14	1 x 15 µmol	590.00	649.00	543.00	-	
3'-Protected Biotin Serinol CPG	20-2993-41	4 x 1 µmol	262.00	288.00	241.00	-	
3'-Protected Biotin Serinol CPG	20-2993-42	4 x 0.2 µmol	157.00	173.00	144.00	-	
3'-Protected BiotinLC Serinol CPG	20-2995-01	0.1 g	157.00	173.00	144.00	-	
3'-Protected BiotinLC Serinol CPG	20-2995-10	1 g	1,304.00	1,434.00	1,200.00	-	
3'-Protected BiotinLC Serinol CPG	20-2995-13	1 x 10 µmol	393.00	432.00	362.00	-	
3'-Protected BiotinLC Serinol CPG	20-2995-14	1 x 15 µmol	590.00	649.00	543.00	-	
3'-Protected BiotinLC Serinol CPG	20-2995-41	4 x 1 µmol	262.00	288.00	241.00	-	
3'-Protected BiotinLC Serinol CPG	20-2995-42	4 x 0.2 µmol	157.00	173.00	144.00	-	
3'-Protected BiotinLC Serinol CPG, Expedite	20-2995-42E	4 x 0.2 µmol	157.00	173.00	144.00	-	
3'-PT-Amino-Modifier C3 CPG	20-2954-01	0.1 g	124.00	136.00	114.00	-	
3'-PT-Amino-Modifier C3 CPG	20-2954-10	1 g	885.00	974.00	814.00	-	
3'-PT-Amino-Modifier C3 CPG	20-2954-13	1 x 10 µmol	328.00	361.00	302.00	-	
3'-PT-Amino-Modifier C3 CPG	20-2954-14	1 x 15 µmol	491.00	540.00	452.00	-	
3'-PT-Amino-Modifier C3 CPG	20-2954-41	4 x 1 µmol	183.00	201.00	168.00	-	
3'-PT-Amino-Modifier C3 CPG	20-2954-42	4 x 0.2 µmol	111.00	122.00	102.00	-	
3'-PT-Amino-Modifier C6 CPG	20-2956-01	0.1 g	124.00	136.00	114.00	-	
3'-PT-Amino-Modifier C6 CPG	20-2956-10	1 g	885.00	974.00	814.00	-	
3'-PT-Amino-Modifier C6 CPG	20-2956-13	1 x 10 µmol	328.00	361.00	302.00	-	
3'-PT-Amino-Modifier C6 CPG	20-2956-14	1 x 15 µmol	491.00	540.00	452.00	-	
3'-PT-Amino-Modifier C6 CPG	20-2956-41	4 x 1 µmol	183.00	201.00	168.00	-	
3'-PT-Amino-Modifier C6 CPG	20-2956-42	4 x 0.2 µmol	111.00	122.00	102.00	-	
3'-PT-Amino-Modifier C6 CPG,Expedite	20-2956-42E	4 x 0.2 µmol	111.00	122.00	102.00	-	
3'-PT-Amino-Modifier C6 PS	26-2956-10	1 g	1,343.00	1,477.00	1,236.00	-	
3'-PT-Amino-Modifier C6 PS	26-2956-52	10 x 200 nmol	288.00	317.00	265.00	-	
3'-PT-Amino-Modifier C6 PS	26-2956-55	10 x 40 nmol	288.00	317.00	265.00	-	
3'-Spacer C3 CPG	20-2913-01	0.1 g	91.70	100.90	84.40	-	
3'-Spacer C3 CPG	20-2913-10	1 g	629.00	692.00	579.00	-	

Description	Reference	Scale/Quantity	Price List 2021				Ice Shipping
			EUR	CHF	GBP	USD	
3'-Spacer C3 CPG	20-2913-13	1 x 10 µmol	236.00	260.00	217.00	-	
3'-Spacer C3 CPG	20-2913-14	1 x 15 µmol	367.00	404.00	338.00	-	
3'-Spacer C3 CPG	20-2913-41	4 x 1 µmol	131.00	144.00	121.00	-	
3'-Spacer C3 CPG	20-2913-42	4 x 0.2 µmol	78.60	86.50	72.30	-	
3'-TAMRA CPG	20-5910-01	0.1 g	145.00	160.00	133.00	-	
3'-TAMRA CPG	20-5910-10	1 g	1,204.00	1,324.00	1,108.00	-	
3'-TAMRA CPG	20-5910-41	4 x 1 µmol	202.00	222.00	186.00	-	
3'-TAMRA CPG	20-5910-42	4 x 0.2 µmol	96.80	106.50	89.10	-	
3'-TAMRA PS	26-5910-52	10 x 200 nmol	393.00	432.00	362.00	-	
3'-TAMRA PS	26-5910-55	10 x 40 nmol	393.00	432.00	362.00	-	
3'-Thiol-Modifier 6 S-S CPG	20-2938-41E	4 x 1 µmol	164.00	180.00	151.00	-	
3'-Thiol-Modifier 6 S-S CPG	20-2938-42	4 x 0.2 µmol	98.30	108.10	90.40	-	
3'-Thiol-Modifier 6 S-S CPG	20-2938-42E	4 x 0.2 µmol	98.30	108.10	90.40	-	
3'-Thiol-Modifier C3 S-S CPG	20-2933-01	0.1 g	103.00	113.00	95.00	-	
3'-Thiol-Modifier C3 S-S CPG	20-2933-10	1 g	726.00	799.00	668.00	-	
3'-Thiol-Modifier C3 S-S CPG	20-2933-13	1 x 10 µmol	272.00	299.00	250.00	-	
3'-Thiol-Modifier C3 S-S CPG	20-2933-14	1 x 15 µmol	423.00	465.00	389.00	-	
3'-Thiol-Modifier C3 S-S CPG	20-2933-41	4 x 1 µmol	139.00	153.00	128.00	-	
3'-Thiol-Modifier C3 S-S CPG	20-2933-42	4 x 0.2 µmol	75.60	83.20	69.60	-	
3'-Thiol-Modifier C3 S-S CPG, Beckman	20-2933-41B		139.00	153.00	128.00	-	
3'-Thiol-Modifier C3 S-S CPG, Pharmacia	20-2933-41P		139.00	153.00	128.00	-	
3'-Thiol-Modifier C3 S-SCPG 1µmol 4 pc M	20-2933-41M		139.00	153.00	128.00	-	
3'-Thiol-Modifier C6 S-S CPG	20-2936-01	0.1 g	122.00	134.00	112.00	-	
3'-Thiol-Modifier C6 S-S CPG	20-2936-10	1 g	859.00	945.00	790.00	-	
3'-Thiol-Modifier C6 S-S CPG	20-2936-13	1 x 10 µmol	322.00	354.00	296.00	-	
3'-Thiol-Modifier C6 S-S CPG	20-2936-14	1 x 15 µmol	501.00	551.00	461.00	-	
3'-Thiol-Modifier C6 S-S CPG	20-2936-42	4 x 0.2 µmol	107.00	118.00	98.00	-	
3'-Thiol-Modifier C6 S-S CPG	20-2938-41	4x1.0µm	164.00	180.00	151.00	-	
3'-Thiol-Modifier C3 S-SCPG-0.2µm(1pk/4)M	20-2933-42M	4 x 0.2 µmol	75.60	83.20	69.60	-	
3'-Uaq Cap CPG	20-2980-01	0.1 g	236.00	260.00	217.00	-	
4-Thio-dT-CE Phosphoramidite	10-1034-02	0.25 g	885.00	974.00	814.00	-	
4-Thio-dT-CE Phosphoramidite	10-1034-90	100 µmol	387.00	426.00	356.00	-	
4-Thio-dT-CE Phosphoramidite	10-1034-95	50 µmol	216.00	238.00	199.00	-	
4-Thio-dU-CE Phosphoramidite	10-1052-02	0.25 g	1,147.00	1,262.00	1,055.00	-	
4-Thio-dU-CE Phosphoramidite	10-1052-90	100 µmol	393.00	432.00	362.00	-	
4-Thio-dU-CE Phosphoramidite	10-1052-95	50 µmol	216.00	238.00	199.00	-	
5' - DMS(O)MT-Amino-Modifier C6, Expedite	10-1907-02E	0.25 g	295.00	325.00	271.00	-	
5' Stearyl Phosphoramidite	10-1979-90	100 µmol	59.00	64.90	54.30	-	✓
5,6-Dihydro-dT-CE Phosphoramidite	10-1530-02	0.25 g	780.00	858.00	718.00	-	
5,6-Dihydro-dT-CE Phosphoramidite	10-1530-90	100 µmol	256.00	282.00	236.00	-	
5,6-Dihydro-dU-CE Phosphoramidite	10-1550-02	0.25 g	780.00	858.00	718.00	-	
5,6-Dihydro-dU-CE Phosphoramidite	10-1550-90	100 µmol	256.00	282.00	236.00	-	
5'-Aldehyde-Modifier C2 Phosphoramidite	10-1933-02	0.25 g	426.00	469.00	392.00	-	
5'-Aldehyde-Modifier C2 Phosphoramidite	10-1933-90	100 µmol	111.00	122.00	102.00	-	
5'-Aldehyde-Modifier C2 Phosphoramidite, Expedite	10-1933-90E	100 µmol	111.00	122.00	102.00	-	
5'-Amino-dT-CE Phosphoramidite	10-1932-02	0.25 g	524.00	576.00	482.00	-	
5'-Amino-dT-CE Phosphoramidite	10-1932-90	100 µmol	197.00	217.00	181.00	-	
5'-Amino-Modifier 5	10-1905-02	0.25 g	262.00	288.00	241.00	-	
5'-Amino-Modifier 5	10-1905-90	100 µmol	78.60	86.50	72.30	-	
5'-Amino-Modifier C12	10-1912-02	0.25 g	393.00	432.00	362.00	-	
5'-Amino-Modifier C12	10-1912-90	100 µmol	118.00	130.00	109.00	-	
5'-Amino-Modifier C3-TFA	10-1923-02	0.25 g	229.00	252.00	211.00	-	✓
5'-Amino-Modifier C3-TFA	10-1923-90	100 µmol	65.50	72.10	60.30	-	
5'-Amino-Modifier C6	10-1906-02	0.25 g	161.00	177.00	148.00	-	
5'-Amino-Modifier C6	10-1906-90	100 µmol	48.40	53.20	44.50	-	
5'-Amino-Modifier C6, Beckman	10-1906-90B	100 µmol	48.40	53.20	44.50	-	
5'-Amino-Modifier C6-PDA	10-1947-02	0.25 g	131.00	144.00	121.00	-	
5'-Amino-Modifier C6-PDA	10-1947-02	0.25 g	131.00	144.00	121.00	-	
5'-Amino-Modifier C6-TFA	10-1916-02	0.25 g	80.60	88.70	74.20	-	✓
5'-Amino-Modifier C6-TFA	10-1916-90	100 µmol	24.20	26.60	22.30	-	✓
5'-Amino-Modifier C6-TFA, Mermade	10-1916-02M	0.25 g	80.60	88.70	74.20	-	
5'-Amino-Modifier TEG CE-Phosphoramidite	10-1917-02	0.25 g	655.00	721.00	603.00	-	
5'-Amino-Modifier TEG CE-Phosphoramidite	10-1917-90	100 µmol	151.00	166.00	139.00	-	
5'-Amino-Modifier TEG CE-Phosphoramidite, Expedite	10-1917-02E	0.25 g	655.00	721.00	603.00	-	
5'-Amino-Modifier TEG CE-Phosphoramidite, Expedite	10-1917-90E	100 µmol	151.00	166.00	139.00	-	
5'-AminoOxy - Modifier 11 - CE - Phosphoramidite	10-1919-95	50 µmol	183.00	201.00	168.00	-	
5'-AminoOxy-Modifier 11-CE-Phosphoramidite	10-1919-02	0.25 g	1,173.00	1,290.00	1,079.00	-	
5'-AminoOxy-Modifier 11-CE-Phosphoramidite	10-1919-90	100 µmol	347.00	382.00	319.00	-	
5'-AquaPhluor® 593 CE Phosphoramidite	10-5923-90	100 µmol	1,042.00	1,146.00	959.00	-	
5'-BBQ-650-CE Phosphoramidite	10-5934-90	100 µmol	511.00	562.00	470.00	-	✓
5'-BHQ-1 Phosphoramidite	10-5931-02	0.25 g	1,376.00	1,514.00	1,266.00	-	
5'-BHQ-1 Phosphoramidite	10-5931-90	100 µmol	367.00	404.00	338.00	-	
5'-BHQ-1 Phosphoramidite	10-5931-95	50 µmol	193.00	212.00	178.00	-	
5'-BHQ-2 Phosphoramidite	10-5932-02	0.25 g	1,376.00	1,514.00	1,266.00	-	
5'-BHQ-2 Phosphoramidite	10-5932-90	100 µmol	393.00	432.00	362.00	-	
5'-BHQ-2 Phosphoramidite	10-5932-95	50 µmol	193.00	212.00	178.00	-	
5'-Biotin Phosphoramidite	10-5950-02	0.25 g	252.00	277.00	232.00	-	
5'-Biotin Phosphoramidite	10-5950-90	100 µmol	90.70	99.80	83.40	-	
5'-Biotin Phosphoramidite	10-5950-95	50 µmol	60.50	66.60	55.70	-	
5'-Biotin Phosphoramidite, Expedite	10-5950-95E	50 µmol	60.50	66.60	55.70	-	
5-Br-dC-CE Phosphoramidite	10-1080-02	0.25 g	210.00	231.00	193.00	-	
5-Br-dC-CE Phosphoramidite	10-1080-90	100 µmol	78.60	86.50	72.30	-	
5-Br-dU-CE Phosphoramidite	10-1090-02	0.25 g	161.00	177.00	148.00	-	
5-Br-dU-CE Phosphoramidite	10-1090-90	100 µmol	60.50	66.60	55.70	-	
5-Br-dU-CE Phosphoramidite, Expedite	10-1090-02E	0.25 g	161.00	177.00	148.00	-	
5-Br-dU-CPG	20-2090-01	0.1 g	65.50	72.10	60.30	-	
5-Br-dU-CPG	20-2090-41	4 x 1 µmol	262.00	288.00	241.00	-	
5-Br-dU-CPG	20-2090-42	4 x 0.2 µmol	157.00	173.00	144.00	-	
5'-Bromohexyl Phosphoramidit, Mermade	10-1946-02M	0.25 g	262.00	288.00	241.00	-	
5'-Bromohexyl Phosphoramidite	10-1946-02	0.25 g	262.00	288.00	241.00	-	
5'-Bromohexyl Phosphoramidite	10-1946-90	100 µmol	78.60	86.50	72.30	-	

Description	Reference	Scale/Quantity	Price List 2021				Ice Shipping
			EUR	CHF	GBP	USD	
5-Carboxy-dC-CE Phosphoramidite	10-1066-02	0.25 g	1,573.00	1,730.00	1,447.00	-	
5-Carboxy-dC-CE Phosphoramidite	10-1066-90	100 µmol	590.00	649.00	543.00	-	✓
5-Carboxy-dC-CE Phosphoramidite	10-1066-95	50 µmol	301.00	331.00	277.00	-	
5'-Carboxy-Modifier C10	10-1935-02	0.25 g	347.00	382.00	319.00	-	
5'-Carboxy-Modifier C10	10-1935-90	100 µmol	85.20	93.70	78.40	-	
5'-CDPI3 MGB™ Phosphoramidite	10-5924-90	100 µmol	1,822.00	2,004.00	1,676.00	-	
5'Cholesteryl-TEG Phospho.	10-1976-02	0.25 g	688.00	757.00	633.00	-	
5'Cholesteryl-TEG Phospho.	10-1976-02	0.25 g	688.00	757.00	633.00	-	
5'Cholesteryl-TEG Phospho.	10-1976-90	100 µmol	229.00	252.00	211.00	-	
5'Cholesteryl-TEG Phospho.	10-1976-95	50 µmol	124.00	136.00	114.00	-	
5'-Dabcyl Phosphoramidite	10-5912-02	0.25 g	852.00	937.00	784.00	-	
5'-Dabcyl Phosphoramidite	10-5912-90	100 µmol	295.00	325.00	271.00	-	
5'-Dabcyl Phosphoramidite	10-5912-95	50 µmol	164.00	180.00	151.00	-	
5'-Dabsyl-dT	10-1532-02	0.25 g	1,016.00	1,118.00	935.00	-	
5'-Dabsyl-dT	10-1532-90	100 µmol	328.00	361.00	302.00	-	
5'-DBCO-TEG Phosphoramidite	10-1941-95	50 µmol	164.00	180.00	151.00	-	
5'-DBCO-TEG Phosphoramidite, Mermade	10-1941-02M	0.25 g	1,016.00	1,118.00	935.00	-	
5'-Dichloro-dimethoxy-Fluorescein Phosphoramidite (Joe)	10-5904-90	100 µmol	487.00	536.00	448.00	-	
