



Price List

2018

Products for DNA research exclusively
distributed by Eurogentec



Description	Reference	Scale/Quantity	Price List 2018				Ice Shipping
			EUR	CHF	GBP	USD	
Glen Research products for oligonucleotide synthesis							EA1
GlenDNA							EA1A1Z
0.2M Saccharin 1-Methylimidazole (SMI) in Acetonitrile	30-3180-52	200 mL	67,00	-	-	-	
0.5M CSO in Anhydrous Acetonitrile	40-4632-62	2 L	1.880,00	-	-	-	
0.5M CSO in Anhydrous Acetonitrile	40-4632-52	200 mL	188,00	-	-	-	
2'-F-U-CE Phosphoramidite	10-3430-05	0.5 g	143,00	-	-	-	✓
3'-Phosphate CPG (High Load)	25-2900-46	4 x 2.5 µmol	172,00	198,00	-	-	
5-aza-5,6-dihydro-dC-CE Phosphoramidite	10-1511-02	0.25 g	1.603,00	1.843,00	-	-	
5-aza-5,6-dihydro-dC-CE Phosphoramidite	10-1511-90	100 µmol	515,00	592,00	-	-	
5-aza-5,6-dihydro-dC-CE Phosphoramidite	10-1511-95	50 µmol	258,00	297,00	-	-	
8-D-dG-CE Phosphoramidite	10-1520-02	0.25 g	344,00	395,00	-	-	
8-D-dG-CE Phosphoramidite	10-1520-90	100 µmol	129,00	148,00	-	-	
AB3900 Ac-dC-CPG 1000A, 1.0µmol	20-2115-61	200 pces	712,00	-	-	-	
AB3900 dA-CPG 1000A, 1.0µmol	20-2101-61	200 pces	712,00	-	-	-	
AB3900 dmf-dG-CPG 1000A, 1.0µmol	20-2129-61	200 pces	712,00	-	-	-	
AB3900 dT-CPG 1000A, 1.0µmol	20-2131-61	200 pces	712,00	-	-	-	
Ac-dC-CPG 1000	20-2115-65	200x40nm	516,00	-	-	-	
dA-CPG 1000	20-2101-65	200x40nm	516,00	-	-	-	
dA-Thiophosphoramidite	10-1700-02	0.25 g	515,00	592,00	-	-	✓
dA-Thiophosphoramidite	10-1700-90	100 µmol	215,00	247,00	-	-	✓
dC-Thiophosphoramidite	10-1710-02	0.25 g	515,00	592,00	-	-	✓
dC-Thiophosphoramidite	10-1710-90	100 µmol	215,00	247,00	-	-	✓
dDs-CE Phosphoramidite	10-1521-02	0.25 g	601,00	691,00	-	-	
dDs-CE Phosphoramidite	10-1521-90	100 µmol	208,00	239,00	-	-	
dG-High Load CPG	25-2120-46M	4x2,5µm	66,00	75,90	-	-	
dG-Thiophosphoramidite	10-1720-02	100 µmol	215,00	247,00	-	-	✓
dG-Thiophosphoramidite	10-1720-02	0.25 g	515,00	592,00	-	-	✓
dmf-dG-CPG	20-2129-65	200x40nm	516,00	-	-	-	
dPa-CE Phosphoramidite	10-1523-02	0.25 g	601,00	691,00	-	-	
dPa-CE Phosphoramidite	10-1523-90	100 µmol	186,00	214,00	-	-	
dT-CE Phosphoramidite, Expedite	10-1030-05E	0.5 g	7,80	-	-	-	
dT-CPG 1000	20-2131-65	200x40nm	516,00	-	-	-	
dT-Thiophosphoramidite	10-1730-02	0.25 g	515,00	592,00	-	-	✓
dT-Thiophosphoramidite	10-1730-90	10 µmol	215,00	247,00	-	-	✓
F-DMT-dC-CE Phosphoramidite	10-1410-02	0.25 g	200,00	230,00	-	-	
F-DMT-dC-CE Phosphoramidite	10-1410-90	100 µmol	107,00	123,00	-	-	
F-DMT-dG-CE Phosphoramidite	10-1420-02	0.25 g	200,00	230,00	-	-	
F-DMT-dG-CE Phosphoramidite	10-1420-90	100 µmol	107,00	123,00	-	-	
F-DMT-dT-CE Phosphoramidite	10-1430-02	0.25 g	200,00	230,00	-	-	
F-DMT-dT-CE Phosphoramidite	10-1430-90	100 µmol	107,00	123,00	-	-	
Glen Gel-Pak™ 0.2 Desalting Column	61-5002-05		43,10	49,50	-	-	
Glen Gel-Pak™ 1.0 Desalting Column	61-5010-05		50,00	57,50	-	-	
Glen Gel-Pak™ 2.5 Desalting Column	61-5025-05	pk/5	65,00	74,70	-	-	
Glen reagent 1-Me-A-CE Phosphoramidite	10-3501-02	0.25 g	1.396,00	1.605,00	-	-	
Glen reagent 1-Me-A-CE Phosphoramidite	10-3501-95	50 µmol	272,00	313,00	-	-	
Glen reagent 4,5-Dicyanoimidazole (DCI), Crystalline	30-3050-10	1 g	27,00	31,00	-	-	
Glen reagent 4,5-Dicyanoimidazole (DCI), Crystalline	30-3050-25	25 g	385,00	443,00	-	-	
Glen reagent 5-Ethylthio-1H-Tetrazole	30-3040-10	1 g	27,00	31,00	-	-	
Glen reagent 5-Ethylthio-1H-Tetrazole	30-3040-20	2 g	46,20	53,10	-	-	
Glen reagent 5-Ethylthio-1H-Tetrazole	30-3040-25	25 g	385,00	443,00	-	-	
Glen reagent 5-I-U-CE Phosphoramidite	10-3091-95	50 µmol	140,00	161,00	-	-	
Glen reagent 5-Me-dC Brancher Phosphoramidite	10-1018-02	0.25 g	723,00	831,00	-	-	
Glen reagent 5-Me-dC Brancher Phosphoramidite	10-1018-90	100 µmol	293,00	337,00	-	-	
Glen reagent 5-Me-U-CE Phosphoramidite	10-3050-95	50 µmol	136,00	156,00	-	-	
Glen reagent AAA Trimer Phosphoramidite	13-1000-90	100 µmol	1.002,00	1.152,00	-	-	
Glen reagent AAA Trimer Phosphoramidite	13-1000-95	50 µmol	501,00	576,00	-	-	
Glen reagent AAC Trimer Phosphoramidite	13-1001-90	100 µmol	1.002,00	1.152,00	-	-	
Glen reagent AAC Trimer Phosphoramidite	13-1001-95	50 µmol	501,00	576,00	-	-	
Glen reagent Ac-dC-CE Phosphoramidite	10-1015-02	0.25 g	4,10	4,71	-	-	
Glen reagent Ac-dC-CE Phosphoramidite	10-1015-02M	0.25 g	4,10	4,71	-	-	
Glen reagent Ac-dC-CE Phosphoramidite	10-1015-05	0.5 g	8,30	9,54	-	-	
Glen reagent Ac-dC-CE Phosphoramidite	10-1015-05M	0.5 g	8,30	9,54	-	-	
Glen reagent Ac-dC-CE Phosphoramidite	10-1015-10	1 g	16,50	19,00	-	-	
Glen reagent Ac-dC-CE Phosphoramidite	10-1015-10M	0.1 g	16,50	19,00	-	-	
Glen reagent Ac-dC-CE Phosphoramidite	10-1015-1B	1 g	16,50	19,00	-	-	
Glen reagent Ac-dC-CE Phosphoramidite	10-1015-1C	1 g	16,50	19,00	-	-	
Glen reagent Ac-dC-CE Phosphoramidite	10-1015-1P	1 g	16,50	19,00	-	-	
Glen reagent Ac-dC-CE Phosphoramidite	10-1015-1S	10 g	165,00	190,00	-	-	
Glen reagent Ac-dC-CE Phosphoramidite	10-1015-20	2 g	33,00	37,90	-	-	
Glen reagent Ac-dC-CE Phosphoramidite	10-1015-2C	2 g	33,00	37,90	-	-	
Glen reagent Ac-dC-CE Phosphoramidite	10-1015-40	4 g	66,00	75,90	-	-	
Glen reagent Ac-dC-CE Phosphoramidite	10-1015-50	5 g	83,00	95,40	-	-	
Glen reagent Ac-dC-CE Phosphoramidite	10-1015-B2	0.25 g	4,10	4,71	-	-	
Glen reagent Ac-dC-CE Phosphoramidite	10-1015-B5	0.5 g	8,30	9,54	-	-	
Glen reagent Ac-dC-CE Phosphoramidite	10-1015-C2	0.25 g	4,10	4,71	-	-	
Glen reagent Ac-dC-CE Phosphoramidite	10-1015-C5	0.5 g	8,30	9,54	-	-	
Glen reagent Ac-dC-CE Phosphoramidite	10-1015-P4	0.4 g	6,60	7,59	-	-	
Glen reagent Ac-dC-CPG 1000	20-2015-01	0.1 g	6,30	7,24	-	-	
Glen reagent Ac-dC-CPG 1000	20-2015-02	0.25 g	14,10	16,20	-	-	
Glen reagent Ac-dC-CPG 1000	20-2015-10	1 g	53,00	60,90	-	-	
Glen reagent Ac-dC-CPG 1000	20-2015-61	48 x 1 µmol	264,00	303,00	-	-	
Glen reagent Ac-dC-CPG 1000	20-2015-62	200 x 0.2 µmol	661,00	760,00	-	-	
Glen reagent Ac-dC-CPG 1000	20-2115-13	1 x 10 µmol	88,00	101,10	-	-	
Glen reagent Ac-dC-CPG 1000	20-2115-41	4 x 1 µmol	30,80	35,40	-	-	
Glen reagent Ac-dC-CPG 1000	20-2115-42	4 x 0.2 µmol	30,80	35,40	-	-	
Glen reagent Ac-dC-CPG 1000	20-2115-45	4 x 40 nmol	30,80	35,40	-	-	
Glen reagent Ac-dC-CPG 1000	20-2215-14	1 x 15 µmol	132,00	152,00	-	-	
Glen reagent Ac-dC-CPG 1000	20-2215-41	4 x 1 µmol	22,00	25,30	-	-	
Glen reagent Ac-dC-CPG 1000	20-2215-42	4 x 0.2 µmol	15,40	17,70	-	-	
Glen reagent Ac-dC-CPG 1000	20-2215-45	4 x 40 nmol	15,40	17,70	-	-	
Glen reagent Ac-dC-CPG 1000	20-2515-P1	10 x 1.3 µmol	106,00	122,00	-	-	
Glen reagent Ac-dC-CPG 1000	20-2515-P2	10 x 0.2 µmol	106,00	122,00	-	-	

