

## EUROGENTEC ANTIBODIES PRICE LIST

Description	Reference	Qty	EUR	CHF	GBP
5-Methylcytidine • purified	BI-MECY-0100	0.1 mg	450.00	702.00	321.00
5-Methylcytidine • purified	BI-MECY-0500	0.5 mg	1575.00	2457.00	1125.00
5-Methylcytidine • purified	BI-MECY-1000	1.0 mg	2205.00	3440.00	1575.00
Actin (alpha smooth muscle)	MB-0100	0.1 mg	338.00	528.00	231.00
ALDP • 1AL-2B4 • mouse • IgG1 K	IGALD-2B4	0.1 mg	300.00	450.00	201.00
ALDP • 2AL-1D6 • mouse • IgG1 K	IGALD-1D6	0.1 mg	300.00	450.00	201.00
Alpha-Interneixin /NF66	MB-0101	0.1 mg	378.00	591.00	258.00
BrdU	MB-0200	0.1 mg	284.00	444.00	194.00
Cadherin-1 (E-cadherin)	MB-0300	0.1 mg	294.00	460.00	201.00
Cadherin-1 (E-cadherin)	MB-0301	0.1 mg	340.00	531.00	232.00
Cadherin-1 (E-cadherin)	MB-0302	0.1 mg	352.00	550.00	240.00
Cadherin-1 (E-cadherin)	MB-0303	0.1 mg	294.00	460.00	201.00
Cadherin-1 (E-cadherin)	MB-0304	0.1 mg	294.00	460.00	201.00
Cadherin-11(OB-cadherin)	MB-0306	0.1 mg	352.00	550.00	240.00
Cadherin-6 (K-cadherin)	MB-0305	0.1 mg	352.00	550.00	240.00
Cardiotin	MB-0307	0.1 mg	294.00	460.00	201.00
Cardiotin	MB-0308	0.1 mg	294.00	460.00	201.00
Cardiotin	MB-0309	0.1 mg	294.00	460.00	201.00
Cardiotin	MB-0310	0.1 mg	294.00	460.00	201.00
Catenin (alphaT-)	MB-0311	0.1 mg	312.00	488.00	213.00
Cre-Recombinase • 2cre 2D8 • mouse • IgG1 K	IGCRE-2D8	0.1 mg	300.00	450.00	201.00
Cytokeratin (basal cell)	MB-0312	0.1 mg	244.00	382.00	167.00
Cytokeratin 10 (DE-K10)	MB-0320	0.1 mg	250.00	391.00	171.00
Cytokeratin 10 (RKSE60)	MB-0319	0.1 mg	244.00	382.00	167.00
Cytokeratin 10 and 13	MB-0321	0.1 mg	250.00	391.00	171.00
Cytokeratin 13	MB-0322	0.1 mg	300.00	469.00	205.00
Cytokeratin 14	MB-0323	0.1 mg	244.00	382.00	167.00
Cytokeratin 14	MB-0324	0.1 mg	248.00	388.00	170.00
Cytokeratin 17	MB-0325	0.1 mg	300.00	469.00	205.00
Cytokeratin 18	MB-0326	0.1 mg	244.00	382.00	167.00
Cytokeratin 18	MB-0327	0.1 mg	244.00	382.00	167.00
Cytokeratin 18	MB-0328	0.1 mg	250.00	391.00	171.00
Cytokeratin 19	MB-0329	0.1 mg	244.00	382.00	167.00
Cytokeratin 4	MB-0313	0.1 mg	270.00	422.00	185.00
Cytokeratin 5 and 8	MB-0314	0.1 mg	244.00	382.00	167.00
Cytokeratin 7	MB-0315	0.1 mg	244.00	382.00	167.00
Cytokeratin 7	MB-0316	0.1 mg	294.00	460.00	201.00
Cytokeratin 8	MB-0317	0.1 mg	304.00	475.00	208.00