5'-Dichloro-dimethoxy-Fluorescein Phosphoramidite (Joe)	10-5904-95	50 µmol	251.00	276.00	231.00	-	
5'-Dichloro-dimethoxy-Fluorescein Phosphoramidite II ( JOE II)	10-5906-02	0.25 g	649.00	714.00	597.00	-	
5'-Dichloro-dimethoxy-Fluorescein Phosphoramidite II ( JOE II)	10-5906-90	100 µmol	259.00	285.00	238.00	-	
5'-Dimethoxytrityl-N-benzoyl-5-cyanoethoxy- methyl-2'-deoxyCytidine,3'-[[2-cyanoethyl]- (N,N-diisopropyl)]-phosphoramidite	10-1062-02	0.25 g	2,195.00	2,415.00	2,019.00	-	✓
5'-Dimethoxytrityl-N-benzoyl-5-cyanoethoxy- methyl-2'-deoxyCytidine,3'-[[2-cyanoethyl]- (N,N-diisopropyl)]-phosphoramidite	10-1062-90	100 µmol	852.00	937.00	784.00	-	
5'-Dimethoxytrityl-N-benzoyl-5-cyanoethoxy- methyl-2'-deoxyCytidine,3'-[[2-cyanoethyl]- (N,N-diisopropyl)]-phosphoramidite	10-1062-95	50 µmol	439.00	483.00	404.00	-	
5'-DMS(O)MT-Amino-Modifier C6	10-1907-02	0.25 g	295.00	325.00	271.00	-	
5'-DMS(O)MT-Amino-Modifier C6	10-1907-90	100 µmol	105.00	116.00	97.00	-	
5-Ethynyl-dU-CE Phosphoramidite	10-1554-02	0.25 g	1,016.00	1,118.00	935.00	-	
5-Ethynyl-dU-CE Phosphoramidite	10-1554-90	100 µmol	321.00	353.00	295.00	-	
5-F-dU-CE Phosphoramidite	10-1092-02	0.25 g	465.00	512.00	428.00	-	
5-F-dU-CE Phosphoramidite	10-1092-90	100 µmol	177.00	195.00	163.00	-	
5'-Fluorescein Phosphoramidite	10-5901-02	0.25 g	192.00	211.00	177.00	-	
5'-Fluorescein Phosphoramidite	10-5901-90	100 µmol	75.60	83.20	69.60	-	
5'-Fluorescein Phosphoramidite	10-5901-90P	100 µmol	75.60	83.20	69.60	-	
5'-Fluorescein Phosphoramidite	10-5901-95	50 µmol	50.40	55.40	46.40	-	
5'-Fluorescein Phosphoramidite, Expedite	10-5901-02E	0.25 g	192.00	211.00	177.00	-	
5'-Fluorescein Phosphoramidite, Mermade	10-5901-02M	0.25 g	192.00	211.00	177.00	-	
5-Formyl dC III CE Phosphoramidite	10-1564-90	100 µmol	1,442.00	1,586.00	1,327.00	-	
5-Formyl-dC-CE Phosphoramidite	10-1514-02	0.25 g	4,226.00	4,649.00	3,888.00	-	
5-Formyl-dC-CE Phosphoramidite	10-1514-95	50 µmol	799.00	879.00	735.00	-	
5-Formylindole-CE Phosphoramidite	10-1934-02	0.25 g	491.00	540.00	452.00	-	
5-Formylindole-CE Phosphoramidite	10-1934-90	100 µmol	151.00	166.00	139.00	-	
5-Formylindole-CE Phosphoramidite, Expedite	10-1934-90E	100 µmol	151.00	166.00	139.00	-	
5'-GalNAc C3 Phosphoramidite	10-1974-02	0.25 g	655.00	721.00	603.00	-	
5'-GalNAc C3 Phosphoramidite	10-1974-90	100 µmol	334.00	367.00	307.00	-	
5'-GalNAc C3 Phosphoramidite, Mermade	10-1974-90M	100 µmol	334.00	367.00	307.00	-	
5'-Hexachloro-Fluorescein Phosphoramidite	10-5901-BULK	1 g	795.00	875.00	731.00	-	
5'-Hexachloro-Fluorescein Phosphoramidite	10-5902-02	0.25 g	756.00	832.00	696.00	-	
5'-Hexachloro-Fluorescein Phosphoramidite	10-5902-90	100 µmol	333.00	366.00	306.00	-	
5'-Hexachloro-Fluorescein Phosphoramidite	10-5902-95	50 µmol	171.00	188.00	157.00	-	
5'-Hexynyl Phosphoramidite	10-1908-02	0.25 g	262.00	288.00	241.00	-	
5'-Hexynyl Phosphoramidite	10-1908-90	100 µmol	78.60	86.50	72.30	-	
5'-Hexynyl Phosphoramidite, Expedite	10-1908-02E	0.25 g	262.00	288.00	241.00	-	
5-Hydroxymethyl dC II CE Phosphoramidite	10-1510-90	100 µmol	878.00	966.00	808.00	-	
5-Hydroxymethyl-dC II-CE Phosphoramidite	10-1510-95	50 µmol	452.00	497.00	416.00	-	✓
5-Hydroxymethyl-dU-CE Phosphoramidite	10-1093-02	0.25 g	885.00	974.00	814.00	-	
5-Hydroxymethyl-dU-CE Phosphoramidite	10-1093-90	100 µmol	295.00	325.00	271.00	-	
5-I-dC-CE Phosphoramidite	10-1081-02	0.25 g	465.00	512.00	428.00	-	
5-I-dC-CE Phosphoramidite	10-1081-90	100 µmol	177.00	195.00	163.00	-	
5-I-dT-CE Phosphoramidite	10-1931-02	0.25 g	387.00	426.00	356.00	-	
5-I-dT-CE Phosphoramidite	10-1931-90	100 µmol	111.00	122.00	102.00	-	
5-I-dU-CE Phosphoramidite	10-1091-02	0.25 g	210.00	231.00	193.00	-	
5-I-dU-CE Phosphoramidite	10-1091-90	100 µmol	78.60	86.50	72.30	-	
5-I-dU-CE Phosphoramidite, Expedite	10-1091-02E	0.25 g	210.00	231.00	193.00	-	
5-I-dU-CE Phosphoramidite, Expedite	10-1091-90E	100 µmol	78.60	86.50	72.30	-	
5'-Maleimide-Modifier Phosphoramidite	10-1938-02	0.25 g	439.00	483.00	404.00	-	
5'-Maleimide-Modifier-Phosphoramidite	10-1938-90	100 µmol	91.70	100.90	84.40	-	
5-Me-2'-deoxyZebularine-CE Phosphoramidite	10-1061-02	0.25 g	1,310.00	1,441.00	1,205.00	-	
5-Me-2'-deoxyZebularine-CE Phosphoramidite	10-1061-90	100 µmol	524.00	576.00	482.00	-	
5-Me-2'-deoxyZebularine-CE Phosphoramidite	10-1061-95	50 µmol	275.00	303.00	253.00	-	
5-Me-dC-CE Phosphoramidite	10-1060-02	0.25 g	121.00	133.00	111.00	-	
5-Me-dC-CE Phosphoramidite	10-1060-90	100 µmol	50.40	55.40	46.40	-	
5-Me-dC-CPG	20-2060-01	0.1 g	72.00	79.20	66.20	-	
5-Me-dC-CPG	20-2160-42	4 x 0.2 µmol	172.00	189.00	158.00	-	
5-Methyl-Uridine a-thiotriphosphate (10mM) (5-Me-UTTPaS)	80-3093-01	0.1 mL	143.00	157.00	132.00	-	
5-Nitroindole-CE Phosphoramidite	10-1044-02	0.25 g	426.00	469.00	392.00	-	
5-Nitroindole-CE Phosphoramidite	10-1044-90	100 µmol	164.00	180.00	151.00	-	
5-O-Dimethoxytrityl-1-O-tert-butylidimethylsilyl-2-deoxyribose- 3'-[[2-cyanoethyl]- (N,N-diisopropyl)]-phosphoramidite	10-1927-02	250 mg	622.00	684.00	572.00	-	
5-O-Dimethoxytrityl-1-O-tert-butylidimethylsilyl-2-deoxyribose- 3'-[[2-cyanoethyl]- (N,N-diisopropyl)]-phosphoramidite	10-1927-90	100 µmol	197.00	217.00	181.00	-	
5-O-Dimethoxytrityl-1-O-tert-butylidimethylsilyl-2-deoxyribose- 3'-[[2-cyanoethyl]- (N,N-diisopropyl)]-phosphoramidite	10-1927-95	50 µmol	105.00	116.00	97.00	-	
5-OH-dC-CE Phosphoramidite	10-1063-02	0.25 g	1,147.00	1,262.00	1,055.00	-	
5-OH-dC-CE Phosphoramidite	10-1063-90	100 µmol	426.00	469.00	392.00	-	
5-OH-dU-CE Phosphoramidite	10-1053-02	0.25 g	1,147.00	1,262.00	1,055.00	-	
5-OH-dU-CE Phosphoramidite	10-1053-90	100 µmol	393.00	432.00	362.00	-	
5'-OMe-dT-CE Phosphoramidite	10-1031-02	0.25 g	465.00	512.00	428.00	-	
5'-OMe-dT-CE Phosphoramidite	10-1031-90	100 µmol	177.00	195.00	163.00	-	
5'-Pyrene Cap Phosphoramidite	10-1987-02	0.25 g	649.00	714.00	597.00	-	
5'-Pyrene Cap Phosphoramidite	10-1987-90	100 µmol	256.00	282.00	236.00	-	
5'-Tetrachloro-Fluorescein Phosphoramidite	10-5903-02	0.25 g	897.00	987.00	825.00	-	

Description	Reference	Scale/Quantity	Price List 2021				Ice Shipping
			EUR	CHF	GBP	USD	
5'-Tetrachloro-Fluorescein Phosphoramidite	10-5903-90	100 µmol	373.00	410.00	343.00	-	
5'-Tetrachloro-Fluorescein Phosphoramidite	10-5903-95	50 µmol	192.00	211.00	177.00	-	
5'-Thiol-Modifier C6	10-1926-02	0.25 g	262.00	288.00	241.00	-	
5'-Thiol-Modifier C6	10-1926-90	100 µmol	78.60	86.50	72.30	-	
5'-Thiol-Modifier C6	10-1926-90E	100 µmol	78.60	86.50	72.30	-	
5'-Trimethoxystilbene Cap Phosphoramidite	10-1986-02	0.25 g	655.00	721.00	603.00	-	
5'-Trimethoxystilbene Cap Phosphoramidite	10-1986-90	100 µmol	256.00	282.00	236.00	-	
6-FAM CPG 1000 -1µm (1pk/4)Mermade	20-2961-41M	4 x 1 µmol	222.00	244.00	204.00	-	
6-Fluorescein Phosphoramidite	10-1964-02	0.25 g	780.00	858.00	718.00	-	
6-Fluorescein Phosphoramidite	10-1964-90	100 µmol	387.00	426.00	356.00	-	
6-Fluorescein Phosphoramidite	10-1964-95	50 µmol	216.00	238.00	199.00	-	
6-Fluorescein Serinol Phosphoramidite	10-1994-02	0.25 g	780.00	858.00	718.00	-	
6-Fluorescein Serinol Phosphoramidite	10-1994-90	100 µmol	387.00	426.00	356.00	-	
6-Fluorescein Serinol Phosphoramidite	10-1994-95	50 µmol	216.00	238.00	199.00	-	
6-thio-dG-CE Phosphoramidite	10-1072-02	0.25 g	1,278.00	1,406.00	1,176.00	-	
6-thio-dG-CE Phosphoramidite	10-1072-90	100 µmol	465.00	512.00	428.00	-	
6-thio-dG-CE Phosphoramidite	10-1072-95	50 µmol	233.00	256.00	214.00	-	
6-Thio-G-CE Phosphoramidite	10-3072-02	0.25 g	1,573.00	1,730.00	1,447.00	-	
6-Thio-G-CE Phosphoramidite	10-3072-90	100 µmol	655.00	721.00	603.00	-	
6-Thio-G-CE Phosphoramidite	10-3072-95	50 µmol	328.00	361.00	302.00	-	
7-deaza-8-aza-dA-CE Phosphoramidite	10-1083-02	0.25 g	1,278.00	1,406.00	1,176.00	-	
7-deaza-8-aza-dA-CE Phosphoramidite	10-1083-90	100 µmol	465.00	512.00	428.00	-	
7-deaza-8-aza-dA-CE Phosphoramidite	10-1083-95	50 µmol	233.00	256.00	214.00	-	
7-Deaza-8-aza-dG-CE Phosphoramidite (PPG )	10-1073-02	0.25 g	2,097.00	2,307.00	1,929.00	-	
7-Deaza-8-aza-dG-CE Phosphoramidite (PPG )	10-1073-90	100 µmol	721.00	793.00	663.00	-	
7-Deaza-8-aza-dG-CE Phosphoramidite (PPG )	10-1073-95	50 µmol	373.00	410.00	343.00	-	
7-deaza-A-CE Phosphoramidite	10-3001-02	0.25 g	1,574.00	1,731.00	1,448.00	-	
7-deaza-A-CE Phosphoramidite	10-3001-90	100 µmol	644.00	708.00	592.00	-	
7-deaza-A-CE Phosphoramidite	10-3001-95	50 µmol	322.00	354.00	296.00	-	
7-Deaza-Adenosine a-thiotriphosphate (10mM) (7-DATTPaS)	80-3303-01	0.1 mL	143.00	157.00	132.00	-	
7-Deaza-dA-CE Phosphoramidite	10-1001-02	0.25 g	1,278.00	1,406.00	1,176.00	-	
7-Deaza-dA-CE Phosphoramidite	10-1001-90	100 µmol	465.00	512.00	428.00	-	
7-Deaza-dA-CE Phosphoramidite	10-1001-95	50 µmol	233.00	256.00	214.00	-	
7-deaza-dG-CE Phosphoramidite	10-1021-02	0.25 g	1,278.00	1,406.00	1,176.00	-	
7-deaza-dG-CE Phosphoramidite	10-1021-90	100 µmol	465.00	512.00	428.00	-	
7-deaza-dG-CE Phosphoramidite	10-1021-95	50 µmol	233.00	256.00	214.00	-	
7-deaza-dX-CE Phosphoramidite	10-1076-02	0.25 g	1,396.00	1,536.00	1,284.00	-	
7-deaza-dX-CE Phosphoramidite	10-1076-90	100 µmol	508.00	559.00	467.00	-	
7-deaza-dX-CE Phosphoramidite	10-1076-95	50 µmol	254.00	279.00	234.00	-	
7-deaza-guanosine thiotriphosphate (1.0mM) (7-DGTPPaS)	80-3321-01	0.1 mL	143.00	157.00	132.00	-	
8,5'-Cyclo-dA CE Phosphoramidite	10-1098-95	50 µmol	1,360.00	1,496.00	1,251.00	-	
8-Amino-dA-CE Phosphoramidite	10-1086-02	0.25 g	1,638.00	1,802.00	1,507.00	-	
8-Amino-dA-CE Phosphoramidite	10-1086-90	100 µmol	577.00	635.00	531.00	-	
8-Amino-dA-CE Phosphoramidite	10-1086-95	50 µmol	288.00	317.00	265.00	-	
8-Amino-dG-CE Phosphoramidite	10-1079-02	0.25 g	1,278.00	1,406.00	1,176.00	-	
8-Amino-dG-CE Phosphoramidite	10-1079-90	100 µmol	465.00	512.00	428.00	-	
8-Amino-dG-CE Phosphoramidite	10-1079-95	50 µmol	233.00	256.00	214.00	-	
8-Aza-7-deaza-A-CE Phosphoramidite	10-3083-02	0.25 g	2,147.00	2,362.00	1,975.00	-	
8-Aza-7-deaza-A-CE Phosphoramidite	10-3083-90	100 µmol	859.00	945.00	790.00	-	
8-Aza-7-deaza-A-CE Phosphoramidite	10-3083-95	50 µmol	429.00	472.00	395.00	-	
8-Br-dA-CE Phosphoramidite	10-1007-02	0.25 g	387.00	426.00	356.00	-	
8-Br-dA-CE Phosphoramidite	10-1007-90	100 µmol	151.00	166.00	139.00	-	
8-Br-dG-CE Phosphoramidite	10-1027-02	0.25 g	334.00	367.00	307.00	-	