Description	Reference	Scale/Quantity	Price List 2018				Ice Shipping
			EUR	CHF	GBP	USD	
Glen reagent Ac-dC-CPG 500	20-2013-01	0.1 g	6,30	7,24	-	-	
Glen reagent Ac-dC-CPG 500	20-2013-02	0.25 g	14,10	16,20	-	-	
Glen reagent Ac-dC-CPG 500	20-2113-13	1 x 10 µmol	88,00	101,10	-	-	
Glen reagent Ac-dC-CPG 500	20-2113-41	4 x 1 µmol	30,80	35,40	-	-	
Glen reagent Ac-dC-CPG 500	20-2113-42	4 x 0.2 µmol	30,80	35,40	-	-	
Glen reagent Ac-dC-CPG 500	20-2213-14	1 x 15 µmol	132,00	152,00	-	-	
Glen reagent Ac-dC-CPG 500	20-2213-41	4 x 1 µmol	22,00	25,30	-	-	
Glen reagent Ac-dC-CPG 500	20-2213-42	4 x 0.2 µmol	15,40	17,70	-	-	
Glen reagent Ac-dC-PACE Phosphoramidite	10-1150-02	0.25 g	143,00	164,00	-	-	
Glen reagent Ac-dC-PACE Phosphoramidite	10-1150-05	0.5 g	286,00	329,00	-	-	
Glen reagent Ac-dC-PACE Phosphoramidite	10-1150-10	1 g	573,00	659,00	-	-	
Glen reagent Ac-G-CE Phosphoramidite	10-3025-02	0.25 g	35,20	40,50	-	-	
Glen reagent Ac-G-CE Phosphoramidite	10-3025-10	1 g	99,00	113,80	-	-	
Glen reagent ACT Trimer Phosphoramidite	13-1013-90	100 µmol	1,002,00	1,152,00	-	-	
Glen reagent ACT Trimer Phosphoramidite	13-1013-95	50 µmol	501,00	576,00	-	-	
Glen reagent ATC Trimer Phosphoramidite	13-1031-90	100 µmol	1,002,00	1,152,00	-	-	
Glen reagent ATC Trimer Phosphoramidite	13-1031-95	50 µmol	501,00	576,00	-	-	
Glen reagent ATG Trimer Phosphoramidite	13-1032-90	100 µmol	1,002,00	1,152,00	-	-	
Glen reagent ATG Trimer Phosphoramidite	13-1032-95	50 µmol	501,00	576,00	-	-	
Glen reagent Br-U-CE Phosphoramidite	10-3090-95	50 µmol	140,00	161,00	-	-	
Glen reagent CAG Trimer Phosphoramidite	13-1102-90	100 µmol	1,002,00	1,152,00	-	-	
Glen reagent CAG Trimer Phosphoramidite	13-1102-95	50 µmol	501,00	576,00	-	-	
Glen reagent CAT Trimer Phosphoramidite	13-1103-90	100 µmol	1,002,00	1,152,00	-	-	
Glen reagent CAT Trimer Phosphoramidite	13-1103-95	50 µmol	501,00	576,00	-	-	
Glen reagent CCG Trimer Phosphoramidite	13-1112-90	100 µmol	1,002,00	1,152,00	-	-	
Glen reagent CCG Trimer Phosphoramidite	13-1112-95	50 µmol	501,00	576,00	-	-	
Glen reagent CGT Trimer Phosphoramidite	13-1123-90	100 µmol	1,002,00	1,152,00	-	-	
Glen reagent CGT Trimer Phosphoramidite	13-1123-95	50 µmol	501,00	576,00	-	-	
Glen reagent CTG Trimer Phosphoramidite	13-1132-90	100 µmol	1,002,00	1,152,00	-	-	
Glen reagent CTG Trimer Phosphoramidite	13-1132-95	50 µmol	501,00	576,00	-	-	
Glen reagent dA LV-PS	26-2100-42	4 x 0.2 µmol	42,50	48,90	-	-	
Glen reagent dA LV-PS	26-2100-45	4 x 40nmol	42,50	48,90	-	-	
Glen reagent dA PS	26-2600-62	200 x 200 nmol	650,00	747,00	-	-	
Glen reagent dA PS	26-2600-65	200 x 40 nmol	650,00	747,00	-	-	
Glen reagent dA, dC, dG, dT-CPG 1000	20-2241-41	4 x 1 µmol	22,00	25,30	-	-	
Glen reagent dA, dC, dG, dT-CPG 1000	20-2241-42	4 x 0.2 µmol	15,40	17,70	-	-	
Glen reagent dA, dC, dG, dT-CPG 1000	20-2241-45	4 x 40 nmol	15,40	17,70	-	-	
Glen reagent dA,dC,dG,dT CPG 1000	20-2141-41	4 x 1 µmol	30,80	35,40	-	-	
Glen reagent dA,dC,dG,dT CPG 1000	20-2141-42	4 x 0.2 µmol	30,80	35,40	-	-	
Glen reagent dA,dC,dG,dT CPG 1000	20-2141-45	4 x 40nmol	30,80	35,40	-	-	
Glen reagent dA,dC,dG,dT- LV-PS 40nm Columns	26-2140-45	4 x 40 nmol	42,50	48,90	-	-	
Glen reagent dA,dC,dG,dT- LV-PS, Columns	26-2140-42	4 x 0.2 µmol	42,50	48,90	-	-	
Glen reagent dA,dC,dG,dT-CPG 2000	20-2142-42	4 x 0.2 µmol	30,80	35,40	-	-	
Glen reagent dA,dC,dG,dT-CPG 2000	20-2242-42	4 x 0.2 µmol	15,40	17,70	-	-	
Glen reagent dA,dC,dG,dT-CPG 500	20-2140-41	4 x 1 µmol	30,80	35,40	-	-	
Glen reagent dA,dC,dG,dT-CPG 500	20-2140-42	4 x 0.2 µmol	30,80	35,40	-	-	
Glen reagent dA,dC,dG,dT-CPG 500	20-2240-41	4 x 1 µmol	22,00	25,30	-	-	
Glen reagent dA,dC,dG,dT-CPG 500	20-2240-42	4 x 0.2 µmol	15,40	17,70	-	-	
Glen reagent dA-CE Phosphoramidite	10-1000-02	0.25 g	3,90	4,48	-	-	
Glen reagent dA-CE Phosphoramidite	10-1000-02M	0.25 g	3,90	4,48	-	-	
Glen reagent dA-CE Phosphoramidite	10-1000-05	0.5 g	7,70	8,85	-	-	
Glen reagent dA-CE Phosphoramidite	10-1000-05M	1.0 g	7,70	8,85	-	-	
Glen reagent dA-CE Phosphoramidite	10-1000-10	1.0 g	15,40	17,70	-	-	
Glen reagent dA-CE Phosphoramidite	10-1000-10M	1.0 g	15,40	17,70	-	-	
Glen reagent dA-CE Phosphoramidite	10-1000-1B	1.0 g	15,40	17,70	-	-	
Glen reagent dA-CE Phosphoramidite	10-1000-1C	1.0 g	15,40	17,70	-	-	
Glen reagent dA-CE Phosphoramidite	10-1000-1P	1.0 g	15,40	17,70	-	-	
Glen reagent dA-CE Phosphoramidite	10-1000-1S	10.0 g	154,00	177,00	-	-	
Glen reagent dA-CE Phosphoramidite	10-1000-20	2.0 g	30,80	35,40	-	-	
Glen reagent dA-CE Phosphoramidite	10-1000-2C	2.0 g	30,80	35,40	-	-	
Glen reagent dA-CE Phosphoramidite	10-1000-40	4.0 g	62,00	71,30	-	-	
Glen reagent dA-CE Phosphoramidite	10-1000-50	5.0 g	77,00	88,50	-	-	
Glen reagent dA-CE Phosphoramidite	10-1000-5S	5.0 g	77,00	88,50	-	-	
Glen reagent dA-CE Phosphoramidite	10-1000-B2	0.25 g	3,90	4,48	-	-	
Glen reagent dA-CE Phosphoramidite	10-1000-C5	0.5 g	7,70	8,85	-	-	
Glen reagent dA-CE Phosphoramidite	10-1000-P4	0.4 g	6,20	7,13	-	-	
Glen reagent dA-CPG 1000	20-2001-01	0.1 g	6,30	7,24	-	-	
Glen reagent dA-CPG 1000	20-2001-02	0.25 g	14,10	16,20	-	-	
Glen reagent dA-CPG 1000	20-2001-10	1 g	53,00	60,90	-	-	
Glen reagent dA-CPG 1000	20-2001-61	48 x 1 µmol	264,00	303,00	-	-	
Glen reagent dA-CPG 1000	20-2001-62	200 x 0.2 µmol	661,00	760,00	-	-	
Glen reagent dA-CPG 1000	20-2101-13	1 x 10 µmol	88,00	101,10	-	-	
Glen reagent dA-CPG 1000	20-2101-41	4 x 1 µmol	30,80	35,40	-	-	
Glen reagent dA-CPG 1000	20-2101-42	4 x 0.2 µmol	30,80	35,40	-	-	
Glen reagent dA-CPG 1000	20-2101-45	4 x 40 nmol	30,80	35,40	-	-	
Glen reagent dA-CPG 1000	20-2201-14	1 x 15 µmol	132,00	152,00	-	-	
Glen reagent dA-CPG 1000	20-2201-41	4 x 1 µmol	22,00	25,30	-	-	
Glen reagent dA-CPG 1000	20-2201-42	4 x 0.2 µmol	15,40	17,70	-	-	
Glen reagent dA-CPG 1000	20-2201-45	4 x 40 nmol	15,40	17,70	-	-	
Glen reagent dA-CPG 1000	20-2202-42	4 x 0.2 µmol	15,40	17,70	-	-	
Glen reagent dA-CPG 1000	20-2210-14	1 x 15 µmol	132,00	152,00	-	-	
Glen reagent dA-CPG 2000	20-2002-01	0.1 g	16,50	19,00	-	-	
Glen reagent dA-CPG 2000	20-2002-02	0.25 g	33,00	37,90	-	-	
Glen reagent dA-CPG 2000	20-2002-10	1 g	116,00	133,00	-	-	
Glen reagent dA-CPG 2000	20-2102-42	4 x 0.2 µmol	30,80	35,40	-	-	
Glen reagent dA-CPG 500	20-2000-01	0.1 g	6,30	7,24	-	-	
Glen reagent dA-CPG 500	20-2000-02	0.25 g	14,10	16,20	-	-	
Glen reagent dA-CPG 500	20-2000-10	1 g	53,00	60,90	-	-	
Glen reagent dA-CPG 500	20-2100-13	1 x 10 µmol	88,00	101,10	-	-	
Glen reagent dA-CPG 500	20-2100-41	4 x 1 µmol	30,80	35,40	-	-	
Glen reagent dA-CPG 500	20-2100-42	4 x 0.2 µmol	30,80	35,40	-	-	

Description	Reference	Scale/Quantity	Price List 2018				Ice Shipping
			EUR	CHF	GBP	USD	
Glen reagent dA-CPG 500	20-2200-14	1 x 15 µmol	132,00	152,00	-	-	
Glen reagent dA-CPG 500	20-2200-41	4 x 1 µmol	22,00	25,30	-	-	
Glen reagent dA-CPG 500	20-2200-42	4 x 0.2 µmol	15,40	17,70	-	-	
Glen reagent dA-CPG 500	20-2500-P2	10 x 0.2 µmol	106,00	122,00	-	-	
Glen reagent dA-CPG 500	20-2510-P2	10 x 0.2 µmol	106,00	122,00	-	-	
Glen reagent dA-High Load-CPG	25-2000-02	0.25 g	17,60	20,20	-	-	
Glen reagent dA-High Load-CPG	25-2000-10	1 g	63,00	72,40	-	-	
Glen reagent dA-High Load-CPG	25-2100-17	1 x 25 µmol	110,00	126,00	-	-	
Glen reagent dA-High Load-CPG	25-2100-46	4 x 2.5 µmol	66,00	75,90	-	-	
Glen reagent dA-High Load-CPG	25-2200-18	1 x 35 µmol	163,00	187,00	-	-	
Glen reagent dA-High Load-CPG	25-2200-46	4 x 2.5 µmol	66,00	75,90	-	-	
Glen reagent dA-PACE Phosphoramidite	10-1140-02	0.25 g	143,00	164,00	-	-	
Glen reagent dA-PACE Phosphoramidite	10-1140-05	0.5 g	286,00	329,00	-	-	
Glen reagent dA-PACE Phosphoramidite	10-1140-10	1 g	573,00	659,00	-	-	
Glen reagent dC LV-PS	26-2110-42	4 x 0.2 µmol	42,50	48,90	-	-	
Glen reagent dC LV-PS	26-2110-45	4 x 40nmol	42,50	48,90	-	-	
Glen reagent dC PS	26-2610-62	200 x 200 nmol	650,00	747,00	-	-	
Glen reagent dC PS	26-2610-65	200 x 40 nmol	650,00	747,00	-	-	
Glen reagent dC-CE Phosphoramidite	10-1010-02	0.25 g	3,90	4,48	-	-	
Glen reagent dC-CE Phosphoramidite	10-1010-02M	0.25 g	3,90	4,48	-	-	
Glen reagent dC-CE Phosphoramidite	10-1010-05	0.5 g	7,70	8,85	-	-	
Glen reagent dC-CE Phosphoramidite	10-1010-05M	0.5 g	7,70	8,85	-	-	
Glen reagent dC-CE Phosphoramidite	10-1010-10	1.0 g	15,40	17,70	-	-	
Glen reagent dC-CE Phosphoramidite	10-1010-10M	1.0 g	15,40	17,70	-	-	
Glen reagent dC-CE Phosphoramidite	10-1010-1B	1.0 g	15,40	17,70	-	-	
Glen reagent dC-CE Phosphoramidite	10-1010-1C	1.0 g	15,40	17,70	-	-	
Glen reagent dC-CE Phosphoramidite	10-1010-1P	1.0 g	15,40	17,70	-	-	
Glen reagent dC-CE Phosphoramidite	10-1010-1S	10.0 g	154,00	177,00	-	-	
Glen reagent dC-CE Phosphoramidite	10-1010-20	2.0 g	30,80	35,40	-	-	
Glen reagent dC-CE Phosphoramidite	10-1010-2C	2.0 g	30,80	35,40	-	-	
Glen reagent dC-CE Phosphoramidite	10-1010-40	4.0 g	62,00	71,30	-	-	
Glen reagent dC-CE Phosphoramidite	10-1010-50	5.0 g	77,00	88,50	-	-	
Glen reagent dC-CE Phosphoramidite	10-1010-5S	5.0 g	77,00	88,50	-	-	
Glen reagent dC-CE Phosphoramidite	10-1010-B2	0.25 g	3,90	4,48	-	-	
Glen reagent dC-CE Phosphoramidite	10-1010-B5	0.5 g	7,70	8,85	-	-	
Glen reagent dC-CE Phosphoramidite	10-1010-C2	0.25 g	3,90	4,48	-	-	
Glen reagent dC-CE Phosphoramidite	10-1010-C5	0.5 g	7,70	8,85	-	-	
Glen reagent dC-CE Phosphoramidite	10-1010-P4	0.4 g	6,20	7,13	-	-	
Glen reagent dC-CPG 1000	20-2011-01	0.1 g	6,30	7,24	-	-	
Glen reagent dC-CPG 1000	20-2011-02	0.25 g	14,10	16,20	-	-	
Glen reagent dC-CPG 1000	20-2011-10	1 g	53,00	60,90	-	-	
Glen reagent dC-CPG 1000	20-2011-61	48 x1 µm	225,00	259,00	-	-	
Glen reagent dC-CPG 1000	20-2111-13	1 x 10 µmol	88,00	101,10	-	-	
Glen reagent dC-CPG 1000	20-2111-41	4 x 1 µmol	30,80	35,40	-	-	
Glen reagent dC-CPG 1000	20-2111-42	4 x 0.2 µmol	30,80	35,40	-	-	
Glen reagent dC-CPG 1000	20-2111-45	4 x 40 nmol	30,80	35,40	-	-	
Glen reagent dC-CPG 1000	20-2211-14	1 x 15 µmol	132,00	152,00	-	-	
Glen reagent dC-CPG 1000	20-2211-41	4 x 1 µmol	22,00	25,30	-	-	
Glen reagent dC-CPG 1000	20-2211-42	4 x 0.2 µmol	15,40	17,70	-	-	
Glen reagent dC-CPG 1000	20-2211-45	4 x 40 nmol	15,40	17,70	-	-	
Glen reagent dC-CPG 2000	20-2012-01	0.1 g	16,50	19,00	-	-	
Glen reagent dC-CPG 2000	20-2012-02	0.25 g	33,00	37,90	-	-	
Glen reagent dC-CPG 2000	20-2012-10	1 g	116,00	133,00	-	-	
Glen reagent dC-CPG 2000	20-2112-42	4 x 0.2 µmol	30,80	35,40	-	-	
Glen reagent dC-CPG 2000	20-2212-42	4 x 0.2 µmol	15,40	17,70	-	-	
Glen reagent dC-CPG 500	20-2010-01	0.1 g	6,30	7,24	-	-	
Glen reagent dC-CPG 500	20-2010-02	0.25 g	14,10	16,20	-	-	
Glen reagent dC-CPG 500	20-2010-10	1 g	53,00	60,90	-	-	
Glen reagent dC-CPG 500	20-2110-13	1 x 10 µmol	88,00	101,10	-	-	
Glen reagent dC-CPG 500	20-2110-41	4 x 1 µmol	30,80	35,40	-	-	
Glen reagent dC-CPG 500	20-2110-42	4 x 0.2 µmol	30,80	35,40	-	-	
Glen reagent dC-CPG 500	20-2210-41	4 x 1 µmol	22,00	25,30	-	-	
Glen reagent dC-CPG 500	20-2210-42	4 x 0.2 µmol	15,40	17,70	-	-	
Glen reagent dC-High Load-CPG	25-2010-02	0.25 g	17,60	20,20	-	-	
Glen reagent dC-High Load-CPG	25-2010-10	1 g	63,00	72,40	-	-	
Glen reagent dC-High Load-CPG	25-2110-17	1 x 25 µmol	110,00	126,00	-	-	
Glen reagent dC-High Load-CPG	25-2110-46	4 x 2.5 µmol	66,00	75,90	-	-	
Glen reagent dC-High Load-CPG	25-2210-18	1 x 35 µmol	163,00	187,00	-	-	
Glen reagent dC-High Load-CPG	25-2210-46	4 x 2.5 µmol	66,00	75,90	-	-	
Glen reagent dG LV-PS	26-2120-42	4 x 0.2 µmol	42,50	48,90	-	-	
Glen reagent dG LV-PS	26-2120-45	4 x 40 nmol	42,50	48,90	-	-	
Glen reagent dG-CE Phosphoramidite	10-1020-02	0.25 g	3,90	4,48	-	-	
Glen reagent dG-CE Phosphoramidite	10-1020-02M	0.25 g	3,90	4,48	-	-	
Glen reagent dG-CE Phosphoramidite	10-1020-05	0.5 g	7,70	8,85	-	-	
Glen reagent dG-CE Phosphoramidite	10-1020-05M	0.5 g	7,70	8,85	-	-	
Glen reagent dG-CE Phosphoramidite	10-1020-10	1 g	15,40	17,70	-	-	
Glen reagent dG-CE Phosphoramidite	10-1020-10M	1 g	15,40	17,70	-	-	
Glen reagent dG-CE Phosphoramidite	10-1020-1B	1 g	15,40	17,70	-	-	
Glen reagent dG-CE Phosphoramidite	10-1020-1C	1 g	15,40	17,70	-	-	
Glen reagent dG-CE Phosphoramidite	10-1020-1P	1 g	15,40	17,70	-	-	
Glen reagent dG-CE Phosphoramidite	10-1020-1S	10 g	154,00	177,00	-	-	
Glen reagent dG-CE Phosphoramidite	10-1020-20	2 g	30,80	35,40	-	-	
Glen reagent dG-CE Phosphoramidite	10-1020-2C	2 g	30,80	35,40	-	-	
Glen reagent dG-CE Phosphoramidite	10-1020-40	4 g	62,00	71,30	-	-	
Glen reagent dG-CE Phosphoramidite	10-1020-50	5 g	77,00	88,50	-	-	
Glen reagent dG-CE Phosphoramidite	10-1020-5S	5 g	77,00	88,50	-	-	
Glen reagent dG-CE Phosphoramidite	10-1020-B2	0.25 g	3,90	4,48	-	-	
Glen reagent dG-CE Phosphoramidite	10-1020-B5	0.5 g	7,70	8,85	-	-	
Glen reagent dG-CE Phosphoramidite	10-1020-C2	0.25 g	3,90	4,48	-	-	
Glen reagent dG-CE Phosphoramidite	10-1020-C5	0.5 g	7,70	8,85	-	-	
Glen reagent dG-CE Phosphoramidite	10-1020-P4	0.4 g	6,20	7,13	-	-	
Glen reagent dG-CPG 1000	20-2021-01	0.1 g	6,30	7,24	-	-	
Glen reagent dG-CPG 1000	20-2021-02	0.25 g	14,10	16,20	-	-	