Cytokeratin 8	MB-0318	0.1 mg	244.00	382.00	167.00
Desmin	MB-0400	0.1 mg	286.00	447.00	195.00
Desmin	MB-0401	0.1 mg	316.00	494.00	216.00
EAPII	MB-0500	0.1 mg	378.00	591.00	258.00
EGFP • 1GFP-2A3 • mouse • IgG1 K	IGGFP-2A3	0.1 mg	300.00	450.00	201.00
EGFP • 1GFP-2A5 • mouse • IgG1 K	IGGFP-2A5	0.1 mg	300.00	450.00	201.00
Entactin	MB-0501	0.1 mg	332.00	519.00	227.00
Ep-CAM	MB-0502	0.1 mg	338.00	528.00	231.00
Episialin	MB-0503	0.1 mg	332.00	519.00	227.00
E-selectin	MB-0504	0.1 mg	344.00	538.00	235.00
E-selectin	MB-0505	0.1 mg	344.00	538.00	235.00
E-selectin	MB-0506	0.1 mg	344.00	538.00	235.00
E-selectin	MB-0507	0.1 mg	344.00	538.00	235.00
Estrogen receptor a (b region) • B10 • mouse • IgG2a K	IGER-B10	0.1 mg	300.00	450.00	201.00
Estrogen receptor a (F region) • F3 • mouse • IgG1 K	IGER-F3	0.1 mg	300.00	450.00	201.00
Ets1 (C-terminal region) • 3ET 1C10 • mouse	IGETS-1C10	0.1 mg	300.00	450.00	201.00
Ets1 (DBD) • 1ET 7F2 • mouse	IGETS-7F2	0.1 mg	300.00	450.00	201.00
FBG3	MB-0600	0.1 mg	378.00	591.00	258.00
FBG4	MB-0601	0.1 mg	378.00	591.00	258.00
Fibronectin	MB-0602	0.1 mg	332.00	519.00	227.00
Filamentous phage coat protein 98p	MB-0603	0.1 mg	276.00	432.00	189.00
Filamentous phage coat protein 98p	MB-0604	0.1 mg	276.00	432.00	189.00
Filamentous phage coat protein 98p	MB-0605	0.1 mg	276.00	432.00	189.00
FMRP • 1C3-1a • mouse • IgG1 K	IGFMR	0.1 mg	300.00	450.00	201.00
Frataxin • 2FRA-1G2 • mouse • IgG1 K	IGFRA-1G2	0.1 mg	300.00	450.00	201.00
Gal 4 • 2GV-10 • mouse • IgG1 K	IGGV-10	0.1 mg	300.00	450.00	201.00
Gal 4 • 2GV-3 • mouse • IgG1 K	IGGV-3	0.1 mg	300.00	450.00	201.00
Gal 4 • 3GV-2 • mouse • IgG1 K	IGGV-2	0.1 mg	300.00	450.00	201.00
GFAP	MB-0700	0.1 mg	294.00	460.00	201.00
GFAP	MB-0701	0.1 mg	326.00	510.00	223.00
GSK-3 / Shaggy Protein Kinase Family • 4G-1E • mouse • IgG1 K	IGGSK-4G	0.1 mg	300.00	450.00	201.00
GSK-3 / Shaggy Protein Kinase Family • 7G-2G5 • mouse • IgG1 K	IGGSK-7G.2	0.1 mg	300.00	450.00	201.00
GST • 1D10 • mouse • IgG1 K	IGGST	0.1 mg	300.00	450.00	201.00
Heparan Sulphate Proteoglycan	MB-0801	0.1 mg	332.00	519.00	227.00
Heterochromatin protein-1 a • 2HP-2G9 • mouse • IgG1 K	IGHP-2G9	0.1 mg	300.00	450.00	201.00

## EUROGENTEC ANTIBODIES PRICE LIST

Description	Reference	Qty	EUR	CHF	GBP