8-Br-dG-CE Phosphoramidite	10-1027-90	100 µmol	138.00	152.00	127.00	-	
8-Br-dG-CE Phosphoramidite, Expedite	10-1027-02E	0.25 g	334.00	367.00	307.00	-	
8-OMe-dG-CE Phosphoramidite	10-1075-02	0.25 g	1,396.00	1,536.00	1,284.00	-	
8-OMe-dG-CE Phosphoramidite	10-1075-90	100 µmol	508.00	559.00	467.00	-	
8-OMe-dG-CE Phosphoramidite	10-1075-95	50 µmol	254.00	279.00	234.00	-	
8-oxo-dA-CE Phosphoramidite	10-1008-02	0.25 g	465.00	512.00	428.00	-	
8-oxo-dA-CE Phosphoramidite	10-1008-90	100 µmol	177.00	195.00	163.00	-	
8-oxo-dG-CE Phosphoramidite	10-1028-02	0.25 g	1,278.00	1,406.00	1,176.00	-	
8-oxo-dG-CE Phosphoramidite	10-1028-90	100 µmol	465.00	512.00	428.00	-	
8-oxo-dG-CE Phosphoramidite	10-1028-95	50 µmol	233.00	256.00	214.00	-	
ABI (240mL capacity)	40-4036-A1	each	7.70	8.47	7.08	-	
ABI (450mL capacity)	40-4036-A2	each	7.70	8.47	7.08	-	
Ac-5-Me-dC-CE Phosphoramidite	10-1560-02	0.25 g	157.00	173.00	144.00	-	
Ac-5-Me-dC-CE Phosphoramidite	10-1560-90	100 µmol	65.50	72.10	60.30	-	
Ac-C-CE Phosphoramidite, Pharmacia	10-3015-05P	0.5 g	48.40	53.20	44.50	-	
Ac-dC-CE Phosphoramidite	10-1015-5S	5 g	70.60	77.70	65.00	-	
Ac-dC-CPG 500	20-2013-10	1 g	48.40	53.20	44.50	-	
Ac-dC-Me Phosphoramidite	10-1115-02	0.25 g	295.00	325.00	271.00	-	
Ac-dC-Me Phosphoramidite	10-1115-05	0.5 g	590.00	649.00	543.00	-	
Ac-dC-Me Phosphoramidite	10-1315-02	0.25 g	32.80	36.10	30.20	-	
Ac-dC-Me Phosphoramidite	10-1315-05	0.5 g	65.50	72.10	60.30	-	
Ac-dC-Me Phosphoramidite	10-1315-10	1 g	131.00	144.00	121.00	-	
Ac-dC-Q-CPG 500	21-2013-01	0.1 g	15.10	16.60	13.90	-	
Ac-dC-Q-CPG 500	21-2013-10	1 g	126.00	139.00	116.00	-	
Ac-dC-Q-CPG 500	21-2113-41	4 x 1 µmol	48.40	53.20	44.50	-	
Ac-dC-Q-CPG 500	21-2113-42	4 x 0.2 µmol	32.30	35.50	29.70	-	
Ac-dC-Q-CPG 500	21-2213-41	4 x 1 µmol	48.40	53.20	44.50	-	
Ac-dC-Q-CPG 500	21-2213-42	4 x 0.2 µmol	32.30	35.50	29.70	-	
Ac-dC-Q-CPG 500,	21-2013-02	0.25 g	37.80	41.60	34.80	-	
Ac-G-CE Phosphoramidite	10-3025-05	0.5 g	48.40	53.20	44.50	-	
Acridine Phosphoramidite	10-1973-02	0.25 g	885.00	974.00	814.00	-	
Acridine Phosphoramidite	10-1973-90	100 µmol	387.00	426.00	356.00	-	
Acridine Phosphoramidite	10-1973-95	50 µmol	216.00	238.00	199.00	-	
Adenosine a-thiotriphosphate (0.5mM) (ATTPaS)	80-3000-01	0.1 mL	143.00	157.00	132.00	-	
Alkyne-Modifier Serinol Phosphoramidite	10-1992-02	0.25 g	754.00	829.00	694.00	-	
Alkyne-Modifier Serinol Phosphoramidite	10-1992-90	100 µmol	242.00	266.00	223.00	-	
Alkyne-Modifier Serinol Phosphoramidite	10-1992-95	50 µmol	131.00	144.00	121.00	-	
Alkyne-NHS Ester	50-1905-24	23 mg	393.00	432.00	362.00	-	

Description	Reference	Scale/Quantity	Price List 2021				Ice Shipping
			EUR	CHF	GBP	USD	
Alkyne-NHS Ester, 2.3mg	50-1905-23		78.60	86.50	72.30	-	✓
Amino-Modifier C2 dT	10-1037-02	0.25 g	472.00	519.00	434.00	-	
Amino-Modifier C2 dT	10-1037-05	0.5 g	944.00	1,038.00	868.00	-	
Amino-Modifier C2 dT	10-1037-90	100 µmol	236.00	260.00	217.00	-	
Amino-Modifier C6 dA	10-1089-02	0.25 g	596.00	656.00	548.00	-	
Amino-Modifier C6 dA	10-1089-90	100 µmol	269.00	296.00	247.00	-	
Amino-Modifier C6 dC	10-1019-02	0.25 g	721.00	793.00	663.00	-	
Amino-Modifier C6 dC	10-1019-90	100 µmol	308.00	339.00	283.00	-	
Amino-Modifier C6 dT	10-1039-02	0.25 g	327.00	360.00	301.00	-	
Amino-Modifier C6 dT	10-1039-05	0.5 g	653.00	718.00	601.00	-	
Amino-Modifier C6 dT	10-1039-90	100 µmol	163.00	179.00	150.00	-	
Amino-Modifier C6 dT, Expedite	10-1039-90E	100 µmol	163.00	179.00	150.00	-	
Amino-Modifier C6-U Phosphoramidite	10-3039-02	0.25 g	2,182.00	2,400.00	2,007.00	-	
Amino-Modifier C6-U Phosphoramidite	10-3039-90	100 µmol	1,075.00	1,183.00	989.00	-	
Amino-Modifier C6-U Phosphoramidite	10-3039-95	50 µmol	550.00	605.00	506.00	-	
Amino-Modifier Serinol Phosphoramidite	10-1997-02	0.25 g	780.00	858.00	718.00	-	
Amino-Modifier Serinol Phosphoramidite	10-1997-90	100 µmol	295.00	325.00	271.00	-	
Amino-Modifier Serinol Phosphoramidite	10-1997-95	50 µmol	164.00	180.00	151.00	-	
AP-dC-CE Phosphoramidite	10-1097-02	0.25 g	1,540.00	1,694.00	1,417.00	-	
AP-dC-CE Phosphoramidite	10-1097-90	100 µmol	603.00	663.00	555.00	-	
AP-dC-CE Phosphoramidite	10-1097-95	50 µmol	301.00	331.00	277.00	-	
AP-dC-CE Phosphoramidite, Mermade	10-1097-02M	0.25 g	1,540.00	1,694.00	1,417.00	-	
Ara-C-CE Phosphoramidite	10-4010-02	0.25 g	508.00	559.00	467.00	-	
Ara-C-CE Phosphoramidite	10-4010-90	100 µmol	193.00	212.00	178.00	-	
Asymmetric Doubler (Lev) Phosphoramidite, Expedite	10-1981-90E	100 µmol	138.00	152.00	127.00	-	
Asymmetric Doubler Phosphoramidite	10-1921-02	0.25 g	429.00	472.00	395.00	-	
Asymmetric Doubler Phosphoramidite	10-1921-90	100 µmol	258.00	284.00	237.00	-	
a-Tocopherol-TEG Phosphoramidite	10-1977-02	0.25 g	754.00	829.00	694.00	-	✓
a-Tocopherol-TEG Phosphoramidite	10-1977-90	100 µmol	393.00	432.00	362.00	-	✓
a-Tocopherol-TEG Phosphoramidite	10-1977-95	50 µmol	210.00	231.00	193.00	-	✓
Azidobutyrate NHS Ester	50-1904-23	2.3 mg	78.60	86.50	72.30	-	
Azidobutyrate NHS Ester	50-1904-24	23 mg	393.00	432.00	362.00	-	
Azobenzene Phosphoramidite	10-5800-02	0.25 g	721.00	793.00	663.00	-	✓
Azobenzene Phosphoramidite	10-5800-90	100 µmol	262.00	288.00	241.00	-	✓
Azobenzene Phosphoramidite	10-5800-95	50 µmol	138.00	152.00	127.00	-	
Azobenzene Phosphoramidite, Expedite	10-5800-02E	0.25 g	721.00	793.00	663.00	-	✓
baseclick Oligo-Click-M-Reload	50-2100-01	1 kit	168.00	185.00	155.00	-	
BBQ-650® CPG	20-5934-01	0.1 g	360.00	396.00	331.00	-	✓
BBQ-650® CPG, Expedite	20-5934-41E	4 x 1 µmol	406.00	447.00	374.00	-	
BBQ-650®-dT-CE Phosphoramidite	10-5944-02	0.25 g	1,605.00	1,766.00	1,477.00	-	
BBQ-650®-dT-CE Phosphoramidite	10-5944-90	100 µmol	950.00	1,045.00	874.00	-	
Beckman Oligo 1000 (30mL capacity)	40-4036-B1	each	7.70	8.47	7.08	-	
BHQ-1-dT	10-5941-02	0.25 g	1,605.00	1,766.00	1,477.00	-	
BHQ-1-dT	10-5941-90	100 µmol	924.00	1,016.00	850.00	-	
BHQ-1-dT	10-5941-95	50 µmol	475.00	523.00	437.00	-	
BHQ-1-dT, Mermade	10-5941-02M	0.25 g	1,605.00	1,766.00	1,477.00	-	
BHQ-2-dT	10-5942-02	0.25 g	1,605.00	1,766.00	1,477.00	-	
BHQ-2-dT	10-5942-90	100 µmol	930.00	1,023.00	856.00	-	
BHQ-2-dT	10-5942-95	50 µmol	478.00	526.00	440.00	-	
BHQ-2-dT, Mermade	10-5942-02M	0.25 g	1,605.00	1,766.00	1,477.00	-	
Biosearch 8000 (100mL capacity)	40-4036-C1	each	11.60	12.80	10.70	-	
Biosearch 8000 (500mL capacity)	40-4036-C2	each	11.60	12.80	10.70	-	
BioTEG-CPG -0.2µmol col.(1pk/4)Mermade	20-2955-42M	0.2µmol	121.00	133.00	111.00	-	
BioTEG-CPG, column	20-2955-41E	1 µmol	222.00	244.00	204.00	-	
BioTEG-CPG-1.0 µmol col.(1pk/4)Mermade	20-2955-41M	1 µmol	222.00	244.00	204.00	-	
Biotin PaTP (10mM)	81-3525-02	0.025 mL	644.00	708.00	592.00	-	
Biotin Phosphoramidite	10-1953-02	0.25 g	885.00	974.00	814.00	-	
Biotin Phosphoramidite	10-1953-90	100 µmol	387.00	426.00	356.00	-	
Biotin Phosphoramidite	10-1953-95	50 µmol	216.00	238.00	199.00	-	
Biotin Phosphoramidite, Mermade	10-1953-02M	0.25 g	885.00	974.00	814.00	-	
Biotin-dT	10-1038-02	0.25 g	756.00	832.00	696.00	-	
Biotin-dT	10-1038-90	100 µmol	393.00	432.00	362.00	-	
Biotin-dT	10-1038-95	50 µmol	203.00	223.00	187.00	-	
BiotinTEG Azide	50-2000-90	100 µmol	590.00	649.00	543.00	-	✓
BiotinTEG Azide	50-2000-92	25 µmol	197.00	217.00	181.00	-	✓
BiotinTEG Phosphoramidite	10-1955-02	0.25 g	476.00	524.00	438.00	-	
BiotinTEG Phosphoramidite	10-1955-90	100 µmol	208.00	229.00	191.00	-	
BiotinTEG Phosphoramidite	10-1955-95	50 µmol	116.00	128.00	107.00	-	
BiotinTEG Phosphoramidite, Mermade	10-1955-02M	0.25 g	476.00	524.00	438.00	-	
Bz-5-Me-C-L A-CE Phosphoramidite, Expedite	10-2011-10E	1 g	197.00	217.00	181.00	-	
Bz-5-Me-C-LA-CE Phosphoramidite	10-2011-10	1 g	197.00	217.00	181.00	-	
Bz-A-LA Phosphoramidite	10-2000-05	0.5 g	98.30	108.10	90.40	-	
Bz-A-LA-CE Phosphoramidite	10-2000-10	1 g	197.00	217.00	181.00	-	
Bz-A-LA-CE Phosphoramidite, Expedite	10-2000-10E	1 g	197.00	217.00	181.00	-	
C8-Alkyne-dT-CE Phosphoramidite	10-1540-90	100 µmol	413.00	454.00	380.00	-	
C8-TMS-Alkyne-dT-CE Phosphoramidite	10-1545-90	100 µmol	518.00	570.00	477.00	-	
Carboxy-dT	10-1035-02	0.25 g	472.00	519.00	434.00	-	
Carboxy-dT	10-1035-90	100 µmol	236.00	260.00	217.00	-	
CDPI3 MGB™ CPG, Expedite	20-5924-01	0.1 g	282.00	310.00	259.00	-	
Chemical Phosphorylation Reagent	10-1900-02	0.25 g	161.00	177.00	148.00	-	
Chemical Phosphorylation Reagent	10-1900-90	100 µmol	45.40	49.90	41.80	-	
Chemical Phosphorylation Reagent II	10-1901-02	0.25 g	202.00	222.00	186.00	-	
Chemical Phosphorylation Reagent II	10-1901-90	100 µmol	60.50	66.60	55.70	-	
Cholesteryl-TEG Phosphoramidite	10-1975-02	0.25 g	714.00	785.00	657.00	-	
Cholesteryl-TEG Phosphoramidite	10-1975-90	100 µmol	347.00	382.00	319.00	-	
Cholesteryl-TEG Phosphoramidite	10-1975-95	50 µmol	183.00	201.00	168.00	-	
Cis-syn Thymine Dimer Phosphoramidite	11-1330-02	0.25 g	16,381.00	18,019.00	15,071.00	-	✓
Cis-syn Thymine Dimer Phosphoramidite	11-1330-90	100 µmol	6,736.00	7,410.00	6,197.00	-	✓
Cis-syn Thymine Dimer Phosphoramidite	11-1330-95	50 µmol	3,381.00	3,719.00	3,111.00	-	✓
Colonnes MerMade 40 A	20-2001-65	200 x 50 nmol	605.00	666.00	557.00	-	
Colonnes MerMade 40 C	20-2015-65	200 x 50 nmol	605.00	666.00	557.00	-	

Description	Reference	Scale/Quantity	Price List 2021				Ice Shipping
			EUR	CHF	GBP	USD	
Colonne MerMade 40 G	20-2021-65	200 x 50 nmol	605.00	666.00	557.00	-	
Colonne MerMade 40 T	20-2031-65	200 x 50 nmol	605.00	666.00	557.00	-	
Coumarin Azide	50-2004-90	100 µmol	459.00	505.00	422.00	-	✓
Cy3 Phosphoramidite, Cyclone	10-5913-95C	50 µmol	269.00	296.00	247.00	-	
Cy3 Phosphoramidite, Expedite	10-5913-95E	50 µmol	269.00	296.00	247.00	-	✓
Cy3.5* Phosphoramidite	10-5914-02	0.25 g	1,305.00	1,436.00	1,201.00	-	
Cy3.5* Phosphoramidite	10-5914-90	100 µmol	565.00	622.00	520.00	-	
Cy3.5* Phosphoramidite	10-5914-95	50 µmol	310.00	341.00	285.00	-	
Cy3* Phosphoramidite	10-5913-02	0.25 g	1,212.00	1,333.00	1,115.00	-	
Cy3* Phosphoramidite	10-5913-90	100 µmol	491.00	540.00	452.00	-	
Cy3* Phosphoramidite	10-5913-95	50 µmol	269.00	296.00	247.00	-	
Cy5 Phosphoramidite, Pharmacia	10-5915-95P	50 µmol	269.00	296.00	247.00	-	
Cy5.5* Phosphoramidite	10-5916-02	0.25 g	1,305.00	1,436.00	1,201.00	-	
Cy5.5* Phosphoramidite	10-5916-90	100 µmol	635.00	699.00	584.00	-	
Cy5.5* Phosphoramidite	10-5916-90	100 µmol	635.00	699.00	584.00	-	