Description	Reference	Scale/Quantity	Price List 2018				Ice Shipping
			EUR	CHF	GBP	USD	
Glen reagent dG-CPG 1000	20-2021-10	1 g	53,00	60,90	-	-	
Glen reagent dG-CPG 1000	20-2021-61	48 x 1 µmol	264,00	303,00	-	-	
Glen reagent dG-CPG 1000	20-2021-62	200 x 0.2 µmol	661,00	760,00	-	-	
Glen reagent dG-CPG 1000	20-2121-13	1 x 10 µmol	88,00	101,10	-	-	
Glen reagent dG-CPG 1000	20-2121-41	4 x 1 µmol	30,80	35,40	-	-	
Glen reagent dG-CPG 1000	20-2121-42	4 x 0.2 µmol	30,80	35,40	-	-	
Glen reagent dG-CPG 1000	20-2121-45	4 x 40 nmol	30,80	35,40	-	-	
Glen reagent dG-CPG 1000	20-2221-14	1 x 15 µmol	132,00	152,00	-	-	
Glen reagent dG-CPG 1000	20-2221-41	4 x 1 µmol	22,00	25,30	-	-	
Glen reagent dG-CPG 1000	20-2221-42	4 x 0.2 µmol	15,40	17,70	-	-	
Glen reagent dG-CPG 1000	20-2221-45	4 x 40 nmol	15,40	17,70	-	-	
Glen reagent dG-CPG 2000	20-2022-01	0.1 g	16,50	19,00	-	-	
Glen reagent dG-CPG 2000	20-2022-02	0.25 g	33,00	37,90	-	-	
Glen reagent dG-CPG 2000	20-2022-10	1 g	116,00	133,00	-	-	
Glen reagent dG-CPG 2000	20-2122-42	4 x 0.2 µmol	30,80	35,40	-	-	
Glen reagent dG-CPG 2000	20-2222-42	4 x 0.2 µmol	15,40	17,70	-	-	
Glen reagent dG-CPG 500	20-2020-01	0.1 g	6,30	7,24	-	-	
Glen reagent dG-CPG 500	20-2020-02	0.25 g	14,10	16,20	-	-	
Glen reagent dG-CPG 500	20-2020-10	1 g	53,00	60,90	-	-	
Glen reagent dG-CPG 500	20-2120-13	1 x 10 µmol	88,00	101,10	-	-	
Glen reagent dG-CPG 500	20-2120-41	4 x 1 µmol	30,80	35,40	-	-	
Glen reagent dG-CPG 500	20-2120-42	4 x 0.2 µmol	30,80	35,40	-	-	
Glen reagent dG-CPG 500	20-2220-14	1 x 15 µmol	132,00	152,00	-	-	
Glen reagent dG-CPG 500	20-2220-41	4 x 1 µmol	22,00	25,30	-	-	
Glen reagent dG-CPG 500	20-2220-42	4 x 0.2 µmol	15,40	17,70	-	-	
Glen reagent dG-CPG 500	20-2520-P2	10 x 0.2 µmol	106,00	122,00	-	-	
Glen reagent dG-High Load CPG	25-2020-02	0.25 g	17,60	20,20	-	-	
Glen reagent dG-High Load CPG	25-2020-10	1 g	63,00	72,40	-	-	
Glen reagent dG-High Load CPG	25-2120-17	1 x 25 µmol	110,00	126,00	-	-	
Glen reagent dG-High Load CPG	25-2120-46	4 x 2.5 µmol	66,00	75,90	-	-	
Glen reagent dG-High Load CPG	25-2220-18	1 x 35 µmol	163,00	187,00	-	-	
Glen reagent dG-High Load CPG	25-2220-46	4 x 2.5 µmol	66,00	75,90	-	-	
Glen reagent dG-PACE Phosphoramidite	10-1160-02	0.25 g	143,00	164,00	-	-	
Glen reagent dG-PACE Phosphoramidite	10-1160-05	0.5 g	286,00	329,00	-	-	
Glen reagent dG-PACE Phosphoramidite	10-1160-10	1 g	573,00	659,00	-	-	
Glen reagent dmf-dG-CE Phosphoramidite	10-1029-02M	0.25 g	3,90	4,48	-	-	
Glen reagent dmf-dG-CE Phosphoramidite	10-1029-05M	0.5 g	7,70	8,85	-	-	
Glen reagent dmf-dG-CE Phosphoramidite	10-1029-10M	1 g	15,40	17,70	-	-	
Glen reagent dmf-dG-CE Phosphoramidite	10-1029-2C	2 g	30,80	35,40	-	-	
Glen reagent dmf-dG-CE Phosphoramidite	10-1029-40	4 g	62,00	71,30	-	-	
Glen reagent dT LV-PS	26-2130-42	4 x 0.2 µmol	42,50	48,90	-	-	
Glen reagent dT LV-PS	26-2130-45	4 x 40 nmol	42,50	48,90	-	-	
Glen reagent dT-CE Phosphoramidite	10-1030-02	0.25 g	3,90	4,48	-	-	
Glen reagent dT-CE Phosphoramidite	10-1030-02M	0.25 g	3,90	4,48	-	-	
Glen reagent dT-CE Phosphoramidite	10-1030-05	0.5 g	7,70	8,85	-	-	
Glen reagent dT-CE Phosphoramidite	10-1030-05M	0.5 g	7,70	8,85	-	-	
Glen reagent dT-CE Phosphoramidite	10-1030-10	1 g	15,40	17,70	-	-	
Glen reagent dT-CE Phosphoramidite	10-1030-10M	1 g	15,40	17,70	-	-	
Glen reagent dT-CE Phosphoramidite	10-1030-1B	1 g	15,40	17,70	-	-	
Glen reagent dT-CE Phosphoramidite	10-1030-1C	1 g	15,40	17,70	-	-	
Glen reagent dT-CE Phosphoramidite	10-1030-1P	1 g	15,40	17,70	-	-	
Glen reagent dT-CE Phosphoramidite	10-1030-1S	10 g	154,00	177,00	-	-	
Glen reagent dT-CE Phosphoramidite	10-1030-20	2 g	30,80	35,40	-	-	
Glen reagent dT-CE Phosphoramidite	10-1030-2C	2 g	30,80	35,40	-	-	
Glen reagent dT-CE Phosphoramidite	10-1030-40	4 g	62,00	71,30	-	-	
Glen reagent dT-CE Phosphoramidite	10-1030-50	5 g	77,00	88,50	-	-	
Glen reagent dT-CE Phosphoramidite	10-1030-55	5 g	77,00	88,50	-	-	
Glen reagent dT-CE Phosphoramidite	10-1030-82	0.25 g	3,90	4,48	-	-	
Glen reagent dT-CE Phosphoramidite	10-1030-B5	0.5 g	7,70	8,85	-	-	
Glen reagent dT-CE Phosphoramidite	10-1030-C2	0.25 g	3,90	4,48	-	-	
Glen reagent dT-CE Phosphoramidite	10-1030-C5	0.5 g	7,70	8,85	-	-	
Glen reagent dT-CE Phosphoramidite	10-1030-P4	0.4 g	6,20	7,13	-	-	
Glen reagent dT-CE Phosphoramidite, Expedite	10-1030-07	0.75 g	22,50	25,90	-	-	
Glen reagent dT-CPG 1000	20-2031-01	0.1 g	6,30	7,24	-	-	
Glen reagent dT-CPG 1000	20-2031-02	0.25 g	14,10	16,20	-	-	
Glen reagent dT-CPG 1000	20-2031-10	1 g	53,00	60,90	-	-	
Glen reagent dT-CPG 1000	20-2031-61	48 x 1 µmol	264,00	303,00	-	-	
Glen reagent dT-CPG 1000	20-2031-62	200 x 0.2 µmol	661,00	760,00	-	-	
Glen reagent dT-CPG 1000	20-2131-13	1 x 10 µmol	88,00	101,10	-	-	
Glen reagent dT-CPG 1000	20-2131-41	4 x 1.0 µmol	30,80	35,40	-	-	
Glen reagent dT-CPG 1000	20-2131-42	4 x 0.2 µmol	30,80	35,40	-	-	
Glen reagent dT-CPG 1000	20-2131-45	4 x 40 nmol	30,80	35,40	-	-	
Glen reagent dT-CPG 1000	20-2231-14	1 x 15 µmol	132,00	152,00	-	-	
Glen reagent dT-CPG 1000	20-2231-41	4 x 1 µmol	22,00	25,30	-	-	
Glen reagent dT-CPG 1000	20-2231-42	4 x 0.2 µmol	15,40	17,70	-	-	
Glen reagent dT-CPG 1000	20-2231-45	4 x 40 nmol	15,40	17,70	-	-	
Glen reagent dT-CPG 2000	20-2032-01	0.1 g	16,50	19,00	-	-	
Glen reagent dT-CPG 2000	20-2032-02	0.25 g	33,00	37,90	-	-	
Glen reagent dT-CPG 2000	20-2032-10	1 g	116,00	133,00	-	-	
Glen reagent dT-CPG 2000	20-2132-42	4 x 0.2 µmol	30,80	35,40	-	-	
Glen reagent dT-CPG 2000	20-2232-42	4 x 0.2 µmol	15,40	17,70	-	-	
Glen reagent dT-CPG 500	20-2030-01	0.1 g	6,30	7,24	-	-	
Glen reagent dT-CPG 500	20-2030-02	0.25 g	14,10	16,20	-	-	
Glen reagent dT-CPG 500	20-2030-10	1 g	53,00	60,90	-	-	
Glen reagent dT-CPG 500	20-2130-13	1 x 10 µmol	88,00	101,10	-	-	
Glen reagent dT-CPG 500	20-2130-41	4 x 1.0 µmol	30,80	35,40	-	-	
Glen reagent dT-CPG 500	20-2130-42	4 x 0.2 µmol	30,80	35,40	-	-	
Glen reagent dT-CPG 500	20-2230-14	1 x 15 µmol	132,00	152,00	-	-	
Glen reagent dT-CPG 500	20-2230-41	4 x 1 µmol	22,00	25,30	-	-	
Glen reagent dT-CPG 500	20-2230-42	4 x 0.2 µmol	15,40	17,70	-	-	
Glen reagent dT-CPG 500	20-2530-P2	10 x 0.2 µmol	106,00	122,00	-	-	
Glen reagent dT-High Load-CPG	25-2030-02	0.25 g	17,60	20,20	-	-	
Glen reagent dT-High Load-CPG	25-2030-10	1 g	63,00	72,40	-	-	

Description	Reference	Scale/Quantity	Price List 2018				Ice Shipping
			EUR	CHF	GBP	USD	
Glen reagent dT-High Load-CPG	25-2130-17	1 x 25 µmol	110,00	126,00	-	-	
Glen reagent dT-High Load-CPG	25-2130-46	4 x 2.5 µmol	66,00	75,90	-	-	
Glen reagent dT-High Load-CPG	25-2230-18	1 x 35 µmol	163,00	187,00	-	-	
Glen reagent dT-High Load-CPG	25-2230-46	4 x 2.5 µmol	66,00	75,90	-	-	
Glen reagent dT-Icaa-CPG 2000	20-2232-41	4 x 1 µm	31,30	36,00	-	-	
Glen reagent dT-PACE Phosphoramidite	10-1170-02	0.25 g	143,00	164,00	-	-	
Glen reagent dT-PACE Phosphoramidite	10-1170-05	0.5 g	286,00	329,00	-	-	
Glen reagent dT-PACE Phosphoramidite	10-1170-10	1 g	573,00	659,00	-	-	
Glen reagent dT-PS	26-2630-62	200 x 200 nmol	650,00	747,00	-	-	
Glen reagent dT-PS	26-2630-65	200 x 40nmol	650,00	747,00	-	-	
Glen reagent dT-Q-CPG 500	21-2230-42	4 x 0.2 µmol	35,20	40,50	-	-	
Glen reagent Empty MerMade Columns, pk/48	20-0050-05	50 nmol	176,00	202,00	-	-	
Glen reagent Empty Synthesis Columns - TWIST 10um/15um, Pk/10	20-0040-00	Pk / 10	264,00	303,00	-	-	
Glen reagent Empty Synthesis Columns, 1um Expedite Style, Pk/10	20-0021-01	Pk / 10	42,30	48,60	-	-	
Glen reagent Empty Synthesis Columns, 40nm, 0.2um Expedite Style, Pk/10	20-0021-02	Pk / 10	42,30	48,60	-	-	
Glen reagent Empty Synthesis Columns-TWIST 40nm, 0.2um or 1um, Pk/10	20-0030-00	Pk / 10	53,00	60,90	-	-	
Glen reagent F-DMT-dA-CE Phosphoramidite	10-1400-02	0.25 g	200,00	230,00	-	-	
Glen reagent F-DMT-dA-CE Phosphoramidite	10-1400-90	100 µmol	107,00	123,00	-	-	
Glen reagent GAA Trimer Phosphoramidite	13-1200-90	100 µmol	1.002,00	1.152,00	-	-	
Glen reagent GAA Trimer Phosphoramidite	13-1200-95	50 µmol	501,00	576,00	-	-	
Glen reagent GAC Trimer Phosphoramidite	13-1201-90	100 µmol	1.002,00	1.152,00	-	-	
Glen reagent GAC Trimer Phosphoramidite	13-1201-95	50 µmol	501,00	576,00	-	-	
Glen reagent GCT Trimer Phosphoramidite	13-1213-90	100 µmol	1.002,00	1.152,00	-	-	
Glen reagent GCT Trimer Phosphoramidite	13-1213-95	50 µmol	501,00	576,00	-	-	
Glen reagent GGT Trimer Phosphoramidite	13-1223-90	100 µmol	1.002,00	1.152,00	-	-	
Glen reagent GGT Trimer Phosphoramidite	13-1223-95	50 µmol	501,00	576,00	-	-	
Glen reagent GTT Trimer Phosphoramidite	13-1233-90	100 µmol	1.002,00	1.152,00	-	-	
Glen reagent GTT Trimer Phosphoramidite	13-1233-95	50 µmol	501,00	576,00	-	-	
Glen reagent I-CE Phosphoramidite	10-3040-90	100 µmol	272,00	313,00	-	-	
Glen reagent I-CE Phosphoramidite	10-3040-95	50 µmol	136,00	156,00	-	-	
Glen reagent pdU-CE Phosphoramidite	10-1054-90	100 µmol	93,00	106,90	-	-	
Glen reagent TAC Trimer Phosphoramidite	13-1301-90	100 µmol	1.002,00	1.152,00	-	-	
Glen reagent TAC Trimer Phosphoramidite	13-1301-95	50 µmol	501,00	576,00	-	-	
Glen reagent TCT Trimer Phosphoramidite	13-1313-90	100 µmol	1.002,00	1.152,00	-	-	
Glen reagent TCT Trimer Phosphoramidite	13-1313-95	50 µmol	501,00	576,00	-	-	
Glen reagent TGC Trimer Phosphoramidite	13-1321-90	100 µmol	1.002,00	1.152,00	-	-	
Glen reagent TGC Trimer Phosphoramidite	13-1321-95	50 µmol	501,00	576,00	-	-	
Glen reagent TGG Trimer Phosphoramidite	13-1322-90	100 µmol	1.002,00	1.152,00	-	-	
Glen reagent TGG Trimer Phosphoramidite	13-1322-95	50 µmol	501,00	576,00	-	-	
Glen reagent Trimer Phosphoramidite Mix 1	13-1991-90	100 µmol	1.474,00	1.694,00	-	-	
Glen reagent Trimer Phosphoramidite Mix 1	13-1991-95	50 µmol	737,00	847,00	-	-	
Glen reagent TTC Trimer Phosphoramidite	13-1331-90	100 µmol	1.002,00	1.152,00	-	-	
Glen reagent TTC Trimer Phosphoramidite	13-1331-95	50 µmol	501,00	576,00	-	-	
Glen UnySupport 1000	20-5041-01	0.1 g	12,10	13,90	-	-	
Glen UnySupport 1000	20-5041-02	0.25 g	27,50	31,60	-	-	
Glen UnySupport 1000	20-5041-10	1 g	105,00	121,00	-	-	
Glen UnySupport 500	20-5040-01	0.1 g	12,10	13,90	-	-	
Glen UnySupport 500	20-5040-02	0.25 g	27,50	31,60	-	-	
Glen UnySupport 500	20-5040-10	1 g	105,00	121,00	-	-	
Glen UnySupport 500	20-5140-13	1 x 10 µmol	88,00	101,10	-	-	
Glen UnySupport 500	20-5140-41	4 x 1 µmol	53,00	60,90	-	-	
Glen UnySupport 500	20-5140-42	4 x 0.2 µmol	35,20	40,50	-	-	
Glen UnySupport 500	20-5140-45	4 x 40 nmol	35,20	40,50	-	-	
Glen UnySupport 500	20-5140-91	96 x 1 µmol	578,00	664,00	-	-	
Glen UnySupport 500	20-5140-92	96 x 0.2 µmol	385,00	443,00	-	-	
Glen UnySupport 500	20-5140-95	96 x 40 nmol	385,00	443,00	-	-	
Glen UnySupport 500	20-5240-14	1 x 15 µmol	132,00	152,00	-	-	
Glen UnySupport 500	20-5240-42	4 x 0.2 µmol	35,20	40,50	-	-	
Glen UnySupport 500	20-5240-45	4 x 40 nmol	35,20	40,50	-	-	
Glen UnySupport PS	26-5140-51	10 x 1 µmol	132,00	152,00	-	-	
Glen UnySupport PS	26-5140-52	10 x 200 nmol	88,00	101,10	-	-	
Glen UnySupport PS	26-5140-55	10 x 40 nmol	88,00	101,10	-	-	
Glen UnySupport™ 1000	20-5241-42	4 x 0,2µm	36,00	-	-	-	
Glen UnySupport™ FC	22-5141-55	10x40nm	86,00	-	-	-	
Glen UnySupport™ FC	22-5141-95	96x40nm	376,00	-	-	-	
High Load Glen UnySupport	25-5040-10	1 g	109,00	-	-	-	EA1B1Z
Pac-ds-CE Phosphoramidite	10-1522-02	0.25 g	601,00	691,00	-	-	
Pac-ds-CE Phosphoramidite	10-1522-90	100 µmol	243,00	279,00	-	-	
Universal Support II	20-5110-91	96 x 1 µmol	578,00	664,00	-	-	
Universal Support III PS	26-5010-10	1 g	156,00	179,00	-	-	
UnyLinker™ CPG 500	20-5030-01	0.1 g	12,10	13,90	-	-	
UnyLinker™ CPG 500	20-5030-02	0.25 g	27,50	31,60	-	-	
UnyLinker™ CPG 500	20-5030-10	1 g	105,00	121,00	-	-	
UnyLinker™ CPG 500	20-5130-13	1 x 10 µmol	88,00	101,10	-	-	
UnyLinker™ CPG 500	20-5130-41	4 x 1 µmol	53,00	60,90	-	-	
UnyLinker™ CPG 500	20-5130-42	4 x 0.2 µmol	35,20	40,50	-	-	
UnyLinker™ CPG 500	20-5130-45	4 x 40 nmol	35,20	40,50	-	-	
UnyLinker™ CPG 500	20-5230-14	1 x 15 µmol	132,00	152,00	-	-	
UnyLinker™ CPG 500	20-5230-41	4 x 1 µmol	53,00	60,90	-	-	
UnyLinker™ CPG 500	20-5230-42	4 x 0.2 µmol	35,20	40,50	-	-	
UnyLinker™ CPG 500	20-5230-45	4 x 40 nmol	35,20	40,50	-	-	
Glen Solvant							
0.5M CSO in Anhydrous Acetonitrile (Expedite)	40-4632-52E	200 mL	188,00	-	-	-	
3% Dichloroacetic acid in Dichloromethane	40-4040-61	960 mL	58,00	66,70	-	-	
Cap Mix 1L (500ml old ref. 40-4026-75 + 500ml old ref. 40-4027-75) Glen Reagent Acetonitrile/Ac20, 500ml + Glen reagent Acetonitrile/Lutidine	40-4028-71	500 mL	123,00	141,00	-	-	
Conjugation Reagent (TCEP)	50-1003-01	1 g	107,00	123,00	-	-	
Fluorous Chemical Phosphorylation Reagent II	10-1904-02	0.25 g	279,00	321,00	-	-	
Fluorous Chemical Phosphorylation Reagent II	10-1904-90	100 µmol	136,00	156,00	-	-	
Glen reagent 0.02M Iodine in Tetrahydrofuran/ Water/ Pyridine	40-4132-52	200 mL	30,80	35,40	-	-	
Glen reagent 0.02M Iodine in Tetrahydrofuran/ Water/ Pyridine	40-4132-55	240 mL	37,00	42,50	-	-	
Glen reagent 0.02M Iodine in Tetrahydrofuran/ Water/ Pyridine	40-4132-57	450 mL	74,00	85,10	-	-	
Glen reagent 0.02M Iodine in Tetrahydrofuran/ Water/ Pyridine	40-4132-62	2000 mL	251,00	289,00	-	-	