Heterochromatin protein-1 b • 1MOD 1A9 • mouse • IgG1 K	IGHP-1A9	0.1 mg	300.00	450.00	201.00
Heterochromatin protein-1 g • 2MOD 1G6 • mouse • IgG1 K	IGHP-1G6	0.1 mg	300.00	450.00	201.00
His-Tag • 1DM1H7 • mouse • IgG1 K	IGHIS-1H7	0.1 mg	300.00	450.00	201.00
His-Tag • 2HIS-2H5 • mouse • IgG1 K	IGHIS-2H5	0.1 mg	300.00	450.00	201.00
Human (TAF10) • 23TA-1H8 • mouse	IGTAF10-1H8	0.1 mg	300.00	450.00	201.00
Human (TAF10) • 6TA-2B11 • mouse	IGTAF10-2B11	0.1 mg	300.00	450.00	201.00
Human (TAF10) • 6TA-2F4 • mouse	IGTAF10-2F4	0.1 mg	300.00	450.00	201.00
Human (TAF5) • 12TA-2D2 • mouse	IGTAF-2D2	0.1 mg	300.00	450.00	201.00
Human (TAF5) • 1TA-1C2 • mouse	IGTAF5-1C2	0.1 mg	300.00	450.00	201.00
Human (TAF6) • 25TA-2G7 • mouse	IGTAF6-2G7	0.1 mg	300.00	450.00	201.00
Human (TAFII68) • 7TA-2B11 • mouse	IGTAF-2B11	0.1 mg	300.00	450.00	201.00
Human (TAFII68) • 8TA-2B10 • mouse	IGTAF-2B10	0.1 mg	300.00	450.00	201.00
Human (TAFII68) • 9TA-1B11 • mouse	IGTAF-1B11	0.1 mg	300.00	450.00	201.00
Human 14-3-3 / Human Stratifin (WB), Pab	CC-0001-0100	0.1 mg	340.00	503.00	232.00
Human A-FABP, C-term (WB, IF), Pab	CC-0002-0100	0.1 mg	309.00	457.00	211.00
Human A-FABP, N-term (WB), Pab	CC-0003-0100	0.1 mg	309.00	457.00	211.00
Human ataxin-3 • 2SCA-1H9 • mouse • IgG1 K	IGSCA-1H9	0.1 mg	300.00	450.00	201.00
Human EF-1 gamma (200-300 region) (WB), Pab	CC-0004-0100	0.1 mg	361.00	534.00	246.00
Human EF-1 gamma (330-430 region) (WB), Pab	CC-0005-0100	0.1 mg	361.00	534.00	246.00
Human FUS /TLS • 1FU-1D2 • mouse	IGFUS-1D2	0.1 mg	300.00	450.00	201.00
Human HSP27 (WB), Pab	CC-0006-0100	0.1 mg	340.00	503.00	232.00
Human Keratin 15 (WB, IF), Pab	CC-0007-0100	0.1 mg	340.00	503.00	232.00
Human MRP14 (WB), Pab	CC-0008-0100	0.1 mg	309.00	457.00	211.00
Human Nucleophosmin (WB), Pab	CC-0009-0100	0.1 mg	309.00	457.00	211.00
Human p28 alpha (WB), Pab	CC-0010-0100	0.1 mg	361.00	534.00	246.00
Human PAI-2 (WB), Pab	CC-0012-0100	0.1 mg	309.00	457.00	211.00
Human PCNA (WB, IF), Pab	CC-0013-0100	0.1 mg	340.00	503.00	232.00
Human Psoriasin, C-term (WB), Pab	CC-0014-0100	0.1 mg	232.00	343.00	159.00
Human Psoriasin, N-term (WB, IF), Pab	CC-0015-0100	0.1 mg	232.00	343.00	159.00
Human Psoriasis associated -FABP (WB), Pab	CC-0011-0100	0.1 mg	361.00	534.00	246.00