Cy5.5* Phosphoramidite	10-5916-95	50 µmol	346.00	381.00	318.00	-	
Cy5* Phosphoramidite	10-5915-02	0.25 g	1,212.00	1,333.00	1,115.00	-	
Cy5* Phosphoramidite	10-5915-90	100 µmol	491.00	540.00	452.00	-	
Cy5* Phosphoramidite	10-5915-95	50 µmol	269.00	296.00	247.00	-	
Cyanine 3 CPG	20-5913-42	4 x 0.2 µmol	100.00	110.00	92.00	-	
Cyanine 5 Phosphoramidite, Expedite	10-5915-95E	50 µmol	269.00	296.00	247.00	-	
Cytidine a-thiotriphosphate (1.0mM) (CTTPaS)	80-3010-01	0.1 mL	143.00	157.00	132.00	-	
dA+dC+dG+dT-CE Phosphoramidite Mix	10-1023-02	0.25 g	26.20	28.80	24.10	-	
dA+dG-CE Phosphoramidite	10-1002-02	0.25 g	26.20	28.80	24.10	-	
dA-5'-CE Phosphoramidite	10-0001-02	0.25 g	83.20	91.50	76.50	-	
dA-5'-CE Phosphoramidite	10-0001-05	0.5 g	166.00	183.00	153.00	-	
dA-5'-CE Phosphoramidite	10-0001-10	1 g	333.00	366.00	306.00	-	
dA-5'-CE Phosphoramidite, Expedite	10-0001-05E	0.5 g	166.00	183.00	153.00	-	
dA-5'-CPG	20-0002-01	0.1 g	65.50	72.10	60.30	-	
dA-5'-CPG	20-0002-10	1 g	491.00	540.00	452.00	-	
dA-5'-CPG	20-0012-13	1 x 10 µmol	295.00	325.00	271.00	-	
dA-5'-CPG	20-0012-14	1 x 15 µmol	393.00	432.00	362.00	-	
dA-5'-CPG	20-0012-41	4 x 1 µmol	131.00	144.00	121.00	-	
dA-5'-CPG	20-0012-42	4 x 0.2 µmol	98.30	108.10	90.40	-	
Dabcyl-dT	10-1058-02	0.25 g	885.00	974.00	814.00	-	
Dabcyl-dT	10-1058-90	100 µmol	426.00	469.00	392.00	-	
Dabcyl-dT	10-1058-95	50 µmol	236.00	260.00	217.00	-	
dA-CE Phosphoramidite, Expedite	10-1000-CSE	0.5 g	7.06	7.77	6.50	-	
dA-Me Phosphoramidite	10-1100-02	0.25 g	295.00	325.00	271.00	-	
dA-Me Phosphoramidite	10-1100-05	0.5 g	590.00	649.00	543.00	-	
dA-Q-CPG 500	21-2000-01	0.1 g	15.10	16.60	13.90	-	
dA-Q-CPG 500	21-2000-10	1 g	136.00	150.00	125.00	-	
dA-Q-CPG 500	21-2100-41	4 x 1 µmol	48.40	53.20	44.50	-	
dA-Q-CPG 500	21-2100-42	4 x 0.2 µmol	32.30	35.50	29.70	-	
dA-Q-CPG 500	21-2200-41	4 x 1 µmol	48.40	53.20	44.50	-	
dA-Q-CPG 500	21-2200-42	4 x 0.2 µmol	32.30	35.50	29.70	-	
dA-Q-CPG 500,	21-2000-02	0.25 g	37.80	41.60	34.80	-	
DBCO TEG	10-1941-02		1,016.00	1,118.00	935.00	-	
dbco nhs ester	50-1941-23		78.60	86.50	72.30	-	
DBCO-dT-CE Phosphoramidite	10-1539-90	100 µmol	636.00	700.00	585.00	-	
DBCO-dT-CE phosphoramidite	10-1539-95	50 µmol	328.00	361.00	302.00	-	
DBCO-dT-CE Phosphoramidite, Expedite	10-1539-95E	50 µmol	328.00	361.00	302.00	-	
DBCO-PEG4-phosphoramidite	10-1941-90	100 µmol	301.00	331.00	277.00	-	
DBCO-Serinol Phosphoramidite, Mermade	10-1998-02M	0.25 g	1,245.00	1,370.00	1,145.00	-	
dC+dT-CE Phosphoramidite	10-1013-02	0.25 g	26.20	28.80	24.10	-	
dC-5'-CE Phosphoramidite	10-0101-02	0.25 g	83.20	91.50	76.50	-	
dC-5'-CE Phosphoramidite	10-0101-05	0.5 g	166.00	183.00	153.00	-	
dC-5'-CE Phosphoramidite	10-0101-10	1 g	333.00	366.00	306.00	-	
dC-5'-CE Phosphoramidite, Expedite	10-0101-02E	0.25 g	83.20	91.50	76.50	-	
dC-5'-CE Phosphoramidite, Expedite	10-0101-05E	0.5 g	166.00	183.00	153.00	-	
dC-5'-CPG	20-0102-01	0.1 g	65.50	72.10	60.30	-	
dC-5'-CPG	20-0102-10	1 g	491.00	540.00	452.00	-	
dC-5'-CPG	20-0112-13	1 x 10 µmol	295.00	325.00	271.00	-	
dC-5'-CPG	20-0112-14	1 x 15 µmol	393.00	432.00	362.00	-	
dC-5'-CPG	20-0112-41	4 x 1 µmol	131.00	144.00	121.00	-	
dC-5'-CPG	20-0112-42	4 x 0.2 µmol	98.30	108.10	90.40	-	
dC-Q-CPG 500	21-2010-01	0.1 g	15.10	16.60	13.90	-	
dC-Q-CPG 500	21-2010-10	1 g	126.00	139.00	116.00	-	
dC-Q-CPG 500	21-2110-41	4 x 1 µmol	48.40	53.20	44.50	-	
dC-Q-CPG 500	21-2110-42	4 x 0.2 µmol	32.30	35.50	29.70	-	
dC-Q-CPG 500	21-2210-41	4 x 1 µmol	48.40	53.20	44.50	-	
dC-Q-CPG 500	21-2210-42	4 x 0.2 µmol	32.30	35.50	29.70	-	
dC-Q-CPG 500,	21-2010-02	0.25 g	37.80	41.60	34.80	-	
def-dA-CE Phosphoramidite	10-1504-10	1 g	78.60	86.50	72.30	-	
def-dA-CE Phosphoramidite, Expedite	10-1504-10E	1 g	78.60	86.50	72.30	-	
DesthiobiotinTEG Azide	50-2001-90	100 µmol	524.00	576.00	482.00	-	✓
DesthiobiotinTEG Azide	50-2001-92	25 µmol	177.00	195.00	163.00	-	✓
DesthiobiotinTEG Phosphoramidite	10-1952-02	0.25 g	1,016.00	1,118.00	935.00	-	
DesthiobiotinTEG Phosphoramidite	10-1952-90	100 µmol	439.00	483.00	404.00	-	
DesthiobiotinTEG Phosphoramidite	10-1952-95	50 µmol	242.00	266.00	223.00	-	
DesthiobiotinTEG-CPG	20-2952-01	0.1 g	183.00	201.00	168.00	-	
DesthiobiotinTEG-CPG	20-2952-10	1 g	1,507.00	1,658.00	1,386.00	-	
DesthiobiotinTEG-CPG	20-2952-13	1 x 10 µmol	452.00	497.00	416.00	-	
DesthiobiotinTEG-CPG	20-2952-14	1 x 15 µmol	681.00	749.00	627.00	-	
DesthiobiotinTEG-CPG	20-2952-41	4 x 1 µmol	301.00	331.00	277.00	-	
DesthiobiotinTEG-CPG	20-2952-42	4 x 0.2 µmol	183.00	201.00	168.00	-	
dG-5'-CPG	20-0202-01	0.1 g	65.50	72.10	60.30	-	
dG-5'-CPG	20-0202-10	1 g	491.00	540.00	452.00	-	
dG-5'-CPG	20-0212-13	1 x 10 µmol	295.00	325.00	271.00	-	
dG-5'-CPG	20-0212-14	1 x 15 µmol	393.00	432.00	362.00	-	
dG-5'-CPG	20-0212-41	4 x 1 µmol	131.00	144.00	121.00	-	



Description	Reference	Scale/Quantity	Price List 2021				Ice Shipping
			EUR	CHF	GBP	USD	
dG-5'-CPG	20-0212-42	4 x 0.2 µmol	98.30	108.10	90.40	-	
dG-CE Phosphoramidite	10-1020-C5E	0.5 g	7.06	7.77	6.50	-	
dG-Me Phosphoramidite	10-1120-02	0.25 g	295.00	325.00	271.00	-	
dG-Me Phosphoramidite	10-1120-05	0.5 g	590.00	649.00	543.00	-	
dl-CE Phosphoramidite	10-1040-02	0.25 g	109.00	120.00	100.00	-	
dl-CE Phosphoramidite	10-1040-90	100 µmol	45.40	49.90	41.80	-	
dl-CPG 1000	20-2041-01	0.1 g	39.30	43.20	36.20	-	
dl-CPG 1000	20-2191-41	4 x 1 µmol	157.00	173.00	144.00	-	
dl-CPG 1000	20-2191-42	4 x 0.2 µmol	103.00	113.00	95.00	-	
dl-CPG 500	20-2040-01	0.1 g	39.30	43.20	36.20	-	
dl-CPG 500	20-2190-41	4 x 1 µmol	157.00	173.00	144.00	-	
dl-CPG 500	20-2190-42	4 x 0.2 µmol	94.40	103.80	86.80	-	
Dithiol Serinol Phosphoramidite	10-1991-02	0.25 g	767.00	844.00	706.00	-	
Dithiol Serinol Phosphoramidite	10-1991-90	100 µmol	282.00	310.00	259.00	-	
dK-CE Phosphoramidite	10-1048-02	0.25 g	780.00	858.00	718.00	-	
dK-CE Phosphoramidite	10-1048-90	100 µmol	256.00	282.00	236.00	-	
dmf-5-Me-isodC-CE Phosphoramidite	10-1065-02	0.25 g	454.00	499.00	418.00	-	
dmf-5-Me-isodC-CE Phosphoramidite	10-1065-90	100 µmol	145.00	160.00	133.00	-	
dmf-5-Me-isodC-CE Phosphoramidite, Expedite	10-1065-02E	0.25 g	454.00	499.00	418.00	-	
dmf-dG PS	26-2629-62	200 x 200 nmol	650.00	715.00	598.00	-	
dmf-dG PS	26-2629-65	200 x 40 nmol	650.00	715.00	598.00	-	
dmf-dG-5'-CE Phosphoramidite	10-9201-02	0.25 g	83.20	91.50	76.50	-	
dmf-dG-5'-CE Phosphoramidite	10-9201-05	0.5 g	166.00	183.00	153.00	-	
dmf-dG-5'-CE Phosphoramidite	10-9201-10	1 g	333.00	366.00	306.00	-	
dmf-dG-CE Phosphoramidite	10-1029-02	0.25 g	3.53	3.88	3.25	-	
dmf-dG-CE Phosphoramidite	10-1029-05	0.5 g	7.06	7.77	6.50	-	
dmf-dG-CE Phosphoramidite	10-1029-10	1 g	14.10	15.50	13.00	-	
dmf-dG-CE Phosphoramidite	10-1029-18	1 g	15.40	16.90	14.20	-	
dmf-dG-CE Phosphoramidite	10-1029-1C	1 g	14.10	15.50	13.00	-	
dmf-dG-CE Phosphoramidite	10-1029-1P	1 g	15.40	16.90	14.20	-	
dmf-dG-CE Phosphoramidite	10-1029-1S	10 g	141.00	155.00	130.00	-	
dmf-dG-CE Phosphoramidite	10-1029-20	2 g	28.20	31.00	25.90	-	
dmf-dG-CE Phosphoramidite	10-1029-50	5 g	70.60	77.70	65.00	-	
dmf-dG-CE Phosphoramidite	10-1029-5S	5 g	70.60	77.70	65.00	-	
dmf-dG-CE Phosphoramidite	10-1029-B2	0.25 g	3.90	4.29	3.59	-	
dmf-dG-CE Phosphoramidite	10-1029-B5	0.5 g	7.70	8.47	7.08	-	
dmf-dG-CE Phosphoramidite	10-1029-C2	0.25 g	3.53	3.88	3.25	-	
dmf-dG-CE Phosphoramidite	10-1029-C5	0.5 g	7.06	7.77	6.50	-	
dmf-dG-CE Phosphoramidite	10-1029-P4	0.4 g	6.20	6.82	5.70	-	
dmf-dG-CE Phosphoramidite, Expedite	10-1029-02E	0.25 g	3.53	3.88	3.25	-	
dmf-dG-CE Phosphoramidite, Expedite	10-1029-C5E	0.5 g	7.06	7.77	6.50	-	
dmf-dG-CPG	20-2029-01	0.1 g	5.81	6.39	5.35	-	
dmf-dG-CPG	20-2029-02	0.25 g	12.90	14.20	11.90	-	
dmf-dG-CPG	20-2029-10	1 g	48.40	53.20	44.50	-	
dmf-dG-CPG	20-2129-13	1 x 10 µmol	80.60	88.70	74.20	-	
dmf-dG-CPG	20-2129-41	4 x 1 µmol	28.20	31.00	25.90	-	
dmf-dG-CPG	20-2129-42	4 x 0.2 µmol	28.20	31.00	25.90	-	
dmf-dG-CPG	20-2129-45	4 x 40 nmol	28.20	31.00	25.90	-	
dmf-dG-CPG	20-2229-14	1 x 15 µmol	121.00	133.00	111.00	-	
dmf-dG-CPG	20-2229-41	4 x 1 µmol	20.20	22.20	18.60	-	
dmf-dG-CPG	20-2229-42	4 x 0.2 µmol	14.10	15.50	13.00	-	
dmf-dG-CPG	20-2229-45	4 x 40 nmol	14.10	15.50	13.00	-	
dmf-dG-Q-CPG 500	21-2029-01	0.1 g	25.20	27.70	23.20	-	
dmf-dG-Q-CPG 500	21-2029-10	1 g	222.00	244.00	204.00	-	
dmf-dG-Q-CPG 500	21-2129-41	4 x 1 µmol	48.40	53.20	44.50	-	
dmf-dG-Q-CPG 500	21-2129-42	4 x 0.2 µmol	32.30	35.50	29.70	-	
dmf-dG-Q-CPG 500	21-2229-41	4 x 1 µmol	48.40	53.20	44.50	-	
dmf-dG-Q-CPG 500	21-2229-42	4 x 0.2 µmol	32.30	35.50	29.70	-	
dmf-dG-Q-CPG 500,	21-2029-02	0.25 g	55.40	60.90	51.00	-	
dmf-G-LA Phosphoramidite	10-2029-05	0.5 g	98.30	108.10	90.40	-	
dmf-G-LA-CE Phosphoramidite	10-2029-10	1 g	197.00	217.00	181.00	-	
dmf-G-LA-CE Phosphoramidite, Expedite	10-2029-10E	1 g	197.00	217.00	181.00	-	
dmf-isodG-CE Phosphoramidite	10-1078-02	0.25 g	394.00	433.00	362.00	-	
dmf-isodG-CE Phosphoramidite	10-1078-90	100 µmol	183.00	201.00	168.00	-	
DNP-TEG Phosphoramidite	10-1985-02	0.25 g	885.00	974.00	814.00	-	
DNP-TEG Phosphoramidite	10-1985-90	100 µmol	387.00	426.00	356.00	-	
DNP-TEG Phosphoramidite	10-1985-95	50 µmol	216.00	238.00	199.00	-	
dP+dK-CE Phosphoramidite	10-1049-02	0.25 g	780.00	858.00	718.00	-	
dP+dK-CE Phosphoramidite	10-1049-90	100 µmol	256.00	282.00	236.00	-	
dP-CE Phosphoramidite	10-1047-02	0.25 g	780.00	858.00	718.00	-	
dP-CE Phosphoramidite	10-1047-90	100 µmol	256.00	282.00	236.00	-	
dSpacer CE Phosphoramidite	10-1914-02	0.25 g	378.00	416.00	348.00	469.00	
dSpacer CE Phosphoramidite	10-1914-90	100 µmol	116.00	128.00	107.00	144.00	
dSpacer CE Phosphoramidite, Expedite	10-1914-02E	0.25 g	378.00	416.00	348.00	-	
dT-5'-CE Phosphoramidite	10-0301-02	0.25 g	83.20	91.50	76.50	-	
dT-5'-CE Phosphoramidite	10-0301-05	0.5 g	166.00	183.00	153.00	-	
dT-5'-CE Phosphoramidite	10-0301-10	1 g	333.00	366.00	306.00	-	
dT-5'-CE Phosphoramidite, Expedite	10-0301-10E	1 g	333.00	366.00	306.00	-	
dT-5'-CPG	20-0302-01	0.1 g	65.50	72.10	60.30	-	
dT-5'-CPG	20-0302-10	1 g	491.00	540.00	452.00	-	
dT-5'-CPG	20-0312-13	1 x 10 µmol	295.00	325.00	271.00	-	
dT-5'-CPG	20-0312-14	1 x 15 µmol	393.00	432.00	362.00	-	
dT-5'-CPG	20-0312-41	4 x 1 µmol	131.00	144.00	121.00	-	
dT-5'-CPG	20-0312-42	4 x 0.2 µmol	98.30	108.10	90.40	-	
dT-5'-Icaa-CPG 500, column	20-0312-41E	4 x 1 µmol	131.00	144.00	121.00	-	