Description	Reference	Scale/Quantity	Price List 2018				Ice Shipping
			EUR	CHF	GBP	USD	
Glen reagent 0.02M Iodine in Tetrahydrofuran/ Water/ Pyridine	40-4132-66	60 mL	15,40	17,70	-	-	
Glen reagent 0.02M Iodine in Tetrahydrofuran/Pyridine/Water	40-4330-52	200 mL	23,10	26,60	-	-	
Glen reagent 0.02M Iodine in Tetrahydrofuran/Pyridine/Water	40-4330-57	450 mL	56,00	64,40	-	-	
Glen reagent 0.02M Iodine in Tetrahydrofuran/Pyridine/Water	40-4330-61	960 mL	119,00	137,00	-	-	
Glen reagent 0.02M Iodine in Tetrahydrofuran/Pyridine/Water	40-4330-62	2000 mL	251,00	289,00	-	-	
Glen reagent 0.05M I2 in Pyridine/H2O	40-4035-71	1000 mL	174,00	200,00	-	-	
Glen reagent 0.05M Potassium Carbonate in Methanol	60-4600-30	30 mL	23,10	26,60	-	-	
Glen reagent 0.25M 5-Benzylthio-1H-Tetrazole in Acetonitrile	30-3172-52	200 mL	77,00	88,50	-	-	
Glen reagent 0.25M 5-Benzylthio-1H-Tetrazole in Acetonitrile, Expedite	30-3172-52E	200 mL	77,00	88,50	-	-	
Glen reagent 0.25M 5-Benzylthio-1H-tetrazole in Acetonitrile	30-3170-45	45 mL	30,80	35,40	-	-	
Glen reagent 0.25M 5-Benzylthio-1H-tetrazole in Acetonitrile	30-3170-52	200 mL	77,00	88,50	-	-	
Glen reagent 0.25M 5-Benzylthio-1H-tetrazole in Acetonitrile	30-3170-57	450 mL	154,00	177,00	-	-	
Glen reagent 0.25M 5-Benzylthio-1H-tetrazole in Acetonitrile	30-3170-62	2000 mL	586,00	674,00	-	-	
Glen reagent 0.25M 5-Ethylthio-1H-Tetrazole in Anhydrous Acetonitrile	30-3140-45	45 mL	30,80	35,40	-	-	
Glen reagent 0.25M 5-Ethylthio-1H-Tetrazole in Anhydrous Acetonitrile	30-3140-52	200 mL	77,00	88,50	-	-	
Glen reagent 0.25M 5-Ethylthio-1H-Tetrazole in Anhydrous Acetonitrile	30-3140-57	450 mL	154,00	177,00	-	-	
Glen reagent 0.25M 5-Ethylthio-1H-Tetrazole in Anhydrous Acetonitrile	30-3140-61	960 mL	281,00	323,00	-	-	
Glen reagent 0.25M 5-Ethylthio-1H-Tetrazole in Anhydrous Acetonitrile	30-3140-62	2000 mL	586,00	674,00	-	-	
Glen reagent 0.25M 5-Ethylthio-1H-Tetrazole in Anhydrous Acetonitrile	30-3142-52	200 mL	77,00	88,50	-	-	
Glen reagent 0.25M DCI in Anhydrous Acetonitrile	30-3150-45	45 mL	30,80	35,40	-	-	
Glen reagent 0.25M DCI in Anhydrous Acetonitrile	30-3150-52	200 mL	77,00	88,50	-	-	
Glen reagent 0.25M DCI in Anhydrous Acetonitrile	30-3150-57	450 mL	154,00	177,00	-	-	
Glen reagent 0.25M DCI in Anhydrous Acetonitrile	30-3150-61	960 mL	281,00	323,00	-	-	
Glen reagent 0.25M DCI in Anhydrous Acetonitrile	30-3150-62	2000 mL	586,00	674,00	-	-	
Glen reagent 0.25M DCI in Anhydrous Acetonitrile	30-3152-52	200 mL	77,00	88,50	-	-	
Glen reagent 0.40M Tetrazole in Acetonitrile	30-3105-71	1000 mL	443,00	509,00	-	-	
Glen reagent 10% 1-Methylimidazole in Tetrahydrofuran	40-4120-52	200 mL	23,10	26,60	-	-	
Glen reagent 10% 1-Methylimidazole in Tetrahydrofuran	40-4120-57	450 mL	56,00	64,40	-	-	
Glen reagent 10% 1-Methylimidazole in Tetrahydrofuran	40-4120-62	2000 mL	251,00	289,00	-	-	
Glen reagent 10% 1-Methylimidazole in Tetrahydrofuran/Pyridine	40-4122-51	150 mL	23,10	26,60	-	-	
Glen reagent 10% 1-Methylimidazole in Tetrahydrofuran/Pyridine	40-4122-52	200 mL	30,80	35,40	-	-	
Glen reagent 10% 1-Methylimidazole in Tetrahydrofuran/Pyridine	40-4122-57	450 mL	74,00	85,10	-	-	
Glen reagent 10% 1-Methylimidazole in Tetrahydrofuran/Pyridine	40-4122-62	2000 mL	251,00	289,00	-	-	
Glen reagent 10% 1-Methylimidazole in Tetrahydrofuran/Pyridine	40-4122-66	60 mL	15,40	17,70	-	-	
Glen reagent 16% 1-Methylimidazole in Tetrahydrofuran	40-4220-45	45 mL	15,40	17,70	-	-	
Glen reagent 16% 1-Methylimidazole in Tetrahydrofuran	40-4220-52	200 mL	30,80	35,40	-	-	
Glen reagent 16% 1-Methylimidazole in Tetrahydrofuran	40-4220-57	450 mL	74,00	85,10	-	-	
Glen reagent 16% 1-Methylimidazole in Tetrahydrofuran	40-4220-61	960 mL	157,00	180,00	-	-	
Glen reagent 16% 1-Methylimidazole in Tetrahydrofuran	40-4220-62	2000 mL	328,00	377,00	-	-	
Glen reagent 2% Aqueous Trifluoroacetic Acid	60-4040-57	450 mL	31,70	36,40	-	-	
Glen reagent 2.0M Triethylamine Acetate, HPLC grade, pH=7	60-4110-52	200 mL	53,00	60,90	-	-	
Glen reagent 2.0M Triethylamine Acetate, HPLC grade, pH=7	60-4110-62	2000 mL	352,00	405,00	-	-	
Glen reagent 2.0M Triethylamine Acetate, HPLC grade, pH=7	60-4110-57	450 mL	106,00	122,00	-	-	
Glen reagent 2.0M Triethylamine Acetate, HPLC grade, pH=7	60-4110-60	960 mL	176,00	202,00	-	-	
Glen reagent 2.5% Dichloroacetic acid in Dichloromethane	40-4042-57	450 mL	27,70	31,80	-	-	
Glen reagent 2.5% Dichloroacetic acid in Dichloromethane	40-4042-62	2000 mL	111,00	128,00	-	-	
Glen reagent 3% Aqueous Trifluoroacetic Acid	60-4041-57	450 mL	31,70	36,40	-	-	
Glen reagent 3% Dichloroacetic acid in Dichloromethane	40-4040-57	450 mL	27,70	31,80	-	-	
Glen reagent 3% Dichloroacetic acid in Dichloromethane	40-4040-62	2000 mL	111,00	128,00	-	-	
Glen reagent 3% Dichloroacetic acid in Dichloromethane	40-4040-71	1000 mL	62,00	71,30	-	-	
Glen reagent 3% TCA/DCM	40-4140-57	450 mL	27,70	31,80	-	-	
Glen reagent 3% TCA/DCM	40-4140-61	960 mL	58,00	66,70	-	-	
Glen reagent 3% TCA/DCM	40-4140-61E	960 mL	58,00	66,70	-	-	
Glen reagent 3% TCA/DCM	40-4140-62	2000 mL	111,00	128,00	-	-	
Glen reagent 3% TCA/DCM	40-4140-68	180 mL	13,90	16,00	-	-	
Glen reagent 3% TCA/DCM	40-4140-71	1000 mL	62,00	71,30	-	-	
Glen reagent 4% Aqueous Trifluoroacetic Acid	60-4042-57	450 mL	31,70	36,40	-	-	
Glen reagent 5% Phenoxyacetic anhydride in THF	40-4212-52	200 mL	108,00	124,00	-	-	
Glen reagent 5% Phenoxyacetic anhydride in THF	40-4212-57	450 mL	231,00	266,00	-	-	
Glen reagent 5% Phenoxyacetic anhydride in THF, Expedite	40-4212-52E	200 mL	108,00	124,00	-	-	
Glen reagent 5% Phenoxyacetic anhydride in THF/Pyridine	40-4210-52	200 mL	108,00	124,00	-	-	
Glen reagent 5% Phenoxyacetic anhydride in THF/Pyridine	40-4210-57	450 mL	231,00	266,00	-	-	
Glen reagent 5-Benzylthio-1H-tetrazole (BTT)	30-3070-10	1 g	27,00	31,00	-	-	
Glen reagent 5-Benzylthio-1H-tetrazole (BTT)	30-3070-20	2 g	46,20	53,10	-	-	
Glen reagent 5-Benzylthio-1H-tetrazole (BTT)	30-3070-25	25 g	385,00	443,00	-	-	
Glen reagent 6.5% Dimethylaminopyridine in Tetrahydrofuran	40-4020-52	200 mL	32,40	37,20	-	-	
Glen reagent Acetonitrile, anhydrous	40-4050-45	60 mL	9,20	10,57	-	-	
Glen reagent Acetonitrile, anhydrous	40-4050-50	100 mL	12,30	14,10	-	-	
Glen reagent Acetonitrile, anhydrous	40-4050-53	300 mL	30,80	35,40	-	-	
Glen reagent Acetonitrile, anhydrous	40-4050-57	450 mL	38,50	44,30	-	-	
Glen reagent Acetonitrile/Melm	40-4015-71	1000 mL	112,00	129,00	-	-	
Glen reagent Acetonitrile/Pyridine (1:1)	40-4500-51	150 mL	21,60	24,80	-	-	
Glen reagent Acetonitrile/Pyridine (1:1)	40-4500-60	960 mL	93,00	106,90	-	-	
Glen reagent Acetonitrile/Pyridine, anhydrous (1:1)	40-4550-50	100 mL	18,50	21,30	-	-	
Glen reagent Acetonitrile/Pyridine, anhydrous (95:5)	40-4551-50	100 mL	18,50	21,30	-	-	
Glen reagent DCM	40-4100-62	2000 mL	93,00	106,90	-	-	
Glen reagent Dichloroacetic acid (DCA)	40-4044-50	100 mL	38,50	44,30	-	-	
Glen reagent Dichloroacetic acid (DCA)	40-4044-54	120 mL	46,20	53,10	-	-	
Glen reagent Isopropyl Phosphite/Acetonitrile/Pyridine	40-4510-45	45 mL	27,70	31,80	-	-	
Glen reagent Sublimed 1H-Tetrazole in Anhydrous Acetonitrile	30-3100-45	45 mL	30,80	35,40	-	-	
Glen reagent Sublimed 1H-Tetrazole in Anhydrous Acetonitrile	30-3100-52	200 mL	77,00	88,50	-	-	
Glen reagent Sublimed 1H-Tetrazole in Anhydrous Acetonitrile	30-3100-57	450 mL	154,00	177,00	-	-	
Glen reagent Sublimed 1H-Tetrazole in Anhydrous Acetonitrile	30-3100-62	2000 mL	586,00	674,00	-	-	
Glen reagent Sublimed 1H-Tetrazole in Anhydrous Acetonitrile	30-3102-52	200 mL	77,00	88,50	-	-	
Glen reagent Sublimed 1H-Tetrazole in Anhydrous Acetonitrile	30-3102-66	60 mL	38,50	44,30	-	-	
Glen reagent Sublimed 1H-Tetrazole in Anhydrous Acetonitrile, Expedite	30-3102-52E	200 mL	77,00	88,50	-	-	
Glen reagent Sulfurizing Reagent II	40-4037-10	1 g	38,50	44,30	-	-	
Glen reagent Sulfurizing Reagent II	40-4037-20	2 g	77,00	88,50	-	-	
Glen reagent Sulfurizing Solution	40-4137-51	150 mL	77,00	88,50	-	-	
Glen reagent Sulfurizing Solution	40-4137-52	200 mL	154,00	177,00	-	-	
Glen reagent Sulfurizing Solution	40-4137-57	450 mL	347,00	399,00	-	-	
Glen reagent Tetrahydrofuran/ 2,6-Lutidine/ Acetic Anhydride	40-4010-52	200 mL	23,10	26,60	-	-	
Glen reagent Tetrahydrofuran/ 2,6-Lutidine/ Acetic Anhydride	40-4010-57	450 mL	56,00	64,40	-	-	