Human SCCA2 (WB), Pab	CC-0016-0100	0.1 mg	361.00	534.00	246.00
Human TCTP (WB), Pab	CC-0017-0100	0.1 mg	361.00	534.00	246.00
Human Thioredoxin (WB), Pab	CC-0018-0100	0.1 mg	309.00	457.00	211.00
Human Uroplakin 3 (WB, IF), Pab	CC-0019-0100	0.1 mg	309.00	457.00	211.00
Humanin	MB-0800	0.1 mg	378.00	591.00	258.00
Huntingtin (HD) • 1HU-4C8 • mouse • IgG1 K	IGHD-4C8	0.1 mg	300.00	450.00	201.00
Huntingtin (HD) • 2HU-2E8 • mouse • IgG1 K	IGHD-2E8	0.1 mg	300.00	450.00	201.00
Integrin a3A	MB-0900	0.1 mg	294.00	460.00	201.00
Integrin a3A	MB-0901	0.1 mg	294.00	460.00	201.00
Integrin a3B	MB-0902	0.1 mg	294.00	460.00	201.00
Integrin a3B /6B	MB-0903	0.1 mg	294.00	460.00	201.00
Integrin a6A	MB-0904	0.1 mg	294.00	460.00	201.00
Integrin a6B	MB-0905	0.1 mg	294.00	460.00	201.00
Integrin b1D	MB-0906	0.1 mg	294.00	460.00	201.00
Integrin b1D	MB-0907	0.1 mg	294.00	460.00	201.00
Ku80	MB-1000	0.1 mg	378.00	591.00	258.00
Ku80	MB-1001	0.1 mg	378.00	591.00	258.00
Lamin A	MB-1101	0.1 mg	310.00	485.00	212.00
Lamin A and C	MB-1102	0.1 mg	310.00	485.00	212.00
Lamin B1	MB-1103	0.1 mg	310.00	485.00	212.00
Lamin B2	MB-1104	0.1 mg	310.00	485.00	212.00
Lamin C	MB-1105	0.1 mg	310.00	485.00	212.00
Laminin	MB-1100	0.1 mg	332.00	519.00	227.00
MBP- Maltose Binding Protein • 1H4 • mouse • IgG1 K	IGMBP	0.1 mg	300.00	450.00	201.00
Mitotic cells	MB-1200	0.1 mg	284.00	444.00	194.00
NCAM	MB-1300	0.1 mg	312.00	488.00	213.00
NCAM	MB-1301	0.1 mg	332.00	519.00	227.00
Nestin	MB-1302	0.1 mg	332.00	519.00	227.00
Neurofilament 160 kD	MB-1304	0.1 mg	276.00	432.00	189.00
Neurofilament 160 kD	MB-1305	0.1 mg	276.00	432.00	189.00
Neurofilament 200 kD	MB-1306	0.1 mg	276.00	432.00	189.00
Neurofilament 200 kD	MB-1307	0.1 mg	276.00	432.00	189.00
Neurofilament 200 kD	MB-1308	0.1 mg	276.00	432.00	189.00
Neurofilament 70 kD	MB-1303	0.1 mg	326.00	510.00	223.00
NKX3.1	MB-1309	0.1 mg	378.00	591.00	258.00
NSP-1A and 1B	MB-1313	0.1 mg	276.00	432.00	189.00
NSP-1A and 1B	MB-1314	0.1 mg	276.00	432.00	189.00
NSP-A	MB-1310	0.1 mg	286.00	447.00	195.00
NSP-A	MB-1311	0.1 mg	286.00	447.00	195.00
NSP-A	MB-1312	0.1 mg	286.00	447.00	195.00
NSP-C RNL-4	MB-1315	0.1 mg	276.00	432.00	189.00
Ovarian cancer associated antigen	MB-1400	0.1 mg	320.00	500.00	219.00
P53	MB-1500	0.1 mg	332.00	519.00	227.00