dT-Me Phosphoramidite	10-1130-02	0.25 g	295.00	325.00	271.00	-	
dT-Me Phosphoramidite	10-1130-05	0.5 g	590.00	649.00	543.00	-	
dT-Me Phosphoramidite	10-1330-02	0.25 g	32.80	36.10	30.20	-	
dT-Me Phosphoramidite	10-1330-05	0.5 g	65.50	72.10	60.30	-	
dT-Me Phosphoramidite	10-1330-10	1 g	131.00	144.00	121.00	-	
dT-PS	26-2630-61	200 x 1 µmol	1,244.00	1,368.00	1,144.00	-	

Description	Reference	Scale/Quantity	Price List 2021				Ice Shipping
			EUR	CHF	GBP	USD	
dT-Q-CPG 500	21-2030-01	0.1 g	15.10	16.60	13.90	-	
dT-Q-CPG 500	21-2030-10	1 g	126.00	139.00	116.00	-	
dT-Q-CPG 500	21-2130-41	4 x 1 µmol	48.40	53.20	44.50	-	
dT-Q-CPG 500	21-2130-42	4 x 0.2 µmol	32.30	35.50	29.70	-	
dT-Q-CPG 500	21-2230-41	4 x 1 µmol	48.40	53.20	44.50	-	
dT-Q-CPG 500,	21-2030-02	0.25 g	37.80	41.60	34.80	-	
dU-CE Phosphoramidite	10-1050-02	0.25 g	80.60	88.70	74.20	-	
dU-CE Phosphoramidite	10-1050-90	100 µmol	28.20	31.00	25.90	-	
dU-CPG 1000	20-2051-01	0.1 g	65.50	72.10	60.30	-	
dU-CPG 1000	20-2151-41	4 x 1 µmol	262.00	288.00	241.00	-	
dU-CPG 1000	20-2151-42	4 x 0.2 µmol	157.00	173.00	144.00	-	
dU-CPG 500	20-2050-01	0.1 g	39.30	43.20	36.20	-	
dU-CPG 500	20-2150-41	4 x 1 µmol	157.00	173.00	144.00	-	
dU-CPG 500	20-2150-42	4 x 0.2 µmol	94.40	103.80	86.80	-	
dX-CE Phosphoramidite	10-1537-02	0.25 g	917.00	1,009.00	844.00	-	
dX-CE Phosphoramidite	10-1537-95	50 µmol	223.00	245.00	205.00	-	
DyLight DY547 Phosphoramidite	10-5917-90	100 µmol	480.00	528.00	442.00	-	✓
DyLight DY647 Phosphoramidite	10-5918-02	0.25 g	1,424.00	1,566.00	1,310.00	-	✓
DyLight DY647 Phosphoramidite	10-5918-90	100 µmol	480.00	528.00	442.00	-	✓
EDTA-C2-dT-CE Phosphoramidite	10-1059-02	0.25 g	1,278.00	1,406.00	1,176.00	-	
EDTA-C2-dT-CE Phosphoramidite	10-1059-90	100 µmol	649.00	714.00	597.00	-	
EDTA-C2-dT-CE Phosphoramidite	10-1059-95	50 µmol	328.00	361.00	302.00	-	
Empty MerMade Columns (200 nmol and 1 µmol)	20-0050-02	pk / 48	161.00	177.00	148.00	-	
Epoch Eclipse QuencherCPG0.2µm(1pk/4)Mer	20-5925-42M	4 x 0.2 µmol	212.00	233.00	195.00	-	
Epoch Eclipse QuencherCPG1µm(1pk/4)Mer	20-5925-41M	4 x 1 µmol	423.00	465.00	389.00	-	
Epoch Eclipse QuencherCPG1µm(1pk/4)Mer	20-5925-41MD		423.00	465.00	389.00	-	
Epoch Eclipse™ Quencher CPG	20-5925-01	0.1 g	393.00	432.00	362.00	-	
Epoch Eclipse™ Quencher CPG	20-5925-10	1 g	3,276.00	3,604.00	3,014.00	-	
Epoch Eclipse™ Quencher CPG	20-5925-13	1 x 10 µmol	1,304.00	1,434.00	1,200.00	-	
Epoch Eclipse™ Quencher CPG	20-5925-14	1 x 15 µmol	1,959.00	2,155.00	1,802.00	-	
Epoch Eclipse™ Quencher CPG	20-5925-41	4 x 1 µmol	423.00	465.00	389.00	-	
Epoch Eclipse™ Quencher CPG	20-5925-42	4 x 0.2 µmol	212.00	233.00	195.00	-	
Epoch Eclipse™ Quencher Phosphoramidite	10-5925-02	0.25 g	2,143.00	2,357.00	1,972.00	-	
Epoch Eclipse™ Quencher Phosphoramidite	10-5925-90	100 µmol	865.00	952.00	796.00	-	
Epoch Eclipse™ Quencher Phosphoramidite	10-5925-95	50 µmol	432.00	475.00	397.00	-	
Epoch Gig Harbor Green™ Phosphoramidite	10-5922-02	0.25 g	1,138.00	1,252.00	1,047.00	-	
Epoch Gig Harbor Green™ Phosphoramidite	10-5922-90	100 µmol	427.00	470.00	393.00	-	
Epoch Gig Harbor Green™ Phosphoramidite	10-5922-95	50 µmol	226.00	249.00	208.00	-	
Epoch Redmond Red™ CPG	20-5920-01	0.1 g	282.00	310.00	259.00	-	
Epoch Redmond Red™ CPG	20-5920-10	1 g	2,359.00	2,595.00	2,170.00	-	
Epoch Redmond Red™ CPG	20-5920-13	1 x 10 µmol	983.00	1,081.00	904.00	-	
Epoch Redmond Red™ CPG	20-5920-14	1 x 15 µmol	1,474.00	1,621.00	1,356.00	-	
Epoch Redmond Red™ CPG	20-5920-41	4 x 1 µmol	491.00	540.00	452.00	-	
Epoch Redmond Red™ CPG	20-5920-42	4 x 0.2 µmol	197.00	217.00	181.00	-	
Epoch Redmond Red™ Phosphoramidite	10-5920-02	0.25 g	1,874.00	2,061.00	1,724.00	-	
Epoch Redmond Red™ Phosphoramidite	10-5920-90	100 µmol	754.00	829.00	694.00	-	
Epoch Redmond Red™ Phosphoramidite	10-5920-95	50 µmol	390.00	429.00	359.00	-	
Epoch Yakima Yellow™ CPG	20-5921-01	0.1 g	282.00	310.00	259.00	-	
Epoch Yakima Yellow™ CPG	20-5921-10	1 g	2,359.00	2,595.00	2,170.00	-	
Epoch Yakima Yellow™ CPG	20-5921-13	1 x 10 µmol	983.00	1,081.00	904.00	-	
Epoch Yakima Yellow™ CPG	20-5921-14	1 x 15 µmol	1,474.00	1,621.00	1,356.00	-	
Epoch Yakima Yellow™ CPG	20-5921-41	4 x 1 µmol	491.00	540.00	452.00	-	
Epoch Yakima Yellow™ CPG	20-5921-42	4 x 0.2 µmol	197.00	217.00	181.00	-	
Epoch Yakima Yellow™ Phosphoramidite	10-5921-02	0.25 g	1,881.00	2,069.00	1,731.00	-	
Epoch Yakima Yellow™ Phosphoramidite	10-5921-90	100 µmol	799.00	879.00	735.00	-	
Epoch Yakima Yellow™ Phosphoramidite	10-5921-95	50 µmol	413.00	454.00	380.00	-	
Etheno-dA-CE Phosphoramidite	10-1006-02	0.25 g	347.00	382.00	319.00	-	
Etheno-dA-CE Phosphoramidite	10-1006-90	100 µmol	138.00	152.00	127.00	-	
Etheno-dA-CE Phosphoramidite, Expedite	10-1006-02E	0.25 g	347.00	382.00	319.00	-	
Expedite 8905 (150mL capacity)	40-4036-C3	each	7.70	8.47	7.08	-	
Expedite 8909 and Cyclone (240mL capacity)	40-4036-C4	each	7.70	8.47	7.08	-	
F-DMT-dT-CE Phosphoramidite, Pharmacia	10-1430-90P	100 µmol	107.00	118.00	98.00	-	✓
Ferrocene-dT-CE Phosphoramidite	10-1576-90	100 µmol	459.00	505.00	422.00	-	
Fluorescein DT Phosphoramidite	10-1056-02	0.25 g	680.00	748.00	626.00	-	
Fluorescein Phosphoramidite	10-1963-02	0.25 g	780.00	858.00	718.00	-	
Fluorescein Phosphoramidite	10-1963-90	100 µmol	387.00	426.00	356.00	-	
Fluorescein Phosphoramidite	10-1963-95	50 µmol	216.00	238.00	199.00	-	
Fluorescein-dT Phosphoramidite	10-1056-90	100 µmol	408.00	449.00	375.00	506.00	
Fluorescein-dT Phosphoramidite	10-1056-95	50 µmol	214.00	235.00	197.00	265.00	
Fluorescein-dT Phosphoramidite, Expedite	10-1056-02E	0.25 g	680.00	748.00	626.00	-	
Fluorescein-dT Phosphoramidite, Mermade	10-1056-02M	0.25 g	680.00	748.00	626.00	-	
Fluorescein-dUTP-dabsyl (1mM, 25nmoles, 10mM Tris, 1mM EDTA)	88-1056-01	0.1 mL	386.00	425.00	355.00	-	
Fmoc Amino-Modifier C6 dT	10-1536-02	0.25 g	472.00	519.00	434.00	-	
Fmoc Amino-Modifier C6 dT	10-1536-90	100 µmol	236.00	260.00	217.00	-	
Formyl dC	10-1564-02		3,715.00	4,087.00	3,418.00	-	
GalNAc C3 CPG	20-2974-10	1 g	419.00	461.00	385.00	-	
GalNAc C3 CPG	20-2974-41E	4 x 1 µmol	131.00	144.00	121.00	-	
Glen Gel-Pak™ 0.2 Desalting Column	61-5002-05	Pk / 5	59.00	64.90	54.30	-	
Glen Gel-Pak™ 1.0 Desalting Column	61-5010-05	Pk / 5	59.00	64.90	54.30	-	
Glen UnySupport™ 500, columns	20-5240-41	4 x 1 µmol	53.00	58.30	48.80	-	
Glen UnySupport™ HybridCPG™	28-5040-10	1 g	130.00	143.00	120.00	-	
Glen UnySupport™ PS	26-5040-10	1 g	164.00	180.00	151.00	-	
Glen-Pak™ DNA purification cartridge	60-5100-10	Pk / 10	64.50	71.00	59.30	-	
Glen-Pak™ DNA purification cartridge	60-5100-96	Pk / 96	408.00	449.00	375.00	-	
Glen-Pak™ DNA purification cartridge (for use with disposable syringes)	60-5200-10	Pk / 10	101.00	111.00	93.00	-	
Glen-Pak™ RNA purification cartridge	60-6100-10	Pk / 10	76.60	84.30	70.50	-	
Glen-Pak™ RNA purification cartridge	60-6100-96	Pk / 96	464.00	510.00	427.00	-	
Glen-Pak™ RNA purification cartridge (for use with disposable syringes)	60-6200-10	Pk / 10	126.00	139.00	116.00	-	
Guanosine a-thiotriphosphate (0.5mM) (GTPPaS)	80-3020-01	0.1 mL	143.00	157.00	132.00	-	
High Load Glen UnySupport	25-5040-10	1 g	116.00	128.00	107.00	-	
I-CE Phosphoramidite	10-3040-02	0.25 g	655.00	721.00	603.00	-	

Description	Reference	Scale/Quantity	Price List 2021				Ice Shipping
			EUR	CHF	GBP	USD	
Inosine a-thiotriphosphate (4mM) (ITTPaS)	80-3050-01	0.1 mL	143.00	157.00	132.00	-	
iPr-Pac-dG-CE Phosphoramidite	10-1621-02	0.25 g	18.10	19.90	16.70	-	
iPr-Pac-dG-CE Phosphoramidite	10-1621-05	0.5 g	36.30	39.90	33.40	-	
iPr-Pac-dG-CE Phosphoramidite	10-1621-10	1 g	72.60	79.90	66.80	-	
iPr-Pac-dG-CE Phosphoramidite, Expedite	10-1621-02E	0.25 g	18.10	19.90	16.70	-	
iPr-Pac-dG-CE Phosphoramidite, Expedite	10-1621-05E	0.5 g	36.30	39.90	33.40	-	
iPr-Pac-dG-CPG	20-2621-01	0.1 g	23.60	26.00	21.70	-	
iPr-Pac-dG-CPG	20-2621-02	0.25 g	52.40	57.60	48.20	-	
iPr-Pac-dG-CPG	20-2621-10	1 g	197.00	217.00	181.00	-	
iPr-Pac-dG-CPG	20-2721-13	1 x 10 µmol	131.00	144.00	121.00	-	
iPr-Pac-dG-CPG	20-2721-41	4 x 1 µmol	78.60	86.50	72.30	-	
iPr-Pac-dG-CPG	20-2721-42	4 x 0.2 µmol	52.40	57.60	48.20	-	
iPr-Pac-dG-CPG	20-2721-45	4 x 40 nmol	52.40	57.60	48.20	-	
iPr-Pac-dG-CPG	20-2821-14	1 x 15 µmol	197.00	217.00	181.00	-	
iPr-Pac-dG-CPG	20-2821-41	4 x 1 µmol	78.60	86.50	72.30	-	
iPr-Pac-dG-CPG	20-2821-42	4 x 0.2 µmol	52.40	57.60	48.20	-	
iPr-Pac-dG-CPG	20-2821-45	4 x 40 nmol	52.40	57.60	48.20	-	
iPr-Pac-dG-Me Phosphoramidite	10-1321-02	0.25 g	32.80	36.10	30.20	-	
iPr-Pac-dG-Me Phosphoramidite	10-1321-05	0.5 g	65.50	72.10	60.30	-	
iPr-Pac-dG-Me Phosphoramidite	10-1321-10	1 g	131.00	144.00	121.00	-	
iPr-Pac-G-RNA-CPG	20-3321-10	1 g	465.00	512.00	428.00	-	
Long Trebler Phosphoramidite	10-1925-02	0.25 g	393.00	432.00	362.00	-	
Long Trebler Phosphoramidite	10-1925-90	100 µmol	262.00	288.00	241.00	-	
Long Trebler Phosphoramidite, Expedite	10-1925-02E	0.25 g	393.00	432.00	362.00	-	
Methacrylate C6 Phosphoramidite	10-1891-02	0.25 g	852.00	937.00	784.00	-	
Methacrylate C6 Phosphoramidite	10-1891-90	100 µmol	144.00	158.00	132.00	-	
Methylene Blue C3 Phosphoramidite	10-5960-90	100 µmol	760.00	836.00	699.00	-	
Methylene Blue NHS Ester	50-1960-23	5.4 mg	754.00	829.00	694.00	-	
N2-Amino-Modifier C6 dG	10-1529-02	0.25 g	1,442.00	1,586.00	1,327.00	-	
N2-Amino-Modifier C6 dG	10-1529-90	100 µmol	629.00	692.00	579.00	-	
N3-Cyanoethyl-dT	10-1531-02	0.25 g	786.00	865.00	723.00	-	
N3-Cyanoethyl-dT	10-1531-90	100 µmol	262.00	288.00	241.00	-	
N4-Et-dC-CE Phosphoramidite	10-1068-02	0.25 g	885.00	974.00	814.00	-	
N4-Et-dC-CE Phosphoramidite	10-1068-90	100 µmol	295.00	325.00	271.00	-	
N4-Et-dC-CE Phosphoramidite	10-1068-95	50 µmol	164.00	180.00	151.00	-	
N4-Et-dC-CE Phosphoramidite	10-1068-SP	1 g	1,670.00	1,837.00	1,536.00	-	
N6-Ac-N6-Me-dA-CE Phosphoramidite	10-1503-02	0.25 g	649.00	714.00	597.00	-	
N6-Ac-N6-Me-dA-CE Phosphoramidite	10-1503-90	100 µmol	213.00	234.00	196.00	-	
N6-Me-Adenosine a-thiotriphosphate (4mM) (N6-Me-ATTPaS)	80-3302-01	0.1 mL	143.00	157.00	132.00	-	
N6-Me-dA-CE Phosphoramidite	10-1003-02	0.25 g	649.00	714.00	597.00	-	
N6-Me-dA-CE Phosphoramidite	10-1003-90	100 µmol	213.00	234.00	196.00	-	
N6-Methyl-A-CE Phosphoramidite	10-3005-90	0.25 g	721.00	793.00	663.00	-	