Description	Reference	Scale/Quantity	Price List 2018				Ice Shipping
			EUR	CHF	GBP	USD	
Glen reagent Tetrahydrofuran/ 2,6-Lutidine/ Acetic Anhydride	40-4010-61	960 mL	119,00	137,00	-	-	
Glen reagent Tetrahydrofuran/ 2,6-Lutidine/ Acetic Anhydride	40-4010-62	2000 mL	251,00	289,00	-	-	
Glen reagent Tetrahydrofuran/ Acetic Anhydride	40-4012-50	100 mL	15,40	17,70	-	-	
Glen reagent Tetrahydrofuran/ Acetic Anhydride	40-4012-52	200 mL	23,10	26,60	-	-	
Glen reagent Tetrahydrofuran/ Acetic Anhydride	40-4012-57	450 mL	56,00	64,40	-	-	
Glen reagent Tetrahydrofuran/ Acetic Anhydride	40-4012-62	2000 mL	212,00	244,00	-	-	
Glen reagent Tetrahydrofuran/ Acetic Anhydride	40-4012-66	60 mL	11,60	13,30	-	-	
Glen reagent Tetrahydrofuran/ Pyridine/ Acetic Anhydride	40-4110-45	45 mL	12,30	14,10	-	-	
Glen reagent Tetrahydrofuran/ Pyridine/ Acetic Anhydride	40-4110-52	200 mL	23,10	26,60	-	-	
Glen reagent Tetrahydrofuran/ Pyridine/ Acetic Anhydride	40-4110-57	450 mL	56,00	64,40	-	-	
Glen reagent Tetrahydrofuran/ Pyridine/ Acetic Anhydride	40-4110-62	2000 mL	251,00	289,00	-	-	
Loading Buffer	61-4120-52		50,00	57,50	-	-	
Oligo Modifying Reagent	50-1001-01	150 µmol	415,00	477,00	-	-	
Peptide Modifying Reagent	50-1002-01	1 g	358,00	411,00	-	-	
Saccharin 1-Methylimidazole (SMI)	30-3080-10	1 gr	23,00	-	-	-	
Glen Purifications							EA1C1Z
Adapter Rack for 96 well manifolds, Pk/1	60-0010-01		8,80	10,11	-	-	
Fluoro-Pak™ columns (with Luer adaptor), Pk/10	61-2100-10A		143,00	164,00	-	-	
Fluoro-Pak™ columns, Pk/10	61-2100-10		143,00	164,00	-	-	
Fluoro-Pak™ II columns (with Luer adaptor), Pk/10	61-4100-10A		193,00	222,00	-	-	
Fluoro-Pak™ II columns, Pk/10	61-4100-10		193,00	222,00	-	-	
Glen reagent Poly-Pak Cartridges, Pk/1	60-1100-01		7,00	8,05	-	-	
Glen reagent Poly-Pak Cartridges, Pk/10	60-1100-10		71,00	81,60	-	-	
Glen reagent Poly-Pak II Cartridge, Pk/1	60-3100-01		10,60	12,20	-	-	
Glen reagent Poly-Pak II Cartridge, Pk/10	60-3100-10		106,00	122,00	-	-	
Glen reagent Poly-Pak Packing	60-1000-05	5 g	108,00	-	-	-	
Glen reagent Poly-Pak Packing	60-1000-25	25 g	542,00	-	-	-	
Glen-Pak™ DNA 30mg 96-Well Plate	60-5400-01	Pk/1	409,00	470,00	-	-	
Glen-Pak™ 50mg DNA Purification Cartridge (For use in vacuum manifolds and high-throughput devices)	60-5000-96	1 pack of 96 pce	367,00	-	-	-	
Seal for unused wells on Adapter Rack, Pk/1	60-0020-01		17,60	20,20	-	-	
Glen Research reagents for cartridge purification							EA1D1Z
1-Methyl-PseudoUridine Phosphoramidite	10-3056-95	50µmoles	607,00	-	-	-	
2-F-dl-CE Phosphoramidite (Expedite)	10-1082-95E	50 µM	258,00	-	-	-	
3'-(6-FAM) PS	26-2961-01	0,1 grams	182,00	209,00	-	-	
3'-(6-FAM) PS	26-2961-10	1 gram	1.306,00	-	-	-	
3'-DTPA CPG	20-2937-01	0.1 g	179,00	206,00	-	-	
3'-DTPA CPG	20-2937-10	1 g	1.281,00	1.472,00	-	-	
3'-DTPA CPG	20-2937-13	1 x 10 µm	480,00	552,00	-	-	
3'-DTPA CPG	20-2937-14	1 x 15 µmol	744,00	855,00	-	-	
3'-DTPA CPG	20-2937-41	4 x 1 µmol	265,00	305,00	-	-	
3'-DTPA CPG	20-2937-42	4 x 0.2 µmol	157,00	180,00	-	-	
3'-Thiol-Modifier 6 S-S CPG	20-2938-01	0.1 g	122,00	140,00	-	-	
3'-Thiol-Modifier C6 S-S CPG	20-2936-01	0.1 g	122,00	140,00	-	-	
3'-Thiol-Modifier C6 S-S CPG	20-2936-10	1 g	859,00	987,00	-	-	
3'-Thiol-Modifier C6 S-S CPG	20-2936-13	1 x 10 µmol	322,00	370,00	-	-	
3'-Thiol-Modifier C6 S-S CPG	20-2936-14	1 x 15 µmol	501,00	576,00	-	-	
3'-Thiol-Modifier C6 S-S CPG	20-2936-41	4 x 1 µmol	179,00	206,00	-	-	
3'-Thiol-Modifier C6 S-S CPG	20-2936-42	4 x 0.2 µmol	107,00	123,00	-	-	
3'-Uaq Cap CPG	20-2980-01	0.1 g	258,00	297,00	-	-	
3'-6-Fluorescein Serinol CPG	20-2994-01	0.1 g	172,00	198,00	-	-	
3'-6-Fluorescein Serinol CPG	20-2994-10	1 g	1.424,00	1.637,00	-	-	
3'-6-Fluorescein Serinol CPG	20-2994-13	1x10 µmol	429,00	493,00	-	-	
3'-6-Fluorescein Serinol CPG	20-2994-14	1 x 15 µmol	644,00	740,00	-	-	
3'-6-Fluorescein Serinol CPG	20-2994-41	4 x 1 µmol	286,00	329,00	-	-	
3'-6-Fluorescein Serinol CPG	20-2994-42	4 x 0.2 µmol	172,00	198,00	-	-	
3'-Alkyne-Modifier Serinol CPG	20-2992-42	4x0.2µm	140,00	-	-	-	
3'-Alkyne-Modifier Serinol CPG	20-2992-10	1 gram	989,00	-	-	-	
3'-Alkyne-Modifier Serinol CPG	20-2992-41	4 x 1 µmol	217,00	-	-	-	
3'-Alkyne-Modifier Serinol CPG	20-2992-01	0.1 g	109,00	-	-	-	
3'-Alkyne-Modifier Serinol CPG	20-2992-42	4x0.2µm	140,00	-	-	-	
3'-Amino-Modifier Serinol CPG	20-2997-01	0.1 g	136,00	156,00	-	-	
3'-Amino-Modifier Serinol CPG	20-2997-10	1 g	966,00	1.110,00	-	-	
3'-Amino-Modifier Serinol CPG	20-2997-13	1 x 10 µmol	358,00	411,00	-	-	
3'-Amino-Modifier Serinol CPG	20-2997-14	1 x 15 µmol	537,00	617,00	-	-	
3'-Amino-Modifier Serinol CPG	20-2997-41	4 x 1 µmol	200,00	230,00	-	-	
3'-Amino-Modifier Serinol CPG	20-2997-42	4 x 0.2 µmol	122,00	140,00	-	-	
3'-BBQ-650® CPG	20-5934-42	4 x 0.2 µmol	115,00	-	-	-	
3'-CPR II CPG	20-2903-10	1 g	670,00	770,00	-	-	
3'-CPR II CPG	20-2903-01	0,1g	85,00	-	-	-	
3'-Propargyl-5-Me-dC CPG	20-2982-01	0,1 g	223,00	-	-	-	
3'-Protected Biotin Serinol CPG	20-2993-01	0.1 g	172,00	198,00	-	-	
3'-Protected Biotin Serinol CPG	20-2993-10	1 g	1.424,00	1.637,00	-	-	
3'-Protected Biotin Serinol CPG	20-2993-13	1 x 10 µmol	429,00	493,00	-	-	
3'-Protected Biotin Serinol CPG	20-2993-14	1 x 15 µmol	644,00	740,00	-	-	
3'-Protected Biotin Serinol CPG	20-2993-41	4 x 1 µmol	286,00	329,00	-	-	
3'-Protected Biotin Serinol CPG	20-2993-42	4 x 0.2 µmol	172,00	198,00	-	-	
3'-Protected BiotinLC Serinol CPG	20-2995-01	0.1 g	172,00	198,00	-	-	
3'-Protected BiotinLC Serinol CPG	20-2995-10	1 g	1.424,00	1.637,00	-	-	
3'-Protected BiotinLC Serinol CPG	20-2995-13	1 x 10 µmol	429,00	493,00	-	-	
3'-Protected BiotinLC Serinol CPG	20-2995-14	1 x 15 µmol	644,00	740,00	-	-	
3'-Protected BiotinLC Serinol CPG	20-2995-41	4 x 1 µmol	286,00	329,00	-	-	
3'-Protected BiotinLC Serinol CPG	20-2995-42	4 x 0.2 µmol	172,00	198,00	-	-	
3'-Protected BiotinLC Serinol CPG (Expedite)	20-2995-42E	4 x 0.2 µm	172,00	-	-	-	
3'-PT-Amino-Modifier C6 CPG, Expedite	20-2956-42E	4x0.2µm	123,00	-	-	-	
3'-PT-Amino-Modifier C6 PS	26-2956-10	1.0gram	1.281,00	-	-	-	
3'-Thiol-Modifier 6 S-S CPG	20-2938-10	1 g	862,00	991,00	-	-	
3'-Thiol-Modifier 6 S-S CPG	20-2938-42	4 x 0.2µm	108,00	-	-	-	
3'-Thiol-Modifier 6 S-S CPG	20-2938-42E	4 x 0.2µm	113,00	130,00	-	-	
3'-Thiol-Modifier 6 S-S CPG, Expedite	20-2938-41E	4x1,0µm	181,00	-	-	-	
5' - DMS(O)MT-Amino-Modifier C6, Expedite	10-1907-02E	0,25 grams	300,00	345,00	-	-	
5' Stearyl Phosphoramidite	10-1979-90	100 µM	75,00	-	-	-	
5'-Amino-Modifier C6-PDA	10-1947-02	0.25 g	124,00	-	-	-	