## EUROGENTEC ANTIBODIES PRICE LIST

Description	Reference	Qty	EUR	CHF	GBP
PARK7 (DJ-1)	MB-1501	0.1 mg	378.00	591.00	258.00
Phospho (Tyr 214)-GSK-3 / Shaggy Protein Kinase Family • 5G-2F	IGGSK-5G	0.1 mg	300.00	450.00	201.00
Phospho RSK-2 • 2RS-1F1 • mouse • IgG1 K	IGRSK2-1F1	0.1 mg	300.00	450.00	201.00
Phospho-GSK-3 (ser9) / Shaggy Protein Kinase Family • 7G-1F2	IGGSK-7G.1	0.1 mg	300.00	450.00	201.00
Plakophilin-3	MB-1502	0.1 mg	340.00	531.00	232.00
Plakophilin-3	MB-1503	0.1 mg	340.00	531.00	232.00
Polyglutamine • 5TF1-1C2-17.2 • mouse • IgG1 K	IGPQ-1C2	0.1 mg	300.00	450.00	201.00
Progesterone receptor • 6A • mouse • IgG2a K	IGPR-6A	0.1 mg	300.00	450.00	201.00
Rat anti-rabbit heavy chain, FITC label	SO-290F-0500	0.5 ml	153.00	230.00	106.00
Rat anti-rabbit heavy chain, HRP label	SO-291H-0500	0.5 ml	287.00	431.00	198.00
Rattin (Humanin-like)	MB-1600	0.1 mg	378.00	591.00	258.00
RNA Polymerase B • 1PB7C2 • mouse • IgG1 K	IGPB-7C2	0.1 mg	300.00	450.00	201.00
RNA Polymerase B • 1PB7E11 • mouse • IgG1 K	IGPB-7E11	0.1 mg	300.00	450.00	201.00
RNA Polymerase II • 10PB2F9 • mouse • IgG3k	IGPB-2F9	0.1 mg	300.00	450.00	201.00
RNA Polymerase II • 2PB4H7 • mouse • IgG1 K	IGPB-4H7	0.1 mg	300.00	450.00	201.00
RNA Polymerase II • 4PB2A7 • mouse • IgG1 K	IGPB-2A7	0.1 mg	300.00	450.00	201.00
RNA Polymerase II • 5PB1D3 • mouse • IgG1 K	IGPB-1D3	0.1 mg	300.00	450.00	201.00
Smoothelin	MB-1700	0.1 mg	316.00	494.00	216.00
SNF2b • 2SNF 2E12 • mouse	IGSNF-2E12	0.1 mg	300.00	450.00	201.00
SUMO 2 /3 (WB, IF), Pab	AV-SM23-0100	0.1 mg	304.00	450.00	208.00
TBP • 3TF1-3G3 • mouse	IGTBP-3G3	0.1 mg	300.00	450.00	201.00
TBP • 5TF1-2C1 • mouse	IGTBP-2C1	0.1 mg	300.00	450.00	201.00
Tet Repressor	MB-1800	0.1 mg	378.00	591.00	258.00
Trp E-Tryptophan Synthetase • A21 • mouse • IgG1 K	IGTRP	0.1 mg	300.00	450.00	201.00
Trp-p8	MB-1801	0.1 mg	378.00	591.00	258.00
Vimentin	MB-1900	0.1 mg	316.00	494.00	216.00
Vimentin	MB-1901	0.1 mg	316.00	494.00	216.00
Vimentin	MB-1902	0.1 mg	344.00	538.00	235.00
VP16 • 2GV-4 • mouse • IgG1 K	IGVP-4	0.1 mg	300.00	450.00	201.00
VP16 • 5GV-2 • mouse • IgG1 K	IGVP-2	0.1 mg	300.00	450.00	201.00
VP16 • 5GV-7 • mouse • IgG1 K	IGVP-7	0.1 mg	300.00	450.00	201.00
WWdomain containing oxidoreductase	MB-2000	0.1 mg	378.00	591.00	258.00