N6-Methyl-A-CE Phosphoramidite	10-3005-95	50 µmol	373.00	410.00	343.00	-	
Nebularine-CE Phosphoramidite	10-3041-02	0.25 g	1,789.00	1,968.00	1,646.00	-	
Nebularine-CE Phosphoramidite	10-3041-90	100 µmol	644.00	708.00	592.00	-	
Nebularine-CE Phosphoramidite	10-3041-95	50 µmol	322.00	354.00	296.00	-	
NHS-Carboxy-dT	10-1535-02	0.25 g	721.00	793.00	663.00	-	
NHS-Carboxy-dT	10-1535-90	100 µmol	275.00	303.00	253.00	-	
N-POM Caged-dT-CE Phosphoramidite	10-1534-02	0.25 g	1,173.00	1,290.00	1,079.00	-	
N-POM Caged-dT-CE Phosphoramidite	10-1534-90	100 µmol	465.00	512.00	428.00	-	
O4-Me-dT-CE Phosphoramidite	10-1032-02	0.25 g	508.00	559.00	467.00	-	
O4-Me-dT-CE Phosphoramidite	10-1032-90	100 µmol	193.00	212.00	178.00	-	
O4-Triazolyl-dU-CE Phosphoramidite	10-1051-02	0.25 g	622.00	684.00	572.00	-	
O4-Triazolyl-dU-CE Phosphoramidite	10-1051-90	100 µmol	197.00	217.00	181.00	-	
O6-Me-dG-CE Phosphoramidite	10-1070-02	0.25 g	334.00	367.00	307.00	-	
O6-Me-dG-CE Phosphoramidite	10-1070-90	100 µmol	138.00	152.00	127.00	-	
O6-Phenyl-dl-CE Phosphoramidite	10-1042-02	0.25 g	465.00	512.00	428.00	-	
O6-Phenyl-dl-CE Phosphoramidite	10-1042-90	100 µmol	177.00	195.00	163.00	-	
Oligo-Affinity Support (PS)	26-4001-01	0.1 g	236.00	260.00	217.00	-	
Oligo-Affinity Support (PS)	26-4001-02	0.25 g	557.00	613.00	512.00	-	
Oligo-Affinity Support (PS)	26-4001-10	1 g	2,084.00	2,292.00	1,917.00	-	
Oligo-Affinity Support (PS)	26-4101-41	4 x 1 µmol	393.00	432.00	362.00	-	
Oligo-Affinity Support (PS), Expedite	26-4101-41E	4 x 1 µmol	393.00	432.00	362.00	-	
Oligonucleotide-Peptide Conjugation Kit (OPeC)	50-1000-01	1 g	709.00	780.00	652.00	-	
Pac-A-RNA-CPG	20-3300-10	1 g	465.00	512.00	428.00	-	
Pac-dA-CE Phosphoramidite	10-1601-02	0.25 g	18.10	19.90	16.70	-	
Pac-dA-CE Phosphoramidite	10-1601-05	0.5 g	36.30	39.90	33.40	-	
Pac-dA-CE Phosphoramidite	10-1601-10	1 g	72.60	79.90	66.80	-	
Pac-dA-CE Phosphoramidite, Expedite	10-1601-02E	0.25 g	18.10	19.90	16.70	-	
Pac-dA-CE Phosphoramidite, Expedite	10-1601-05E	0.5 g	36.30	39.90	33.40	-	
Pac-dA-CPG	20-2601-01	0.1 g	23.60	26.00	21.70	-	
Pac-dA-CPG	20-2601-02	0.25 g	52.40	57.60	48.20	-	
Pac-dA-CPG	20-2601-10	1 g	197.00	217.00	181.00	-	
Pac-dA-CPG	20-2701-13	1 x 10 µmol	131.00	144.00	121.00	-	
Pac-dA-CPG	20-2701-41	4 x 1 µmol	78.60	86.50	72.30	-	
Pac-dA-CPG	20-2701-42	4 x 0.2 µmol	52.40	57.60	48.20	-	
Pac-dA-CPG	20-2701-45	4 x 40 nmol	52.40	57.60	48.20	-	
Pac-dA-CPG	20-2801-14	1 x 15 µmol	197.00	217.00	181.00	-	
Pac-dA-CPG	20-2801-41	4 x 1 µmol	78.60	86.50	72.30	-	
Pac-dA-CPG	20-2801-42	4 x 0.2 µmol	52.40	57.60	48.20	-	
Pac-dA-CPG	20-2801-45	4 x 40 nmol	52.40	57.60	48.20	-	
Pac-dA-Me Phosphoramidite	10-1301-02	0.25 g	32.80	36.10	30.20	-	
Pac-dA-Me Phosphoramidite	10-1301-05	0.5 g	65.50	72.10	60.30	-	
Pac-dA-Me Phosphoramidite	10-1301-10	1 g	131.00	144.00	121.00	-	
PC Amino-Modifier Phosphoramidite	10-4906-02	0.25 g	518.00	570.00	477.00	-	
PC Amino-Modifier Phosphoramidite	10-4906-90	100 µmol	177.00	195.00	163.00	-	
PC Biotin Phosphoramidite	10-4950-02	0.25 g	885.00	974.00	814.00	-	
PC Biotin Phosphoramidite	10-4950-90	100 µmol	367.00	404.00	338.00	-	
PC Biotin Phosphoramidite	10-4950-95	50 µmol	190.00	209.00	175.00	-	
PC Linker Phosphoramidite	10-4920-02	0.25 g	1,042.00	1,146.00	959.00	-	
PC Linker Phosphoramidite	10-4920-90	100 µmol	334.00	367.00	307.00	-	

Description	Reference	Scale/Quantity	Price List 2021				Ice Shipping
			EUR	CHF	GBP	USD	
PC Spacer Phosphoramidite	10-4913-02	0.25 g	518.00	570.00	477.00	-	
PC Spacer Phosphoramidite	10-4913-90	100 µmol	177.00	195.00	163.00	-	
pdC-CE Phosphoramidite	10-1014-02	0.25 g	321.00	353.00	295.00	-	
pdC-CE Phosphoramidite	10-1014-05	0.5 g	642.00	706.00	591.00	-	
pdC-CE Phosphoramidite	10-1014-90	100 µmol	144.00	158.00	132.00	-	
pdC-CE Phosphoramidite, Mermade	10-1014-02M	0.25 g	321.00	353.00	295.00	-	
pdU-CE Phosphoramidite	10-1054-02	0.25 g	256.00	282.00	236.00	-	
pdU-CE Phosphoramidite	10-1054-05	0.5 g	511.00	562.00	470.00	-	
pdU-CE Phosphoramidite, Mermade	10-1054-02M	0.25 g	256.00	282.00	236.00	-	
Perylene-dU	10-1591-02	0.25 g	983.00	1,081.00	904.00	-	
Perylene-dU	10-1591-90	100 µmol	419.00	461.00	385.00	-	
Perylene-dU	10-1591-95	50 µmol	223.00	245.00	205.00	-	
Perylene-dU-CE Phosphoramidite, Expedite	10-1591-90E	100 µmol	419.00	461.00	385.00	-	
Protected Biotin Serinol Phosphoramidite	10-1993-02	0.25 g	885.00	974.00	814.00	-	
Protected Biotin Serinol Phosphoramidite	10-1993-90	100 µmol	387.00	426.00	356.00	-	
Protected Biotin Serinol Phosphoramidite	10-1993-95	50 µmol	216.00	238.00	199.00	-	
Protected BiotinLC Serinol Phosphoramidite	10-1995-02	0.25 g	885.00	974.00	814.00	-	
Protected BiotinLC Serinol Phosphoramidite	10-1995-90	100 µmol	478.00	526.00	440.00	-	
Protected BiotinLC Serinol Phosphoramidite	10-1995-95	50 µmol	269.00	296.00	247.00	-	
Protected BiotinLC Serinol Phosphoramidite, Mermade	10-1995-02M	0.25 g	885.00	974.00	814.00	-	
PseudoUridine-CE Phosphoramidite	10-3055-02	0.25 g	1,428.00	1,571.00	1,314.00	-	
PseudoUridine-CE Phosphoramidite	10-3055-90	100 µmol	518.00	570.00	477.00	-	
PseudoUridine-CE Phosphoramidite	10-3055-95	50 µmol	272.00	299.00	250.00	-	
Psoralen Azide	50-2009-92	25 µmol	151.00	166.00	139.00	-	
Psoralen C2 Phosphoramidite	10-1982-02	0.25 g	649.00	714.00	597.00	-	
Psoralen C2 Phosphoramidite	10-1982-90	100 µmol	256.00	282.00	236.00	-	
Psoralen C6 Phosphoramidite	10-1983-02	0.25 g	649.00	714.00	597.00	-	
Psoralen C6 Phosphoramidite	10-1983-90	100 µmol	256.00	282.00	236.00	-	
Purine riboside a-thiotriphosphate (20 mM) (PR-TTPaS)	80-3301-01	0.1 mL	143.00	157.00	132.00	-	
Puromycin-CPG	20-4040-01	0.1 g	157.00	173.00	144.00	-	
Puromycin-CPG	20-4040-10	1 g	1,304.00	1,434.00	1,200.00	-	
Puromycin-CPG	20-4140-13	1 x 10 µmol	472.00	519.00	434.00	-	
Puromycin-CPG	20-4140-14	1 x 15 µmol	708.00	779.00	651.00	-	
Puromycin-CPG	20-4140-41	4 x 1 µmol	262.00	288.00	241.00	-	
Puromycin-CPG	20-4140-42	4 x 0.2 µmol	157.00	173.00	144.00	-	
Pyrene-dU	10-1590-02	0.25 g	740.00	814.00	681.00	-	
Pyrene-dU	10-1590-90	100 µmol	308.00	339.00	283.00	-	
Pyrene-dU	10-1590-95	50 µmol	167.00	184.00	154.00	-	
Pyrene-dU-CE Phosphoramidite, Expedite	10-1590-90E	100 µmol	308.00	339.00	283.00	-	
Pyridin-2-one-CE Phosphoramidite	10-3012-02	0.25 g	1,783.00	1,961.00	1,640.00	-	
Pyridin-2-one-CE Phosphoramidite	10-3012-90	100 µmol	614.00	675.00	565.00	-	
Pyridin-2-one-CE Phosphoramidite	10-3012-95	50 µmol	307.00	338.00	282.00	-	
Pyrrolidine CE Phosphoramidite	10-1915-02	0.25 g	1,422.00	1,564.00	1,308.00	-	
Pyrrolidine CE Phosphoramidite	10-1915-90	100 µmol	498.00	548.00	458.00	-	
Pyrrolidine CE Phosphoramidite	10-1915-95	50 µmol	249.00	274.00	229.00	-	
Pyrrolo-CTP (10mM)	81-3017-01	0.1 mL	386.00	425.00	355.00	-	
Pyrrolo-dC-CE Phosphoramidite	10-1017-02	0.25 g	885.00	974.00	814.00	-	
Pyrrolo-dC-CE Phosphoramidite	10-1017-90	100 µmol	288.00	317.00	265.00	-	
Pyrrolo-dC-CE Phosphoramidite	10-1017-95	50 µmol	144.00	158.00	132.00	-	
Pyrrolo-dCTP (10mM)	81-1017-01	0.1 mL	215.00	237.00	198.00	-	
Quenching buffer for Glen-Pak RNA purification	60-4120-80	1 l	220.00	242.00	202.00	-	
Quenching buffer for Glen-Pak RNA purification	60-4120-82	250 mL	64.50	71.00	59.30	-	
Replacement Filters-Expedite	20-0021-0F	Pk / 20	16.10	17.70	14.80	-	
Replacement Frits - TWIST 10 µmol / 15 µmol	20-0040-0F	Pk / 20	24.20	26.60	22.30	-	
rSpacer CE Phosphoramidite	10-3914-02	0.25 g	709.00	780.00	652.00	-	
rSpacer CE Phosphoramidite	10-3914-90	100 µmol	258.00	284.00	237.00	-	
rSpacer CE Phosphoramidite	10-3914-95	50 µmol	129.00	142.00	119.00	-	
rSpacer TBDMS CE Phosphoramidite	10-3915-02	0.25 g	518.00	570.00	477.00	-	
S-Bz-Thiol-Modifier C6-dT	10-1538-02	0.25 g	721.00	793.00	663.00	-	
S-Bz-Thiol-Modifier C6-dT	10-1538-90	100 µmol	321.00	353.00	295.00	-	
S-Bz-Thiol-Modifier C6-dT	10-1538-95	50 µmol	170.00	187.00	156.00	-	
SIMA (HEX) Phosphoramidite	10-5905-90	100 µmol	229.00	252.00	211.00	-	
SIMA (HEX) Phosphoramidite,	10-5905-02	0.25 g	524.00	576.00	482.00	-	
SIMA (HEX)-dT Phosphoramidite	10-5945-90	100 µmol	885.00	974.00	814.00	-	
SolidChemical Phosphor Reagent II,	10-1902-02	0.25 g	262.00	288.00	241.00	-	
SolidChemicalPhosphor Reagent II,	10-1902-90	100 µmol	78.60	86.50	72.30	-	
Spacer C12 CE Phosphoramidite	10-1928-02	0.25 g	315.00	347.00	290.00	-	
Spacer C12 CE Phosphoramidite	10-1928-90	100 µmol	124.00	136.00	114.00	-	
Spacer Phosphoramidite 18	10-1918-02	0.25 g	242.00	266.00	223.00	-	
Spacer Phosphoramidite 18	10-1918-90	100 µmol	95.80	105.40	88.10	-	
Spacer Phosphoramidite 9	10-1909-02	0.25 g	315.00	347.00	290.00	-	
Spacer Phosphoramidite 9	10-1909-90	100 µmol	98.30	108.10	90.40	-	
Spacer Phosphoramidite C3	10-1913-02	0.25 g	315.00	347.00	290.00	-	
Spacer Phosphoramidite C3	10-1913-90	100 µmol	98.30	108.10	90.40	-	
Spacer Phosphoramidite C3, Expedite	10-1913-90E	100 µmol	98.30	108.10	90.40	-	
Spermine Phosphoramidite	10-1939-02	0.25 g	704.00	774.00	648.00	-	
Spermine Phosphoramidite	10-1939-90	100 µmol	354.00	389.00	326.00	-	
Spermine Phosphoramidite	10-1939-95	50 µmol	190.00	209.00	175.00	-	
Stearyl Phosphoramidite	10-1979-02	0.25 g	236.00	260.00	217.00	-	
Stearyl Phosphoramidite, Expedite	10-1979-02E	0.25 g	236.00	260.00	217.00	-	
sTP (10mM)	81-3522-02	0.025 mL	501.00	551.00	461.00	-	
Sulfurizing Reagent	40-4036-10	1 g	38.50	42.40	35.40	-	
Sulfurizing Reagent	40-4036-20	2 g	77.00	84.70	70.80	-	
Symmetric Doubler Phosphoramidite	10-1920-02	0.25 g	315.00	347.00	290.00	-	
Symmetric Doubler Phosphoramidite	10-1920-90	100 µmol	197.00	217.00	181.00	-	
Symmetric Doubler Phosphoramidite, Expedite	10-1920-90E	100 µmol	197.00	217.00	181.00	-	
TAMRA NHS Ester	50-5910-66	60 µL	315.00	347.00	290.00	-	
TAMRA-dT	10-1057-02	0.25 g	1,278.00	1,406.00	1,176.00	-	
TAMRA-dT	10-1057-90	100 µmol	649.00	714.00	597.00	-	
TAMRA-dT	10-1057-95	50 µmol	328.00	361.00	302.00	-	
tC <sup>+</sup> -CE Phosphoramidite	10-1517-02	250 mg	1,913.00	2,104.00	1,760.00	-	

Description	Reference	Scale/Quantity	Price List 2021				Ice Shipping
			EUR	CHF	GBP	USD	
tC <sup>+</sup> -CE Phosphoramidite	10-1517-90	100 µmol	642.00	706.00	591.00	-	
tC <sup>-</sup> -CE Phosphoramidite	10-1517-95	50 µmol	328.00	361.00	302.00	-	✓
tC-CE Phosphoramidite	10-1516-95	50 µmol	313.00	344.00	288.00	-	