Description	Reference	Scale/Quantity	Price List 2018				Ice Shipping
			EUR	CHF	GBP	USD	
5'-Amino-Modifier C6-PDA	10-1947-02	0.25 g	124,00	-	-	-	
5'-Amino-Modifier TEG CE-Phosphoramidite	10-1917-02	0.25 g	719,00	826,00	-	-	
5'-Amino-Modifier TEG CE-Phosphoramidite	10-1917-90	100 µmol	165,00	190,00	-	-	
5'-AminoOxy - Modifier 11 - CE - Phosphoramidite	10-1919-95	50 µmoles	203,00	-	-	-	
5'-AminoOxy - Modifier 11 - CE Phosphoramidite	10-1919-90	100 µmoles	325,00	-	-	-	
5'-AminoOxy-Modifier-11-CE Phosphoramidite	10-1919-02	0.25 grams	1.105,00	-	-	-	
5'-BBQ-650 [®] -CE Phosphoramidite	10-5934-90	100 µmol	438,00	-	-	-	✓
5-Carboxy-dC-CE Phosphoramidite	10-1066-02	0.25 g	1.725,00	1.983,00	-	-	
5-Carboxy-dC-CE Phosphoramidite	10-1066-90	100 µmol	647,00	-	-	-	✓
5'-Dichloro-dimethoxy-Fluorescein Phosphoramidite II	10-5906-90		248,00	-	-	-	
5'-Dimethoxytrityl-N-benzoyl-5-cyanoethoxy-methyl-2'-deoxyCytidine,3'-[(2-cyanoethyl)-(N,N-diisopropyl)]-phosphoramidite	10-1062-95	50 µmol	480,00	552,00	421,00	-	
5'-Dimethoxytrityl-N-benzoyl-5-cyanoethoxy-methyl-2'-deoxyCytidine,3'-[(2-cyanoethyl)-(N,N-diisopropyl)]-phosphoramidite	10-1062-02	0.25 g	2.397,00	2.755,00	2.103,00	-	✓
5'-DMS(O)MT-Amino-Modifier C6, Expedite	10-19077-02e	0,25 grams	289,00	-	-	-	
5-Ethynyl-dU-CE Phosphoramidite	10-1554-02	0.25 g	960,00	-	-	-	
5-Ethynyl-dU-CE Phosphoramidite 100µmoles	10-1554-90		305,00	-	-	-	
5-Formyl dC III CE Phosphoramidite	10-1564-90	100 µmoles	875,00	-	-	-	
5-Formyl-dC-CE Phosphoramidite	10-1514-02	0.25 g	4.635,00	5.328,00	-	-	
5-Hydroxymethyl-dC II-CE Phosphoramidite	10-1510-90	100 µmoles	838,00	-	-	-	
5-Hydroxymethyl-dC II-CE Phosphoramidite	10-1510-95	50 µmol	496,00	-	-	-	✓
5-Hydroxymethyl-dC-CE Ph.amidite	10-1062-90	100 µmol	930,00	1.069,00	-	-	
5-I-dU-CE Phosphoramidite (Expedite)	10-1091-90E	100 µM	86,00	-	-	-	
5'-Maleimide-Modifier Phosphoramidite	10-1938-02	0,25 grams	419,00	482,00	-	-	
5'-Maleimide-Modifier-Phosphoramidite	10-1938-90	100 µmoles	101,00	-	-	-	
5'-O-Dimethoxytrityl-1-O-tert-butylidimethylsilyl-2-deoxyribose-3'-[(2-cyanoethyl)-(N,N-diisopropyl)]-phosphoramidite	10-1927-02	250 mg	680,00	782,00	-	-	
5'-O-Dimethoxytrityl-1-O-tert-butylidimethylsilyl-2-deoxyribose-3'-[(2-cyanoethyl)-(N,N-diisopropyl)]-phosphoramidite	10-1927-90	100 µmol	215,00	247,00	-	-	
5'-O-Dimethoxytrityl-1-O-tert-butylidimethylsilyl-2-deoxyribose-3'-[(2-cyanoethyl)-(N,N-diisopropyl)]-phosphoramidite	10-1927-95	50 µmol	115,00	132,00	-	-	
6-Fluorescein Serinol Phosphoramidite	10-1994-02	0.25 g	852,00	979,00	-	-	
6-Fluorescein Serinol Phosphoramidite	10-1994-90	100 µmol	422,00	485,00	-	-	
6-Fluorescein Serinol Phosphoramidite	10-1994-95	50 µmol	236,00	271,00	-	-	
Ac-5-Me-dC-CE Phosphoramidite	10-1560-90	100 µmoles	63,00	-	-	-	
Alkyne-Modifier Serinol Phosphoramidite	10-1992-02	0.25 g	711,00	-	-	-	
Alkyne-NHS Ester	50-1905-24	23 mg	429,00	-	-	-	✓
Amino-Modifier Serinol Phosphoramidite	10-1997-02	0.25 g	852,00	979,00	-	-	
Amino-Modifier Serinol Phosphoramidite	10-1997-90	100 µmol	322,00	370,00	-	-	
Amino-Modifier Serinol Phosphoramidite	10-1997-95	50 µmol	179,00	206,00	-	-	
Azidobutyrate NHS Ester	50-1904-23	2.3 mg	86,00	98,90	-	-	
Azidobutyrate NHS Ester	50-1904-24	23 mg	429,00	493,00	-	-	
Azobenzene Phosphoramidite	10-5800-02	0.25grams	769,00	-	-	-	✓
Azobenzene Phosphoramidite	10-5800-90	100 µM	288,00	-	-	-	✓
baseclick Oligo-Click-M-Reload	50-2100-01	1 kit	168,00	-	-	-	
BBQ-650 [®] CPG	20-5934-01	0.1 g	235,00	-	-	-	✓
BBQ-650 [®] -dT-CE Phosphoramidite	10-5944-90	100 µmol	799,00	-	-	-	
Biotin PaTP (10mM)	81-3525-02	0.025 mL	644,00	740,00	-	-	
BiotinTEG Azide	50-2000-90	100 µmol	647,00	744,00	-	-	✓
BiotinTEG Azide	50-2000-92	25 µmol	216,00	248,00	-	-	✓
C8-Alkyne-dT-CE Phosphoramidite	10-1540-90	100µM	470,00	540,00	-	-	
Coumarin Azide	50-2004-90	100 µmol	503,00	-	-	-	✓
Cy3 [™] Phosphoramidite - Expedite	10-5913-95E	50 µmol	321,00	369,00	-	-	✓
dA-CE Phosphoramidite, Expedite	10-1000-C5E	0,5 grams	7,80	-	-	-	
DBCO-dT-CE Phosphoramidite	10-1539-90	100 µMol	589,00	-	-	-	
DBCO-PEG4-phosphoramidite	10-1941-90	100 µmol	298,00	-	-	-	
DesthiobiotinTEG Azide	50-2001-90	100 µmol	575,00	661,00	-	-	✓
DesthiobiotinTEG Azide	50-2001-92	25 µmol	194,00	223,00	-	-	✓
dG-CE Phosphoramidite, Expedite	10-1020-C5E	0,5 grams	7,80	-	-	-	
Dithiol Serinol Phosphoramidite	10-1991-02	0.25 grams	818,00	-	-	-	
Dithiol Serinol Phosphoramidite	10-1991-90	100 µmoles	269,00	-	-	-	
dmf-dG-5'-CE Phosphoramidite	10-9201-02	0.25 g	91,00	104,60	-	-	
dmf-dG-5'-CE Phosphoramidite	10-9201-05	0.5 g	182,00	209,00	-	-	
dmf-dG-5'-CE Phosphoramidite	10-9201-10	1 g	363,00	417,00	-	-	
dmf-dG-CE Phosphoramidite, Expedite	10-1029-C5E	0,5 grams	7,80	-	-	-	
DyLight DY547 Phosphoramidite	10-5917-90	100 µmol	480,00	552,00	-	-	✓
DyLight DY547 Phosphoramidite	10-5917-95	50 µmol	251,00	289,00	-	-	✓
DyLight DY647 Phosphoramidite	10-5918-02	0.25 g	1.424,00	1.637,00	-	-	✓
DyLight DY647 Phosphoramidite	10-5918-90	100 µmol	480,00	552,00	-	-	✓
DyLight DY647 Phosphoramidite	10-5918-95	50 µmol	251,00	289,00	-	-	✓
F-DMT-dT-CE Phosphoramidite Pharmacia	10-1430-90P	100 µmol	107,00	123,00	-	-	✓
Ferrocene-dT-CE Phosphoramidite	10-1576-90	100 µmoles	461,00	-	-	-	
Glen reagent 1-Me-A-CE Phosphoramidite	10-3501-90	100 µmol	544,00	625,00	-	-	
Glen reagent 1-Me-dA-CE Phosphoramidite	10-1501-02	0.25 g	1.073,00	1.233,00	-	-	
Glen reagent 1-Me-dA-CE Phosphoramidite	10-1501-90	100 µmol	358,00	411,00	-	-	
Glen reagent 1-Me-dA-CE Phosphoramidite	10-1501-95	50 µmol	179,00	206,00	-	-	
Glen reagent 2,2'-Dipicolylamine Phosphoramidite	10-5801-02	0.25 g	895,00	1.029,00	-	-	✓
Glen reagent 2,2'-Dipicolylamine Phosphoramidite	10-5801-90	100 µmol	286,00	329,00	-	-	
Glen reagent 2,2'-Dipicolylamine Phosphoramidite	10-5801-95	50 µmol	150,00	172,00	-	-	✓
Glen reagent 2',3'-ddA-CE Phosphoramidite	10-7001-02	0.25 g	780,00	897,00	-	-	
Glen reagent 2',3'-ddA-CE Phosphoramidite	10-7001-90	100 µmol	186,00	214,00	-	-	
Glen reagent 2',3'-ddC-CE Phosphoramidite	10-7101-02	0.25 g	780,00	897,00	-	-	
Glen reagent 2',3'-ddC-CE Phosphoramidite	10-7101-90	100 µmol	186,00	214,00	-	-	
Glen reagent 2',3'-ddC-CPG	20-2017-01	0.1 g	429,00	493,00	-	-	
Glen reagent 2',3'-ddC-CPG	20-2117-41	4 x 1 µmol	859,00	987,00	-	-	
Glen reagent 2',3'-ddC-CPG	20-2117-42	4 x 0.2 µmol	286,00	329,00	-	-	
Glen reagent 2',3'-ddG-CE Phosphoramidite	10-7201-02	0.25 g	1.396,00	1.605,00	-	-	
Glen reagent 2',3'-ddG-CE Phosphoramidite	10-7201-90	100 µmol	329,00	378,00	-	-	
Glen reagent 2',3'-ddT-CE Phosphoramidite	10-7301-02	0.25 g	780,00	897,00	-	-	
Glen reagent 2',3'-ddT-CE Phosphoramidite	10-7301-90	100 µmol	186,00	214,00	-	-	
Glen reagent 2,6-Diaminopurine riboside a-thiotriphosphate (0.25mM) (DTTPaS)	80-3305-01	0.1 mL	143,00	164,00	-	-	
Glen reagent 2,6-Diaminopurine-TOM-CE Phosphoramidite	10-3085-90	100 µmol	608,00	699,00	-	-	
Glen reagent 2,6-Diaminopurine-TOM-CE Phosphoramidite	10-3085-95	50 µmol	304,00	349,00	-	-	
Glen reagent 2-Amino-dA-CE Phosphoramidite	10-1085-02	0.25 g	358,00	411,00	-	-	
Glen reagent 2-Amino-dA-CE Phosphoramidite	10-1085-90	100 µmol	179,00	206,00	-	-	

Description	Reference	Scale/Quantity	Price List 2018			
			EUR	CHF	GBP	USD
Glen reagent 2-Amino-dA-CE Phosphoramidite	10-1085-95	50 µmol	100,00	115,00	-	-
Glen reagent 2-Aminopurine riboside a-thiotriphosphate (20mM) (2-AP-TTPaS)	80-3304-01	0.1 mL	143,00	164,00	-	-
Glen reagent 2-Aminopurine-CE Phosphoramidite	10-1046-02	0.25 g	508,00	584,00	-	-
Glen reagent 2-Aminopurine-CE Phosphoramidite	10-1046-90	100 µmol	193,00	222,00	-	-
Glen reagent 2'-deoxyAdenosine a-thiotriphosphate (15mM) (2'-dATTPaS)	80-1000-01	0.1 mL	143,00	164,00	-	-
Glen reagent 2'-deoxyCytidine a-thiotriphosphate (15mM) (2'-dCTTPaS)	80-1010-01	0.1 mL	143,00	164,00	-	-
Glen reagent 2'-deoxyGuanosine a-thiotriphosphate (5mM) (2'-dGTTPaS)	80-1020-01	0.1 mL	143,00	164,00	-	-
Glen reagent 2'-DeoxyNebularine-CE Phosphoramidite (Purine)	10-1041-02	0.25 g	365,00	420,00	-	-
Glen reagent 2'-DeoxyNebularine-CE Phosphoramidite (Purine)	10-1041-90	100 µmol	150,00	172,00	-	-
Glen reagent 2'-deoxypseudoU-CE Phosphoramidite	10-1055-02	0.25 g	1.396,00	1.605,00	-	-
Glen reagent 2'-deoxypseudoU-CE Phosphoramidite	10-1055-90	100 µmol	508,00	584,00	-	-
Glen reagent 2'-deoxypseudoU-CE Phosphoramidite	10-1055-95	50 µmol	254,00	292,00	-	-
Glen reagent 2'-deoxyUridine a-thiotriphosphate (5mM) (2'dUTTP)	80-1040-01	0.1 mL	143,00	164,00	-	-
Glen reagent 2'-F-A-CE Phosphoramidite	10-3400-02	0.25 g	179,00	206,00	-	-
Glen reagent 2'-F-A-CE Phosphoramidite	10-3400-90	100 µmol	72,00	82,80	-	-
Glen reagent 2'-F-dl-CE Phosphoramidite	10-1082-02	0.25 g	1.396,00	1.605,00	-	-
Glen reagent 2'-F-dl-CE Phosphoramidite	10-1082-90	100 µmol	515,00	592,00	-	-
Glen reagent 2'-F-dl-CE Phosphoramidite	10-1082-95	50 µmol	258,00	297,00	-	-
Glen reagent 2'-F-G-CE Phosphoramidite	10-3420-02	0.25 g	179,00	206,00	-	-
Glen reagent 2'-F-G-CE Phosphoramidite	10-3420-90	100 µmol	72,00	82,80	-	-
Glen reagent 2'-OMe-A-CE Phosphoramidite	10-3100-SP	2g / 60mL bottle	400,00	460,00	-	-
Glen reagent 2'-OMe-G-CE Phosphor.	10-3121-SP	2g / 60mL bottle	400,00	460,00	-	-
Glen reagent 2'-OMe-Pac-A-CE Phosphoramidite	10-3601-10	1 g	358,00	411,00	-	-
Glen reagent 2'-O-Methyl-Uridine thiotriphosphate (20mM)	80-1141-01	0.1 mL	143,00	164,00	-	-
Glen reagent 2'-OMe-U-RNA-CPG	20-3630-01	0.1 g	57,00	65,50	-	-
Glen reagent 2-Thio-dT-CE Phosphoramidite	10-1036-02	0.25 g	966,00	1.110,00	-	-
Glen reagent 2-Thio-dT-CE Phosphoramidite	10-1036-90	100 µmol	422,00	485,00	-	-
Glen reagent 2-Thio-dT-CE Phosphoramidite	10-1036-95	50 µmol	236,00	271,00	-	-
Glen reagent 3'-(6-FAM) CPG	20-2961-01	0.1 g	172,00	198,00	-	-
Glen reagent 3'-(6-FAM) CPG	20-2961-10	1 g	1.424,00	1.637,00	-	-
Glen reagent 3'-(6-FAM) CPG	20-2961-13	1 x 10 µmol	429,00	493,00	-	-
Glen reagent 3'-(6-FAM) CPG	20-2961-14	1 x 15 µmol	644,00	740,00	-	-
Glen reagent 3'-(6-FAM) CPG	20-2961-41	4 x 1 µmol	286,00	329,00	-	-
Glen reagent 3'-(6-FAM) CPG	20-2961-42	4 x 0.2 µmol	172,00	198,00	-	-
Glen reagent 3'-(6-FAM) PS	26-2961-52	10 x 200 nmol	429,00	493,00	-	-
Glen reagent 3'-(6-FAM) PS	26-2961-55	10 x 40 nmol	429,00	493,00	-	-
Glen reagent 3'-(6-Fluorescein) CPG	20-2964-01	0.1 g	172,00	198,00	-	-
Glen reagent 3'-(6-Fluorescein) CPG	20-2964-10	1 g	1.424,00	1.637,00	-	-
Glen reagent 3'-(6-Fluorescein) CPG	20-2964-13	1 x 10 µmol	429,00	493,00	-	-
Glen reagent 3'-(6-Fluorescein) CPG	20-2964-14	1 x 15 µmol	644,00	740,00	-	-
Glen reagent 3'-(6-Fluorescein) CPG	20-2964-41	4 x 1 µmol	286,00	329,00	-	-
Glen reagent 3'-(6-Fluorescein) CPG	20-2964-42	4 x 0.2 µmol	172,00	198,00	-	-
Glen reagent 3'-Acridine CPG	20-2973-01	0.1 g	172,00	198,00	-	-
Glen reagent 3'-Acridine CPG	20-2973-10	1 g	1.424,00	1.637,00	-	-
Glen reagent 3'-Acridine CPG	20-2973-13	1 x 10 µmol	429,00	493,00	-	-
Glen reagent 3'-Acridine CPG	20-2973-14	1 x 15 µmol	644,00	740,00	-	-
Glen reagent 3'-Acridine CPG	20-2973-41	4 x 1 µmol	286,00	329,00	-	-
Glen reagent 3'-Acridine CPG	20-2973-42	4 x 0.2 µmol	172,00	198,00	-	-
Glen reagent 3'-Amino-Modifier C6 dC CPG	20-2019-01	0.1 g	172,00	198,00	-	-
Glen reagent 3'-Amino-Modifier C6 dC CPG	20-2019-10	1 g	1.424,00	1.637,00	-	-
Glen reagent 3'-Amino-Modifier C6 dC CPG	20-2019-13	1 x 10 µmol	429,00	493,00	-	-
Glen reagent 3'-Amino-Modifier C6 dC CPG	20-2019-14	1 x 15 µmol	644,00	740,00	-	-
Glen reagent 3'-Amino-Modifier C6 dC CPG	20-2019-41	4 x 1 µmol	286,00	329,00	-	-
Glen reagent 3'-Amino-Modifier C6 dC CPG	20-2019-42	4 x 0.2 µmol	172,00	198,00	-	-
Glen reagent 3'-Amino-Modifier C6 dT CPG	20-2039-01	0.1 g	137,00	157,00	-	-
Glen reagent 3'-Amino-Modifier C6 dT CPG	20-2039-10	1 g	1.145,00	1.316,00	-	-
Glen reagent 3'-Amino-Modifier C6 dT CPG	20-2039-13	1 x 10 µmol	344,00	395,00	-	-
Glen reagent 3'-Amino-Modifier C6 dT CPG	20-2039-14	1 x 15 µmol	515,00	592,00	-	-
Glen reagent 3'-Amino-Modifier C6 dT CPG	20-2039-41	4 x 1 µmol	229,00	263,00	-	-
Glen reagent 3'-Amino-Modifier C6 dT CPG	20-2039-42	4 x 0.2 µmol	137,00	157,00	-	-
Glen reagent 3'-Amino-Modifier C7 CPG	20-2957-01	0.1 g	136,00	156,00	-	-
Glen reagent 3'-Amino-Modifier C7 CPG	20-2957-10	1 g	966,00	1.110,00	-	-
Glen reagent 3'-Amino-Modifier C7 CPG	20-2957-13	1 x 10 µmol	358,00	411,00	-	-
Glen reagent 3'-Amino-Modifier C7 CPG	20-2957-14	1 x 15 µmol	537,00	617,00	-	-
Glen reagent 3'-Amino-Modifier C7 CPG	20-2957-41	4 x 1 µmol	200,00	230,00	-	-
Glen reagent 3'-Amino-Modifier C7 CPG	20-2957-42	4 x 0.2 µmol	122,00	140,00	-	-
Glen reagent 3'-Amino-Modifier C7 CPG 1000	20-2958-01	0.1 g	136,00	156,00	-	-
Glen reagent 3'-Amino-Modifier C7 CPG 1000	20-2958-10	1 g	966,00	1.110,00	-	-
Glen reagent 3'-Amino-Modifier C7 CPG 1000	20-2958-13	1 x 10 µmol	358,00	411,00	-	-
Glen reagent 3'-Amino-Modifier C7 CPG 1000	20-2958-14	1 x 15 µmol	537,00	617,00	-	-
Glen reagent 3'-Amino-Modifier C7 CPG 1000	20-2958-41	4 x 1 µmol	200,00	230,00	-	-
Glen reagent 3'-Amino-Modifier C7 CPG 1000	20-2958-42	4 x 0.2 µmol	122,00	140,00	-	-
Glen reagent 3'-BHQ-1 CPG	20-5931-01	0.1 g	272,00	313,00	-	-
Glen reagent 3'-BHQ-1 CPG	20-5931-10	1 g	2.147,00	2.468,00	-	-
Glen reagent 3'-BHQ-1 CPG	20-5931-13	1 x 10 µmol	823,00	946,00	-	-
Glen reagent 3'-BHQ-1 CPG	20-5931-14	1 x 15 µmol	1.181,00	1.357,00	-	-
Glen reagent 3'-BHQ-1 CPG	20-5931-41	4 x 1 µmol	429,00	493,00	-	-
Glen reagent 3'-BHQ-1 CPG	20-5931-42	4 x 0.2 µmol	115,00	132,00	-	-
Glen reagent 3'-BHQ-2 CPG	20-5932-01	0.1 g	272,00	313,00	-	-
Glen reagent 3'-BHQ-2 CPG	20-5932-10	1 g	2.147,00	2.468,00	-	-
Glen reagent 3'-BHQ-2 CPG	20-5932-13	1 x 10 µmol	823,00	946,00	-	-
Glen reagent 3'-BHQ-2 CPG	20-5932-14	1 x 15 µmol	1.181,00	1.357,00	-	-
Glen reagent 3'-BHQ-2 CPG	20-5932-41	4 x 1 µmol	429,00	493,00	-	-
Glen reagent 3'-BHQ-2 CPG	20-5932-42	4 x 0.2 µmol	115,00	132,00	-	-
Glen reagent 3'-BHQ-3 CPG	20-5933-01	0.1 g	272,00	313,00	-	-
Glen reagent 3'-BHQ-3 CPG	20-5933-10	1 g	2.147,00	2.468,00	-	-
Glen reagent 3'-BHQ-3 CPG	20-5933-13	1 x 10 µmol	823,00	946,00	-	-
Glen reagent 3'-BHQ-3 CPG	20-5933-14	1 x 15 µmol	1.181,00	1.357,00	-	-
Glen reagent 3'-BHQ-3 CPG	20-5933-41	4 x 1 µmol	429,00	493,00	-	-
Glen reagent 3'-BHQ-3 CPG	20-5933-42	4 x 0.2 µmol	115,00	132,00	-	-
Glen reagent 3'-BiotinTEG CPG	20-2955-01	0.1 g	260,00	299,00	-	-
Glen reagent 3'-BiotinTEG CPG	20-2955-10	1 g	2.155,00	2.477,00	-	-

Description	Reference	Scale/Quantity	Price List 2018				Ice Shipping
			EUR	CHF	GBP	USD	
Glen reagent 3'-BiotinTEG CPG	20-2955-13	1 x 10 µmol	650,00	747,00	-	-	
Glen reagent 3'-BiotinTEG CPG	20-2955-14	1 x 15 µmol	975,00	1.121,00	-	-	
Glen reagent 3'-BiotinTEG CPG	20-2955-41	4 x 1 µmol	433,00	498,00	-	-	
Glen reagent 3'-BiotinTEG CPG	20-2955-42	4 x 0.2 µmol	260,00	299,00	-	-	
Glen reagent 3'-BiotinTEG PS	26-2955-52	10 x 200 nmol	429,00	493,00	-	-	
Glen reagent 3'-BiotinTEG PS	26-2955-55	10 x 40 nmol	429,00	493,00	-	-	
Glen reagent 3'-Cholesteryl-TEG CPG	20-2975-01	0.1 g	122,00	-	-	-	
Glen reagent 3'-Cholesteryl-TEG CPG	20-2975-10	1 g	1.002,00	1.152,00	-	-	
Glen reagent 3'-Cholesteryl-TEG CPG	20-2975-13	1 x 10 µmol	301,00	-	-	-	
Glen reagent 3'-Cholesteryl-TEG CPG	20-2975-14	1 x 15 µmol	451,00	-	-	-	
Glen reagent 3'-Cholesteryl-TEG CPG	20-2975-41	4 x 1 µmol	200,00	-	-	-	
Glen reagent 3'-Cholesteryl-TEG CPG	20-2975-42	4 x 0.2 µmol	120,00	-	-	-	
Glen reagent 3'-Cholesteryl-TEG CPG	20-2975-10	1 g	1.002,00	1.152,00	-	-	
Glen reagent 3'-Dabcyl PS	26-5912-52	10 x 200 nmol	429,00	493,00	-	-	
Glen reagent 3'-Dabcyl PS	26-5912-55	10 x 40 nmol	429,00	493,00	-	-	
Glen reagent 3'-Dabsyl CPG	20-5911-01	0.1 g	172,00	198,00	-	-	
Glen reagent 3'-Dabsyl CPG	20-5911-10	1 g	1.396,00	1.605,00	-	-	
Glen reagent 3'-Dabsyl CPG	20-5911-13	1 x 10 µmol	501,00	576,00	-	-	
Glen reagent 3'-Dabsyl CPG	20-5911-14	1 x 15 µmol	716,00	823,00	-	-	
Glen reagent 3'-Dabsyl CPG	20-5911-41	4 x 1 µmol	286,00	329,00	-	-	
Glen reagent 3'-Dabsyl CPG	20-5911-42	4 x 0.2 µmol	172,00	198,00	-	-	
Glen reagent 3'-Dabsyl CPG	20-5912-01	0.1 g	172,00	198,00	-	-	
Glen reagent 3'-Dabsyl CPG	20-5912-10	1 g	1.396,00	1.605,00	-	-	
Glen reagent 3'-Dabsyl CPG	20-5912-13	1 x 10 µmol	501,00	576,00	-	-	
Glen reagent 3'-Dabsyl CPG	20-5912-14	1 x 15 µmol	716,00	823,00	-	-	
Glen reagent 3'-Dabsyl CPG	20-5912-41	4 x 1 µmol	286,00	329,00	-	-	
Glen reagent 3'-Dabsyl CPG	20-5912-42	4 x 0.2 µmol	172,00	198,00	-	-	
Glen reagent 3'-Dabsyl CPG	20-5912-SP	4 x 1.3 µmol	333,00	383,00	-	-	
Glen reagent 3'-dA-CE Phosphoramidite	10-1004-02	0.25 g	1.396,00	1.605,00	-	-	
Glen reagent 3'-dA-CE Phosphoramidite	10-1004-90	100 µmol	508,00	584,00	-	-	
Glen reagent 3'-dA-CE Phosphoramidite	10-1004-95	50 µmol	254,00	292,00	-	-	
Glen reagent 3'-dA-CPG	20-2004-01	0.1 g	429,00	493,00	-	-	
Glen reagent 3'-dA-CPG	20-2104-41	4 x 1 µmol	859,00	987,00	-	-	
Glen reagent 3'-dA-CPG	20-2104-42	4 x 0.2 µmol	286,00	329,00	-	-	
Glen reagent 3'-dC-CE Phosphoramidite	10-1064-02	0.25 g	1.396,00	1.605,00	-	-	
Glen reagent 3'-dC-CE Phosphoramidite	10-1064-90	100 µmol	508,00	584,00	-	-	
Glen reagent 3'-dC-CE Phosphoramidite	10-1064-95	50 µmol	254,00	292,00	-	-	
Glen reagent 3'-dC-CPG	20-2064-01	0.1 g	429,00	493,00	-	-	
Glen reagent 3'-dC-CPG	20-2164-41	4 x 1 µmol	859,00	987,00	-	-	
Glen reagent 3'-dC-CPG	20-2164-42	4 x 0.2 µmol	286,00	329,00	-	-	
Glen reagent 3'-deaza-dA-CE Phosphoramidite	10-1088-02	0.25 g	1.396,00	1.605,00	-	-	
Glen reagent 3'-deaza-dA-CE Phosphoramidite	10-1088-90	100 µmol	508,00	584,00	-	-	
Glen reagent 3'-deaza-dA-CE Phosphoramidite	10-1088-95	50 µmol	254,00	292,00	-	-	
Glen reagent 3'-dG-CE Phosphoramidite	10-1074-02	0.25 g	1.396,00	1.605,00	-	-	
Glen reagent 3'-dG-CE Phosphoramidite	10-1074-90	100 µmol	508,00	584,00	-	-	
Glen reagent 3'-dG-CE Phosphoramidite	10-1074-95	50 µmol	254,00	292,00	-	-	
Glen reagent 3'-dG-CPG	20-2074-01	0.1 g	429,00	493,00	-	-	
Glen reagent 3'-dG-CPG	20-2174-41	4 x 1 µmol	859,00	987,00	-	-	
Glen reagent 3'-dG-CPG	20-2174-42	4 x 0.2 µmol	286,00	329,00	-	-	
Glen reagent 3'-dT-CE Phosphoramidite	10-1084-02	0.25 g	1.396,00	1.605,00	-	-	
Glen reagent 3'-dT-CE Phosphoramidite	10-1084-90	100 µmol	508,00	584,00	-	-	
Glen reagent 3'-dT-CE Phosphoramidite	10-1084-95	50 µmol	254,00	292,00	-	-	
Glen reagent 3'-dT-CPG	20-2084-01	0.1 g	429,00	493,00	-	-	
Glen reagent 3'-dT-CPG	20-2184-41	4 x 1 µmol	859,00	987,00	-	-	
Glen reagent 3'-dT-CPG	20-2184-42	4 x 0.2 µmol	286,00	329,00	-	-	
Glen reagent 3'-Fluorescein CPG	20-2963-01	0.1 g	172,00	198,00	-	-	
Glen reagent 3'-Fluorescein CPG	20-2963-10	1 g	1.424,00	1.637,00	-	-	
Glen reagent 3'-Fluorescein CPG	20-2963-13	1 x 10 µmol	429,00	493,00	-	-	
Glen reagent 3'-Fluorescein CPG	20-2963-14	1 x 15 µmol	644,00	740,00	-	-	
Glen reagent 3'-Fluorescein CPG	20-2963-41	4 x 1 µmol	286,00	329,00	-	-	
Glen reagent 3'-Fluorescein CPG	20-2963-42	4 x 0.2 µmol	172,00	198,00	-	-	
Glen reagent 3'-Fluorescein-dT CPG	20-2056-01	0.1 g	172,00	198,00	-	-	
Glen reagent 3'-Fluorescein-dT CPG	20-2056-10	1 g	1.424,00	1.637,00	-	-	
Glen reagent 3'-Fluorescein-dT CPG	20-2056-13	1 x 10 µmol	429,00	493,00	-	-	
Glen reagent 3'-Fluorescein-dT CPG	20-2056-14	1 x 15 µmol	644,00	740,00	-	-	
Glen reagent 3'-Fluorescein-dT CPG	20-2056-41	4 x 1 µmol	286,00	329,00	-	-	
Glen reagent 3'-Fluorescein-dT CPG	20-2056-42	4 x 0.2 µmol	172,00	198,00	-	-	
Glen reagent 3'-Glyceryl CPG	20-2902-01	0.1 g	122,00	140,00	-	-	
Glen reagent 3'-Glyceryl CPG	20-2902-10	1 g	859,00	987,00	-	-	
Glen reagent 3'-Glyceryl CPG	20-2902-13	1 x 10 µmol	322,00	370,00	-	-	
Glen reagent 3'-Glyceryl CPG	20-2902-14	1 x 15 µmol	501,00	576,00	-	-	
Glen reagent 3'-Glyceryl CPG	20-2902-41	4 x 1 µmol	179,00	206,00	-	-	
Glen reagent 3'-Glyceryl CPG	20-2902-42	4 x 0.2 µmol	107,00	123,00	-	-	
Glen reagent 3'Glyceryl CPG col, Pharm.	20-2902-41P	1 µm (4 x 1 pce)	179,00	206,00	-	-	
Glen reagent 3'-Phosphate CPG	20-2900-01	0.1 g	100,00	115,00	-	-	
Glen reagent 3'-Phosphate CPG	20-2900-10	1 g	687,00	790,00	-	-	
Glen reagent 3'-Phosphate CPG	20-2900-13	1 x 10 µmol	258,00	297,00	-	-	
Glen reagent 3'-Phosphate CPG	20-2900-14	1 x 15 µmol	401,00	461,00	-	-	
Glen reagent 3'-Phosphate CPG	20-2900-41	4 x 1 µmol	143,00	164,00	-	-	
Glen reagent 3'-Phosphate CPG	20-2900-42	4 x 0.2 µmol	86,00	98,90	-	-	
Glen reagent 3'-Phosphate CPG (High Load)	25-2900-01	0.1 g	122,00	140,00	-	-	
Glen reagent 3'-Phosphate CPG (High Load)	25-2900-10	1 g	859,00	987,00	-	-	
Glen reagent 3'-Phosphate PS	26-2900-52	10 x 200 nmol	215,00	247,00	-	-	
Glen reagent 3'-Phosphate PS	26-2900-55	10 x 40 nmol	215,00	247,00	-	-	
Glen reagent 3'-PT-Amino-Modifier C3 CPG	20-2954-01	0.1 g	136,00	156,00	-	-	
Glen reagent 3'-PT-Amino-Modifier C3 CPG	20-2954-10	1 g	966,00	1.110,00	-	-	
Glen reagent 3'-PT-Amino-Modifier C3 CPG	20-2954-13	1 x 10 µmol	358,00	411,00	-	-	
Glen reagent 3'-PT-Amino-Modifier C3 CPG	20-2954-14	1 x 15 µmol	537,00	617,00	-	-	
Glen reagent 3'-PT-Amino-Modifier C3 CPG	20-2954-41	4 x 1 µmol	200,00	230,00	-	-	
Glen reagent 3'-PT-Amino-Modifier C3 CPG	20-2954-42	4 x 0.2 µmol	122,00	140,00	-	-	
Glen reagent 3'-PT-Amino-Modifier C6 CPG	20-2956-01	0.1 g	136,00	156,00	-	-	
Glen reagent 3'-PT-Amino-Modifier C6 CPG	20-2956-10	1 g	966,00	1.110,00	-	-	
Glen reagent 3'-PT-Amino-Modifier C6 CPG	20-2956-13	1 x 10 µmol	358,00	411,00	-	-	