tCnitro-CE Phosphoramidite	10-1518-95	50 µmol	347.00	382.00	319.00	-	
TEMPO Azide	50-2007-90	100 µmoles	459.00	505.00	422.00	-	
TEMPO Azide, 25 µmol	50-2007-92	Pk / 1	151.00	166.00	139.00	-	
Thiol-Modifier C6 S-S	10-1936-02	0.25 g	290.00	319.00	267.00	-	
Thiol-Modifier C6 S-S	10-1936-90	100 µmol	121.00	133.00	111.00	-	
Thiol-Modifier C6 S-S, Cyclone	10-1936-90C	100 µmol	121.00	133.00	111.00	-	
Thiol-Modifier C6 S-S, Expedite	10-1936-02E	0.25 g	290.00	319.00	267.00	-	
Thiol-Modifier C6 S-S, Mermade	10-1936-02M	0.25 g	290.00	319.00	267.00	-	
THPTA	50-1004-90	100 µmol	236.00	260.00	217.00	-	
Thymidine a-thiotriphosphate (15mM) (TTPaS)	80-1030-01	0.1 mL	143.00	157.00	132.00	-	
Thymidine Glycol CE Phosphoramidite	10-1096-02	0.25 g	1,278.00	1,406.00	1,176.00	-	
Thymidine Glycol CE Phosphoramidite	10-1096-90	100 µmol	544.00	598.00	500.00	-	
Thymidine Glycol CE Phosphoramidite	10-1096-95	50 µmol	282.00	310.00	259.00	-	
T-LA-CE Phosphoramidite	10-2030-10	1 g	197.00	217.00	181.00	-	
T-LA-CE Phosphoramidite, Expedite	10-2030-10E	1 g	197.00	217.00	181.00	-	
TMP-F-dU-CE Phosphoramidite	10-1016-02	0.25 g	649.00	714.00	597.00	-	
TMP-F-dU-CE Phosphoramidite	10-1016-90	100 µmol	256.00	282.00	236.00	-	
Trebler Phosphoramidite	10-1922-02	0.25 g	491.00	540.00	452.00	-	
Trebler Phosphoramidite	10-1922-90	100 µmol	282.00	310.00	259.00	-	
Trebler Phosphoramidite, Expedite	10-1922-90E	100 µmol	282.00	310.00	259.00	-	
Trimer Phosphoramidite Mix 2	13-1992-90	100 µmol	1,638.00	1,802.00	1,507.00	-	
Trimer Phosphoramidite Mix 2	13-1992-95	50 µmol	819.00	901.00	753.00	-	
Trimer Phosphoramidite Mix 2 - Mix of 19 Trimers	13-1992-90E	100 µmol	1,638.00	1,802.00	1,507.00	-	
UniCap Phosphoramidite	10-4410-02	0.25 g	65.50	72.10	60.30	-	
UniCap Phosphoramidite	10-4410-05	0.5 g	131.00	144.00	121.00	-	
UniCap Phosphoramidite	10-4410-10	1 g	262.00	288.00	241.00	-	
UniCap Phosphoramidite	10-4410-20	2 g	524.00	576.00	482.00	-	
UniCap Phosphoramidite, Mermade	10-4410-10M	1 g	262.00	288.00	241.00	-	
UniCap Phosphoramidite, Mermade	10-4410-20M	2 g	524.00	576.00	482.00	-	
Universal Support II	20-5010-01	0.1 g	12.10	13.30	11.10	-	
Universal Support II	20-5010-02	0.25 g	27.50	30.30	25.30	-	
Universal Support II	20-5110-13	1 x 10 µmol	88.00	96.80	81.00	-	
Universal Support II	20-5110-41	4 x 1 µmol	53.00	58.30	48.80	-	
Universal Support II	20-5110-42	4 x 0.2 µmol	35.20	38.70	32.40	-	
Universal Support II	20-5110-45	4 x 40 nmol	35.20	38.70	32.40	-	
Universal Support II	20-5110-92	96 x 0.2 µmol	385.00	424.00	354.00	-	
Universal Support II	20-5110-95	96 x 40 nmol	385.00	424.00	354.00	-	
Universal Support II	20-5210-14	1 x 15 µmol	132.00	145.00	121.00	-	
Universal Support II	20-5210-41	4 x 1 µmol	53.00	58.30	48.80	-	
Universal Support II	20-5210-42	4 x 0.2 µmol	35.20	38.70	32.40	-	
Universal Support II	20-5210-45	4 x 40 nmol	35.20	38.70	32.40	-	
Universal Support III	20-5010-10	1 g	105.00	116.00	97.00	-	
Universal Support III HybridCPGTM	28-5010-10	1.0 g	139.00	153.00	128.00	-	
Universal Support III PS	26-5110-41	4 x 1 µmol	48.40	53.20	44.50	-	
Universal Support III PS	26-5110-42	4 x 0.2 µmol	32.30	35.50	29.70	-	
Universal Support III PS	26-5110-52	10 x 200 nmol	80.60	88.70	74.20	-	
Universal Support III PS	26-5110-55	10 x 40 nmol	80.60	88.70	74.20	-	
Universal Support III PS	26-5210-41	4 x 1 µmol	48.40	53.20	44.50	-	
Universal Support III PS	27-5010-01	0.1 g	16.50	18.20	15.20	-	
Universal Support III PS	27-5010-02	0.25 g	38.50	42.40	35.40	-	
Universal Support III PS	27-5110-92	96 x 0.2 µmol	385.00	424.00	354.00	-	
Universal Support III PS	27-5110-95	96 x 40 nmol	385.00	424.00	354.00	-	
Universal Support III PS	27-5110-P1	10 x 1.3 µmol	132.00	145.00	121.00	-	
Universal Support III PS, Expedite	26-5210-42	4 x 0.2 µmol	32.30	35.50	29.70	-	
Uridine a-thiotriphosphate (0.5mM) (UTTPaS)	80-3040-01	0.1 mL	143.00	157.00	132.00	-	
Zebularine-CE Phosphoramidite	10-3011-02	0.25 g	852.00	937.00	784.00	-	
Zebularine-CE Phosphoramidite	10-3011-90	100 µmol	328.00	361.00	302.00	-	
Zebularine-CE Phosphoramidite	10-3011-95	50 µmol	164.00	180.00	151.00	-	
<b>GlenRNA Synthesis</b>							<b>EA1E1Z</b>
2'-F-A-CE Phosphoramidite, Expedite	10-3400-02E	0.25 g	98.30	108.10	90.40	-	
2'-F-U-CE Phosphoramidite, Expedite	10-3430-02E	0.25 g	65.50	72.10	60.30	-	
2'-OME-ibu-G-CE Phosphoramidite, Expedite	10-3120-02E	0.25 g	27.70	30.50	25.50	-	
2,6-Diaminopurine-TOM-CE Phosphoramidite	10-3085-02	0.25 g	1,278.00	1,406.00	1,176.00	-	
2-Aminopurine-TBDMS-CE Phosphoramidite	10-3070-02	0.25 g	1,278.00	1,406.00	1,176.00	-	
2-Aminopurine-TBDMS-CE Phosphoramidite	10-3070-90	100 µmol	557.00	613.00	512.00	-	
2-Aminopurine-TBDMS-CE Phosphoramidite	10-3070-95	50 µmol	278.00	306.00	256.00	-	
2'-F-Ac-C-CE Phosphoramidite	10-3415-02	0.25 g	65.50	72.10	60.30	-	
2'-F-Ac-C-CE Phosphoramidite	10-3415-05	0.5 g	131.00	144.00	121.00	-	
2'-F-Ac-C-CE Phosphoramidite	10-3415-90	100 µmol	241.00	265.00	222.00	-	
2'-F-Ac-C-CE Phosphoramidite, Expedite	10-3415-05E	0.5 g	131.00	144.00	121.00	-	
2'-F-Ac-C-CE Phosphoramidite, Expedite	10-3415-02E	0.25 g	65.50	72.10	60.30	-	
2'-Fluoro-Adenosine a-thiotriphosphate (10mM) (2'-F-ATTPaS)	80-1102-01	0.1 mL	143.00	157.00	132.00	-	
2'-F-U-CE Phosphoramidite	10-3430-02	0.25 g	65.50	72.10	60.30	-	
2'-F-U-CE Phosphoramidite	10-3430-05	0.5 g	131.00	144.00	121.00	-	
2'-F-U-CE Phosphoramidite	10-3430-90	100 µmol	241.00	265.00	222.00	-	
2'ome a thiophosphoramidite	10-3170-90		393.00	432.00	362.00	-	
2'-Ome-2,6-Diaminopurine-CE Phosphoramidite	10-3124-02	0.25 g	1,310.00	1,441.00	1,205.00	-	
2'-Ome-2,6-Diaminopurine-CE Phosphoramidite	10-3124-90	100 µmol	511.00	562.00	470.00	-	
2'-Ome-2,6-Diaminopurine-CE Phosphoramidite	10-3124-95	50 µmol	269.00	296.00	247.00	-	
2'-Ome-3-deaza-5-aza-C-CE Phosphoramidite	10-3116-02	0.25 g	1,396.00	1,536.00	1,284.00	-	
2'-Ome-3-deaza-5-aza-C-CE Phosphoramidite	10-3116-90	100 µmol	508.00	559.00	467.00	-	
2'-Ome-3-deaza-5-aza-C-CE Phosphoramidite	10-3116-95	50 µmol	254.00	279.00	234.00	-	
2'-Ome-5-Br-U-CE Phosphoramidite	10-3190-02	0.25 g	1,134.00	1,247.00	1,043.00	-	
2'-Ome-5-Br-U-CE Phosphoramidite	10-3190-90	100 µmol	406.00	447.00	374.00	-	
2'-Ome-5-Br-U-CE Phosphoramidite in bulk quantities > 10 g	10-3190-BULK	1 g	2,690.00	2,959.00	2,475.00	-	✓
2'-Ome-5-F-U-CE Phosphoramidite	10-3132-02	0.25 g	1,427.00	1,570.00	1,313.00	-	
2'-Ome-5-F-U-CE Phosphoramidite	10-3132-90	100 µmol	508.00	559.00	467.00	-	
2'-Ome-5-F-U-CE Phosphoramidite	10-3132-95	50 µmol	254.00	279.00	234.00	-	

Description	Reference	Scale/Quantity	Price List 2021				Ice Shipping
			EUR	CHF	GBP	USD	
2'-O-Me-5-Me-C-CE Phosphoramidite	10-3160-02	0.25 g	885.00	974.00	814.00	-	
2'-O-Me-5-Me-C-CE Phosphoramidite	10-3160-90	100 µmol	315.00	347.00	290.00	-	
2'-O-Me-5-Me-U-CE Phosphoramidite	10-3131-02	0.25 g	472.00	519.00	434.00	-	
2'-O-Me-5-Me-U-CE Phosphoramidite	10-3131-90	100 µmol	197.00	217.00	181.00	-	
2'-O-Me-5-Me-U-CE Phosphoramidite, Mermade	10-3131-02M	0.25 g	472.00	519.00	434.00	-	
2'-O-Me-Ac-C-CE Phosphoramidite	10-3115-02	0.25 g	27.70	30.50	25.50	-	
2'-O-Me-Ac-C-CE Phosphoramidite	10-3115-05	0.5 g	55.40	60.90	51.00	-	
2'-O-Me-Ac-C-CE Phosphoramidite	10-3115-10	1 g	111.00	122.00	102.00	-	
2'-O-Me-Ac-C-CE Phosphoramidite	10-3115-90	100 µmol	18.10	19.90	16.70	-	
2'-O-Me-Ac-C-CE Phosphoramidite, Expedite	10-3115-02E	0.25 g	27.70	30.50	25.50	-	
2'-O-Me-Ac-C-CE Phosphoramidite, Expedite	10-3115-05E	0.5 g	55.40	60.90	51.00	-	
2'-O-Me-Ac-C-RNA-CPG	20-3615-01	0.1 g	52.40	57.60	48.20	-	
2'-O-Me-Ac-C-RNA-CPG	20-3615-02	0.25 g	124.00	136.00	114.00	-	
2'-O-Me-Ac-C-RNA-CPG	20-3615-10	1 g	465.00	512.00	428.00	-	
2'-O-Me-Ac-C-RNA-CPG	20-3715-13	1 x 10 µmol	295.00	325.00	271.00	-	
2'-O-Me-Ac-C-RNA-CPG	20-3715-14	1 x 15 µmol	393.00	432.00	362.00	-	
2'-O-Me-Ac-C-RNA-CPG	20-3715-41	4 x 1 µmol	131.00	144.00	121.00	-	
2'-O-Me-Ac-C-RNA-CPG	20-3715-42	4 x 0.2 µmol	98.30	108.10	90.40	-	
2'-O-Me-A-CE Phosphoramidite	10-3100-02E	0.25 g	27.70	30.50	25.50	-	
2'-O-Me-A-CE Phosphoramidite	10-3100-05	0.5 g	55.40	60.90	51.00	-	
2'-O-Me-A-CE Phosphoramidite	10-3100-10	1 g	111.00	122.00	102.00	-	
2'-O-Me-A-CE Phosphoramidite	10-3100-90	100 µmol	18.10	19.90	16.70	-	
2'-O-Me-A-CE Phosphoramidite, Expedite	10-3100-02	0.25 g	27.70	30.50	25.50	-	
2'-O-Me-A-CE Phosphoramidite, Expedite	10-3100-05E	0.5 g	55.40	60.90	51.00	-	
2'-O-Me-Adenosine a-thiotriphosphate (20mM) (2'-O-Me-ATTPaS)	80-1101-01	0.1 mL	143.00	157.00	132.00	-	
2'-O-Me-A-RNA-CPG	20-3600-01	0.1 g	52.40	57.60	48.20	-	
2'-O-Me-A-RNA-CPG	20-3600-02	0.25 g	124.00	136.00	114.00	-	
2'-O-Me-A-RNA-CPG	20-3600-10	1 g	465.00	512.00	428.00	-	
2'-O-Me-A-RNA-CPG	20-3700-13	1 x 10 µmol	295.00	325.00	271.00	-	
2'-O-Me-A-RNA-CPG	20-3700-14	1 x 15 µmol	393.00	432.00	362.00	-	
2'-O-Me-A-RNA-CPG	20-3700-41	4 x 1 µmol	131.00	144.00	121.00	-	
2'-O-Me-A-RNA-CPG	20-3700-42	4 x 0.2 µmol	98.30	108.10	90.40	-	
2'-O-Me-A-RNA-CPG, Expedite	20-3700-42E	4 x 0.2 µmol	98.30	108.10	90.40	-	
2'-O-Me-A-Thiophosphoramidite	10-3170-02	0.25 g	944.00	1,038.00	868.00	-	
2'-O-Me-C-CE Phosphoramidite	10-3110-02	0.25 g	61.00	67.10	56.10	-	
2'-O-Me-C-CE Phosphoramidite	10-3110-05	0.5 g	121.00	133.00	111.00	-	
2'-O-Me-C-CE Phosphoramidite	10-3110-10	1 g	242.00	266.00	223.00	-	
2'-O-Me-C-CE Phosphoramidite	10-3110-90	100 µmol	24.20	26.60	22.30	-	
2'-O-Me-C-CE Phosphoramidite, Expedite	10-3110-02E	0.25 g	61.00	67.10	56.10	-	
2'-O-Me-C-RNA-CPG	20-3610-01	0.1 g	52.40	57.60	48.20	-	
2'-O-Me-C-RNA-CPG	20-3610-02	0.25 g	124.00	136.00	114.00	-	
2'-O-Me-C-RNA-CPG	20-3610-10	1 g	465.00	512.00	428.00	-	
2'-O-Me-C-RNA-CPG	20-3710-13	1 x 10 µmol	295.00	325.00	271.00	-	
2'-O-Me-C-RNA-CPG	20-3710-14	1 x 15 µmol	393.00	432.00	362.00	-	
2'-O-Me-C-RNA-CPG	20-3710-41	4 x 1 µmol	131.00	144.00	121.00	-	
2'-O-Me-C-RNA-CPG	20-3710-42	4 x 0.2 µmol	98.30	108.10	90.40	-	
2'-O-Me-G-CE Phosphoramidite	10-3121-02	0.25 g	45.40	49.90	41.80	-	
2'-O-Me-G-CE Phosphoramidite	10-3121-05	0.5 g	90.70	99.80	83.40	-	
2'-O-Me-G-CE Phosphoramidite	10-3121-10	1 g	181.00	199.00	167.00	-	
2'-O-Me-G-CE Phosphoramidite	10-3121-90	100 µmol	22.20	24.40	20.40	-	
2'-O-Me-G-CE Phosphoramidite, Expedite	10-3121-02E	0.25 g	45.40	49.90	41.80	-	
2'-O-Me-G-CE Phosphoramidite, Expedite	10-3121-05E	0.5 g	90.70	99.80	83.40	-	
2'-O-Me-G-RNA-CPG	20-3621-01	0.1 g	52.40	57.60	48.20	-	
2'-O-Me-G-RNA-CPG	20-3621-02	0.25 g	124.00	136.00	114.00	-	
2'-O-Me-G-RNA-CPG	20-3621-10	1 g	465.00	512.00	428.00	-	
2'-O-Me-G-RNA-CPG	20-3721-13	1 x 10 µmol	295.00	325.00	271.00	-	
2'-O-Me-G-RNA-CPG	20-3721-14	1 x 15 µmol	393.00	432.00	362.00	-	
2'-O-Me-G-RNA-CPG	20-3721-41	4 x 1 µmol	131.00	144.00	121.00	-	
2'-O-Me-G-RNA-CPG	20-3721-42	4 x 0.2 µmol	98.30	108.10	90.40	-	
2'-O-Me-G-Thiophosphoramidite	10-3172-02	0.25 g	944.00	1,038.00	868.00	-	
2'-O-Me-ibu-G-CE Phosphoramidite	10-3120-02	0.25 g	27.70	30.50	25.50	-	
2'-O-Me-I-CE Phosphoramidite	10-3140-02	0.25 g	472.00	519.00	434.00	-	
2'-O-Me-I-CE Phosphoramidite	10-3140-90	100 µmol	197.00	217.00	181.00	-	
2'-O-Me-iPr-Pac-G-CE Phosphoramidite	10-3621-02	0.25 g	81.90	90.10	75.30	-	
2'-O-Me-iPr-Pac-G-CE Phosphoramidite	10-3621-05	0.5 g	164.00	180.00	151.00	-	
2'-O-Me-iPr-Pac-G-CE Phosphoramidite	10-3621-10	1 g	328.00	361.00	302.00	-	
2'-O-Me-iPr-Pac-G-CE Phosphoramidite	10-3621-90	100 µmol	35.80	39.40	32.90	-	
2'-O-Me-Pac-A-CE Phosphoramidite	10-3601-02	0.25 g	81.90	90.10	75.30	-	
2'-O-Me-Pac-A-CE Phosphoramidite	10-3601-05	0.5 g	164.00	180.00	151.00	-	
2'-O-Me-Pac-A-CE Phosphoramidite	10-3601-90	100 µmol	35.80	39.40	32.90	-	
2'-O-Me-TMP-5-F-U-CE Phosphoramidite	10-3111-02	0.25 g	1,396.00	1,536.00	1,284.00	-	
2'-O-Me-TMP-5-F-U-CE Phosphoramidite	10-3111-90	100 µmol	508.00	559.00	467.00	-	
2'-O-Me-TMP-5-F-U-CE Phosphoramidite	10-3111-95	50 µmol	254.00	279.00	234.00	-	
2'-O-Me-U-CE Phosphoramidite	10-3130-02	0.25 g	27.70	30.50	25.50	-	
2'-O-Me-U-CE Phosphoramidite	10-3130-05	0.5 g	55.40	60.90	51.00	-	
2'-O-Me-U-CE Phosphoramidite	10-3130-10	1 g	111.00	122.00	102.00	-	
2'-O-Me-U-CE Phosphoramidite	10-3130-90	100 µmol	18.10	19.90	16.70	-	
2'-O-Me-U-CE Phosphoramidite, Expedite	10-3130-02E	0.25 g	27.70	30.50	25.50	-	
2'-O-Me-U-RNA-CPG	20-3630-02	0.25 g	124.00	136.00	114.00	-	
2'-O-Me-U-RNA-CPG	20-3630-10	1 g	465.00	512.00	428.00	-	
2'-O-Me-U-RNA-CPG	20-3730-13	1 x 10 µmol	295.00	325.00	271.00	-	
2'-O-Me-U-RNA-CPG	20-3730-14	1 x 15 µmol	393.00	432.00	362.00	-	
2'-O-Me-U-RNA-CPG	20-3730-41	4 x 1 µmol	131.00	144.00	121.00	-	
2'-O-Me-U-RNA-CPG	20-3730-42	4 x 0.2 µmol	98.30	108.10	90.40	-	
2'-O-Me-U-Thiophosphoramidite	10-3173-02	0.25 g	944.00	1,038.00	868.00	-	
2'-Se-Me-U-CE Phosphoramidite	10-3730-02	0.25 g	1,396.00	1,536.00	1,284.00	-	
2'-Se-Me-U-CE Phosphoramidite	10-3730-90	100 µmol	501.00	551.00	461.00	-	
2'-Se-Me-U-CE Phosphoramidite	10-3730-95	50 µmol	251.00	276.00	231.00	-	
4-Thio-U-TOM-CE Phosphoramidite	10-3052-02	0.25 g	1,278.00	1,406.00	1,176.00	-	
4-Thio-U-TOM-CE Phosphoramidite	10-3052-90	100 µmol	557.00	613.00	512.00	-	
4-Thio-U-TOM-CE Phosphoramidite	10-3052-95	50 µmol	278.00	306.00	256.00	-	

Description	Reference	Scale/Quantity	Price List 2021				Ice Shipping
			EUR	CHF	GBP	USD	
5-Br-U-CE Phosphoramidite	10-3090-02	0.25 g	622.00	684.00	572.00	-	
5-Br-U-CE Phosphoramidite	10-3090-90	100 µmol	256.00	282.00	236.00	-	
5-I-U-CE Phosphoramidite	10-3091-02	0.25 g	622.00	684.00	572.00	-	
5-I-U-CE Phosphoramidite	10-3091-90	100 µmol	256.00	282.00	236.00	-	
5-Me-C-TOM-CE Phosphoramidite	10-3064-02	0.25 g	622.00	684.00	572.00	-	
5-Me-C-TOM-CE Phosphoramidite	10-3064-90	100 µmol	249.00	274.00	229.00	-	
5-Me-C-TOM-CE Phosphoramidite	10-3064-95	50 µmol	124.00	136.00	114.00	-	
5-Me-U-CE Phosphoramidite	10-3050-02	0.25 g	622.00	684.00	572.00	-	
5-Me-U-CE Phosphoramidite	10-3050-90	100 µmol	249.00	274.00	229.00	-	
Ac-A-RNA-CPG	20-3304-01	0.1 g	52.40	57.60	48.20	-	
Ac-A-RNA-CPG	20-3304-10	1 g	465.00	512.00	428.00	-	
Ac-C-CE Phosphoramidite	10-3015-02	0.25 g	32.30	35.50	29.70	-	
Ac-C-CE Phosphoramidite	10-3015-05	0.5 g	48.40	53.20	44.50	-	
Ac-C-CE Phosphoramidite	10-3015-10	1 g	90.70	99.80	83.40	-	
Ac-C-CE Phosphoramidite MerMade	10-3015-10M	1 g	90.70	99.80	83.40	-	
Ac-C-CE Phosphoramidite, Expedite	10-3015-05E	0.5 g	48.40	53.20	44.50	-	
Ac-C-RNA 500 Å	20-3415-41A	4x1 µmol	131.00	144.00	121.00	-	
Ac-C-RNA-CPG	20-3315-01	0.1 g	52.40	57.60	48.20	-	
Ac-C-RNA-CPG	20-3315-02	0.25 g	124.00	136.00	114.00	-	
Ac-C-RNA-CPG	20-3315-10	1 g	465.00	512.00	428.00	-	
Ac-C-RNA-CPG	20-3415-13	1 x 10 µmol	295.00	325.00	271.00	-	
Ac-C-RNA-CPG	20-3415-14	1 x 15 µmol	393.00	432.00	362.00	-	
Ac-C-RNA-CPG	20-3415-41	4 x 1 µmol	131.00	144.00	121.00	-	
Ac-C-RNA-CPG	20-3415-42	4 x 0.2 µmol	98.30	108.10	90.40	-	
Ac-G-CE Phosphoramidite MerMade	10-3025-10M	1 g	90.70	99.80	83.40	-	
Ac-G-RNA-CPG	20-3324-02	0.25 g	124.00	136.00	114.00	-	
Ac-G-RNA-CPG	20-3324-10	1 g	465.00	512.00	428.00	-	
A-RNA-TOM-CPG	20-3304-02	0.25 g	124.00	136.00	114.00	-	
A-RNA-TOM-CPG	20-3404-13	1 x 10 µmol	295.00	325.00	271.00	-	
A-RNA-TOM-CPG	20-3404-14	1 x 15 µmol	393.00	432.00	362.00	-	
A-RNA-TOM-CPG	20-3404-41	4 x 1 µmol	131.00	144.00	121.00	-	
A-RNA-TOM-CPG	20-3404-42	4 x 0.2 µmol	98.30	108.10	90.40	-	
A-TOM-CE Phosphoramidite	10-3004-02	0.25 g	75.60	83.20	69.60	-	
A-TOM-CE Phosphoramidite	10-3004-05	0.5 g	151.00	166.00	139.00	-	
A-TOM-CE Phosphoramidite	10-3004-10	1 g	277.00	305.00	255.00	-	
A-TOM-CE Phosphoramidite MerMade	10-3004-10M	1 g	277.00	305.00	255.00	-	
A-TOM-CE Phosphoramidite, Expedite	10-3004-10E	1 g	277.00	305.00	255.00	-	
Bz-A-CE Phosphoramidite	10-3003-02	0.25 g	32.30	35.50	29.70	-	
Bz-A-CE Phosphoramidite	10-3003-05	0.5 g	48.40	53.20	44.50	-	
Bz-A-CE Phosphoramidite	10-3003-10	1 g	90.70	99.80	83.40	-	
Bz-A-CE Phosphoramidite MerMade	10-3003-10M	1 g	90.70	99.80	83.40	-	
Bz-A-RNA-CPG	20-3303-01	0.1 g	52.40	57.60	48.20	-	
Bz-A-RNA-CPG	20-3303-02	0.25 g	124.00	136.00	114.00	-	
Bz-A-RNA-CPG	20-3303-10	1 g	465.00	512.00	428.00	-	
Bz-A-RNA-CPG	20-3403-13	1 x 10 µmol	295.00	325.00	271.00	-	
Bz-A-RNA-CPG	20-3403-14	1 x 15 µmol	393.00	432.00	362.00	-	
Bz-A-RNA-CPG	20-3403-41	4 x 1 µmol	131.00	144.00	121.00	-	
Bz-A-RNA-CPG	20-3403-42	4 x 0.2 µmol	98.30	108.10	90.40	-	
C-TOM-CE Phosphoramidite	10-3014-02	0.25 g	75.60	83.20	69.60	-	
C-TOM-CE Phosphoramidite	10-3014-05	0.5 g	151.00	166.00	139.00	-	
C-TOM-CE Phosphoramidite	10-3014-10	1 g	277.00	305.00	255.00	-	
C-TOM-CE Phosphoramidite MerMade	10-3014-10M	1 g	277.00	305.00	255.00	-	
C-TOM-CE Phosphoramidite, Expedite	10-3014-10E	1 g	277.00	305.00	255.00	-	
dmf-G-RNA-CPG	20-3329-01	0.1 g	57.00	62.70	52.40	-	
dmf-G-RNA-CPG	20-3329-02	0.25 g	136.00	150.00	125.00	-	
dmf-G-RNA-CPG	20-3429-13	1 x 10 µmol	322.00	354.00	296.00	-	
dmf-G-RNA-CPG	20-3429-14	1 x 15 µmol	429.00	472.00	395.00	-	
dmf-G-RNA-CPG	20-3429-41	4 x 1 µmol	143.00	157.00	132.00	-	
dmf-G-RNA-CPG	20-3429-42	4 x 0.2 µmol	107.00	118.00	98.00	-	
G-RNA-TOM-CPG	20-3324-01	0.1 g	52.40	57.60	48.20	-	
G-RNA-TOM-CPG	20-3424-13	1 x 10 µmol	295.00	325.00	271.00	-	
G-RNA-TOM-CPG	20-3424-14	1 x 15 µmol	393.00	432.00	362.00	-	
G-RNA-TOM-CPG	20-3424-41	4 x 1 µmol	131.00	144.00	121.00	-	
G-RNA-TOM-CPG	20-3424-42	4 x 0.2 µmol	98.30	108.10	90.40	-	
G-TOM-CE Phosphoramidite	10-3024-02	0.25 g	75.60	83.20	69.60	-	
G-TOM-CE Phosphoramidite	10-3024-05	0.5 g	151.00	166.00	139.00	-	
G-TOM-CE Phosphoramidite	10-3024-10	1 g	277.00	305.00	255.00	-	
G-TOM-CE Phosphoramidite MerMade	10-3024-10M	1 g	277.00	305.00	255.00	-	
G-TOM-CE Phosphoramidite, Expedite	10-3024-10E	1 g	277.00	305.00	255.00	-	
iPr-Pac-G-CE Phosphoramidite	10-3021-02	0.25 g	105.00	116.00	97.00	-	
iPr-Pac-G-CE Phosphoramidite	10-3021-05	0.5 g	210.00	231.00	193.00	-	
iPr-Pac-G-CE Phosphoramidite	10-3021-10	1 g	419.00	461.00	385.00	-	
iPr-Pac-G-CE Phosphoramidite, Expedite	10-3021-05E	0.5 g	210.00	231.00	193.00	-	
iPr-Pac-G-RNA 500 Å	20-3421-41A	4x1 µmol	131.00	144.00	121.00	-	
iPr-Pac-G-RNA-CPG	20-3321-01	0.1 g	52.40	57.60	48.20	-	
iPr-Pac-G-RNA-CPG	20-3321-02	0.25 g	124.00	136.00	114.00	-	
iPr-Pac-G-RNA-CPG	20-3421-13	1 x 10 µmol	295.00	325.00	271.00	-	
iPr-Pac-G-RNA-CPG	20-3421-14	1 x 15 µmol	393.00	432.00	362.00	-	
iPr-Pac-G-RNA-CPG	20-3421-41	4 x 1 µmol	131.00	144.00	121.00	-	
iPr-Pac-G-RNA-CPG	20-3421-42	4 x 0.2 µmol	98.30	108.10	90.40	-	
N6-Methyl-A-CE Phosphoramidite	10-3005-02	0.25 g	1,697.00	1,867.00	1,561.00	-	
Pac-A-CE Phosphoramidite	10-3000-02	0.25 g	98.30	108.10	90.40	-	
Pac-A-CE Phosphoramidite	10-3000-05	0.5 g	197.00	217.00	181.00	-	
Pac-A-CE Phosphoramidite	10-3000-10	1 g	360.00	396.00	331.00	-	
Pac-A-CE Phosphoramidite, Expedite	10-3000-05E	0.5 g	197.00	217.00	181.00	-	
Pac-A-RNA 500 Å	20-3400-41A	4x1 µmol	131.00	144.00	121.00	-	
Pac-A-RNA-CPG	20-3300-01	0.1 g	52.40	57.60	48.20	-	
Pac-A-RNA-CPG	20-3300-02	0.25 g	124.00	136.00	114.00	-	
Pac-A-RNA-CPG	20-3400-13	1 x 10 µmol	295.00	325.00	271.00	-	
Pac-A-RNA-CPG	20-3400-14	1 x 15 µmol	393.00	432.00	362.00	-	
Pac-A-RNA-CPG	20-3400-41	4 x 1 µmol	131.00	144.00	121.00	-	

Description	Reference	Scale/Quantity	Price List 2021				Ice Shipping
			EUR	CHF	GBP	USD	
Pac-A-RNA-CPG	20-3400-42	4 x 0.2 µmol	98.30	108.10	90.40	-	
Pseudo-Uridine-a-thiotriphosphate (1.0mM)	80-3341-01	0.1 mL	143.00	157.00	132.00	-	
Pyrrolo-C-TOM-CE Phosphoramidite	10-3017-02	0.25 g	1,605.00	1,766.00	1,477.00	-	
Pyrrolo-C-TOM-CE Phosphoramidite	10-3017-90	100 µmol	649.00	714.00	597.00	-	
Pyrrolo-C-TOM-CE Phosphoramidite	10-3017-95	50 µmol	337.00	371.00	310.00	-	
U-CE Phosphoramidite	10-3030-02	0.25 g	32.30	35.50	29.70	-	
U-CE Phosphoramidite	10-3030-05	0.5 g	48.40	53.20	44.50	-	
U-CE Phosphoramidite	10-3030-10	1 g	90.70	99.80	83.40	-	
U-CE Phosphoramidite MerMade	10-3030-10M	1 g	90.70	99.80	83.40	-	
U-CE Phosphoramidite, Expedite	10-3030-05E	0.5 g	48.40	53.20	44.50	-	
U-RNA-CPG	20-3330-01	0.1 g	52.40	57.60	48.20	-	
U-RNA-CPG	20-3330-02	0.25 g	124.00	136.00	114.00	-	
U-RNA-CPG	20-3330-10	1 g	465.00	512.00	428.00	-	
U-RNA-CPG	20-3430-13	1 x 10 µmol	295.00	325.00	271.00	-	
U-RNA-CPG	20-3430-14	1 x 15 µmol	393.00	432.00	362.00	-	
U-RNA-CPG	20-3430-41	4 x 1 µmol	131.00	144.00	121.00	-	
U-RNA-CPG	20-3430-42	4 x 0.2 µmol	98.30	108.10	90.40	-	
U-TOM-CE Phosphoramidite	10-3034-02	0.25 g	75.60	83.20	69.60	-	
U-TOM-CE Phosphoramidite	10-3034-05	0.5 g	151.00	166.00	139.00	-	
U-TOM-CE Phosphoramidite	10-3034-10	1 g	277.00	305.00	255.00	-	
U-TOM-CE Phosphoramidite MerMade	10-3034-10M	1 g	277.00	305.00	255.00	-	

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