Description	Reference	Scale/Quantity	Price List 2018				Ice Shipping
			EUR	CHF	GBP	USD	
Glen reagent 3'-PT-Amino-Modifier C6 CPG	20-2956-14	1 x 15 µmol	537,00	617,00	-	-	
Glen reagent 3'-PT-Amino-Modifier C6 CPG	20-2956-41	4 x 1 µmol	200,00	230,00	-	-	
Glen reagent 3'-PT-Amino-Modifier C6 CPG	20-2956-42	4 x 0.2 µmol	122,00	140,00	-	-	
Glen reagent 3'-PT-Amino-Modifier C6 PS	26-2956-52	10 x 200 nmol	315,00	362,00	-	-	
Glen reagent 3'-PT-Amino-Modifier C6 PS	26-2956-55	10 x 40 nmol	315,00	362,00	-	-	
Glen reagent 3'-Spacer C3 CPG	20-2913-01	0.1 g	100,00	115,00	-	-	
Glen reagent 3'-Spacer C3 CPG	20-2913-10	1 g	687,00	790,00	-	-	
Glen reagent 3'-Spacer C3 CPG	20-2913-13	1 x 10 µmol	258,00	297,00	-	-	
Glen reagent 3'-Spacer C3 CPG	20-2913-14	1 x 15 µmol	401,00	461,00	-	-	
Glen reagent 3'-Spacer C3 CPG	20-2913-41	4 x 1 µmol	143,00	164,00	-	-	
Glen reagent 3'-Spacer C3 CPG	20-2913-42	4 x 0.2 µmol	86,00	98,90	-	-	
Glen reagent 3'-TAMRA CPG	20-5910-01	0.1 g	172,00	198,00	-	-	
Glen reagent 3'-TAMRA CPG	20-5910-10	1 g	1,424,00	1,637,00	-	-	
Glen reagent 3'-TAMRA CPG	20-5910-41	4 x 1 µmol	286,00	329,00	-	-	
Glen reagent 3'-TAMRA CPG	20-5910-42	4 x 0.2 µmol	143,00	-	-	-	
Glen reagent 3'-TAMRA PS	26-5910-52	10 x 200 nmol	330,00	-	-	-	
Glen reagent 3'-TAMRA PS	26-5910-55	10 x 40 nmol	330,00	-	-	-	
Glen reagent 3'-Thiol-Modifier C3 S-S CPG	20-2933-01	0.1 g	122,00	140,00	-	-	
Glen reagent 3'-Thiol-Modifier C3 S-S CPG	20-2933-10	1 g	859,00	987,00	-	-	
Glen reagent 3'-Thiol-Modifier C3 S-S CPG	20-2933-13	1 x 10 µmol	322,00	370,00	-	-	
Glen reagent 3'-Thiol-Modifier C3 S-S CPG	20-2933-14	1 x 15 µmol	501,00	576,00	-	-	
Glen reagent 3'-Thiol-Modifier C3 S-S CPG	20-2933-41	4 x 1 µmol	179,00	206,00	-	-	
Glen reagent 3'-Thiol-Modifier C3 S-S CPG	20-2933-42	4 x 0.2 µmol	107,00	123,00	-	-	
Glen reagent 4-Thio-dT-CE Phosphoramidite	10-1034-02	0.25 g	966,00	1,110,00	-	-	
Glen reagent 4-Thio-dT-CE Phosphoramidite	10-1034-90	100 µmol	422,00	485,00	-	-	
Glen reagent 4-Thio-dT-CE Phosphoramidite	10-1034-95	50 µmol	236,00	271,00	-	-	
Glen reagent 4-Thio-dU-CE Phosphoramidite	10-1052-02	0.25 g	966,00	1,110,00	-	-	
Glen reagent 4-Thio-dU-CE Phosphoramidite	10-1052-90	100 µmol	422,00	485,00	-	-	
Glen reagent 4-Thio-dU-CE Phosphoramidite	10-1052-95	50 µmol	236,00	271,00	-	-	
Glen reagent 5,6-Dihydro-dT-CE Phosphoramidite	10-1530-02	0.25 g	852,00	979,00	-	-	
Glen reagent 5,6-Dihydro-dT-CE Phosphoramidite	10-1530-90	100 µmol	279,00	321,00	-	-	
Glen reagent 5,6-Dihydro-dU-CE Phosphoramidite	10-1550-02	0.25 g	852,00	979,00	-	-	
Glen reagent 5,6-Dihydro-dU-CE Phosphoramidite	10-1550-90	100 µmol	279,00	321,00	-	-	
Glen reagent 5'-Aldehyde-Modifier C2 Phosphoramidite	10-1933-02	0.25 g	465,00	534,00	-	-	
Glen reagent 5'-Aldehyde-Modifier C2 Phosphoramidite	10-1933-90	100 µmol	122,00	140,00	-	-	
Glen reagent 5'-Aldehyde-Modifier C2 Phosphoramidite, Expedite	10-1933-90E	100 µmol	122,00	-	-	-	
Glen reagent 5'-Amino-dT-CE Phosphoramidite	10-1932-02	0.25 g	573,00	659,00	-	-	
Glen reagent 5'-Amino-dT-CE Phosphoramidite	10-1932-90	100 µmol	215,00	247,00	-	-	
Glen reagent 5'-Amino-Modifier 5	10-1905-02	0.25 g	286,00	329,00	-	-	
Glen reagent 5'-Amino-Modifier 5	10-1905-90	100 µmol	86,00	98,90	-	-	
Glen reagent 5'-Amino-Modifier C12	10-1912-02	0.25 g	429,00	493,00	-	-	
Glen reagent 5'-Amino-Modifier C12	10-1912-90	100 µmol	129,00	148,00	-	-	
Glen reagent 5'-Amino-Modifier C3-TFA	10-1923-02	0.25 g	251,00	289,00	-	-	
Glen reagent 5'-Amino-Modifier C3-TFA	10-1923-90	100 µmol	72,00	82,80	-	-	
Glen reagent 5'-Amino-Modifier C6	10-1906-02	0.25 g	220,00	253,00	-	-	
Glen reagent 5'-Amino-Modifier C6	10-1906-90	100 µmol	66,00	75,90	-	-	
Glen reagent 5'-Amino-Modifier C6, Beckman	10-1906-90B	100 µmol	66,00	-	-	-	
Glen reagent 5'-Amino-Modifier C6-TFA	10-1916-02	0.25 g	89,00	102,30	78,10	106,80	
Glen reagent 5'-Amino-Modifier C6-TFA	10-1916-90	100 µmol	29,70	34,10	-	-	
Glen reagent 5'-BHQ-1 Phosphoramidite	10-5931-02	0.25 g	1,002,00	1,152,00	-	-	
Glen reagent 5'-BHQ-1 Phosphoramidite	10-5931-90	100 µmol	286,00	329,00	-	-	
Glen reagent 5'-BHQ-1 Phosphoramidite	10-5931-95	50 µmol	143,00	164,00	-	-	
Glen reagent 5'-BHQ-2 Phosphoramidite	10-5932-02	0.25 g	1,002,00	1,152,00	-	-	
Glen reagent 5'-BHQ-2 Phosphoramidite	10-5932-90	100 µmol	286,00	329,00	-	-	
Glen reagent 5'-BHQ-2 Phosphoramidite	10-5932-95	50 µmol	143,00	164,00	-	-	
Glen reagent 5'-Biotin Phosphoramidite	10-5950-02	0.25 g	573,00	659,00	-	-	
Glen reagent 5'-Biotin Phosphoramidite	10-5950-90	100 µmol	198,00	228,00	-	-	
Glen reagent 5'-Biotin Phosphoramidite	10-5950-95	50 µmol	110,00	126,00	-	-	
Glen reagent 5'-Biotin Phosphoramidite, Expedite	10-5950-95E	50 µmol	110,00	126,00	-	-	
Glen reagent 5-Br-dC-CE Phosphoramidite	10-1080-02	0.25 g	229,00	263,00	-	-	
Glen reagent 5-Br-dC-CE Phosphoramidite	10-1080-90	100 µmol	86,00	98,90	-	-	
Glen reagent 5-Br-dU-CE Phosphoramidite	10-1090-02	0.25 g	176,00	202,00	-	-	
Glen reagent 5-Br-dU-CE Phosphoramidite	10-1090-90	100 µmol	66,00	75,90	-	-	
Glen reagent 5-Br-dU-CPG	20-2090-01	0.1 g	72,00	82,80	-	-	
Glen reagent 5-Br-dU-CPG	20-2090-41	4 x 1 µmol	286,00	329,00	-	-	
Glen reagent 5-Br-dU-CPG	20-2090-42	4 x 0.2 µmol	172,00	198,00	-	-	
Glen reagent 5'-Bromohexyl Phosphoramidite	10-1946-02	0.25 g	286,00	329,00	-	-	
Glen reagent 5'-Bromohexyl Phosphoramidite	10-1946-90	100 µmol	86,00	98,90	-	-	
Glen reagent 5'-Carboxy-Modifier C10	10-1935-02	0.25 g	320,00	368,00	-	-	
Glen reagent 5'-Carboxy-Modifier C11	10-1935-90	100 µmol	94,00	108,00	-	-	
Glen reagent 5'-Cholesteryl-TEG Phosphoramidite	10-1976-02	0.25 g	751,00	863,00	-	-	
Glen reagent 5'-Cholesteryl-TEG Phosphoramidite	10-1976-90	100 µmol	250,00	-	-	-	
Glen reagent 5'-Cholesteryl-TEG Phosphoramidite	10-1976-95	50 µmol	136,00	-	-	-	
Glen reagent 5'-Cholesteryl-TEG Phosphoramidite	10-1976-02	0.25 g	751,00	863,00	-	-	
Glen reagent 5'-DabcyI Phosphoramidite	10-5912-02	0.25 g	930,00	1,069,00	-	-	
Glen reagent 5'-DabcyI Phosphoramidite	10-5912-90	100 µmol	322,00	370,00	-	-	
Glen reagent 5'-DabcyI Phosphoramidite	10-5912-95	50 µmol	179,00	206,00	-	-	
Glen reagent 5'-DabcyI-dT	10-1532-02	0.25 g	1,109,00	1,275,00	-	-	
Glen reagent 5'-DabcyI-dT	10-1532-90	100 µmol	358,00	411,00	-	-	
Glen reagent 5'-Dichloro-dimethoxy-Fluorescein Phosphoramidite (Joe)	10-5904-02	0.25 g	1,109,00	1,275,00	-	-	
Glen reagent 5'-Dichloro-dimethoxy-Fluorescein Phosphoramidite (Joe)	10-5904-90	100 µmol	487,00	560,00	-	-	
Glen reagent 5'-Dichloro-dimethoxy-Fluorescein Phosphoramidite (Joe)	10-5904-95	50 µmol	251,00	289,00	-	-	
Glen reagent 5'-DMS(O)MT-Amino-Modifier C6	10-1907-02	0.25 g	286,00	329,00	-	-	
Glen reagent 5'-DMS(O)MT-Amino-Modifier C6	10-1907-90	100 µmol	86,00	98,90	-	-	
Glen reagent 5'-F-dU-CE Phosphoramidite	10-1092-02	0.25 g	508,00	584,00	-	-	
Glen reagent 5'-F-dU-CE Phosphoramidite	10-1092-90	100 µmol	193,00	222,00	-	-	
Glen reagent 5'-Fluorescein Phosphoramidite	10-5901-90P	100µmol	124,00	143,00	-	-	
Glen reagent 5'-Fluorescein Phosphoramidite	10-5901-02	0.25 g	297,00	341,00	-	-	
Glen reagent 5'-Fluorescein Phosphoramidite	10-5901-90	100 µmol	109,00	125,00	-	-	
Glen reagent 5'-Fluorescein Phosphoramidite	10-5901-95	50 µmol	66,00	75,90	-	-	
Glen reagent 5'-Fluorescein Phosphoramidite (6-FAM)	10-5901-BULK	1 g	795,00	914,00	-	-	
Glen reagent 5'-Formylindole-CE Phosphoramidite	10-1934-02	0.25 g	465,00	534,00	-	-	
Glen reagent 5'-Formylindole-CE Phosphoramidite	10-1934-90	100 µmol	122,00	140,00	-	-	

Description	Reference	Scale/Quantity	Price List 2018				Ice Shipping
			EUR	CHF	GBP	USD	
			Glen reagent 5'-Hexachloro-Fluorescein Phosphoramidite	10-5902-02	0.25 g	870,00	
Glen reagent 5'-Hexachloro-Fluorescein Phosphoramidite	10-5902-90	100 µmol	385,00	443,00	-	-	
Glen reagent 5'-Hexachloro-Fluorescein Phosphoramidite	10-5902-95	50 µmol	198,00	228,00	-	-	
Glen reagent 5'-Hexynyl Phosphoramidite	10-1908-02	0.25 g	286,00	329,00	-	-	
Glen reagent 5'-Hexynyl Phosphoramidite	10-1908-90	100 µmol	86,00	98,90	-	-	
Glen reagent 5-Hydroxymethyl-dU-CE Phosphoramidite	10-1093-02	0.25 g	966,00	1.110,00	-	-	
Glen reagent 5-Hydroxymethyl-dU-CE Phosphoramidite	10-1093-90	100 µmol	322,00	370,00	-	-	
Glen reagent 5-I-dC-CE Phosphoramidite	10-1081-02	0.25 g	508,00	584,00	-	-	
Glen reagent 5-I-dC-CE Phosphoramidite	10-1081-90	100 µmol	193,00	222,00	-	-	
Glen reagent 5'-I-dT-CE Phosphoramidite	10-1931-02	0.25 g	422,00	485,00	-	-	
Glen reagent 5'-I-dT-CE Phosphoramidite	10-1931-90	100 µmol	122,00	140,00	-	-	
Glen reagent 5-I-dU-CE Phosphoramidite	10-1091-02	0.25 g	229,00	263,00	-	-	
Glen reagent 5-I-dU-CE Phosphoramidite	10-1091-90	100 µmol	86,00	98,90	-	-	
Glen reagent 5-I-dU-CE Phosphoramidite, Expedite	10-1091-02E	0.25 g	229,00	-	-	-	
Glen reagent 5-Me-2'-deoxyZebularine-CE Phosphoramidite	10-1061-02	0.25 g	1.396,00	1.605,00	-	-	
Glen reagent 5-Me-2'-deoxyZebularine-CE Phosphoramidite	10-1061-90	100 µmol	573,00	659,00	-	-	
Glen reagent 5-Me-2'-deoxyZebularine-CE Phosphoramidite	10-1061-95	50 µmol	286,00	329,00	-	-	
Glen reagent 5-Me-dC-CE Phosphoramidite	10-1060-02	0.25 g	132,00	152,00	-	-	
Glen reagent 5-Me-dC-CE Phosphoramidite	10-1060-90	100 µmol	55,00	63,20	-	-	
Glen reagent 5-Me-dC-CPG	20-2060-01	0.1 g	72,00	82,80	-	-	
Glen reagent 5-Me-dC-CPG	20-2160-41	4 x 1 µmol	286,00	329,00	-	-	
Glen reagent 5-Me-dC-CPG	20-2160-42	4 x 0.2 µmol	172,00	198,00	-	-	
Glen reagent 5-Methyl-Uridine a-thiotriphosphate (10mM) (5-Me-UTTPaS)	80-3093-01	0.1 mL	143,00	164,00	-	-	
Glen reagent 5-Nitroindole-CE Phosphoramidite	10-1044-02	0.25 g	465,00	534,00	-	-	
Glen reagent 5-Nitroindole-CE Phosphoramidite	10-1044-90	100 µmol	179,00	206,00	-	-	
Glen reagent 5-OH-dC-CE Phosphoramidite	10-1063-02	0.25 g	1.109,00	1.275,00	-	-	
Glen reagent 5-OH-dC-CE Phosphoramidite	10-1063-90	100 µmol	394,00	453,00	-	-	
Glen reagent 5-OH-dU-CE Phosphoramidite	10-1053-02	0.25 g	966,00	1.110,00	-	-	
Glen reagent 5-OH-dU-CE Phosphoramidite	10-1053-90	100 µmol	322,00	370,00	-	-	
Glen reagent 5'-OMe-dT-CE Phosphoramidite	10-1031-02	0.25 g	508,00	584,00	-	-	
Glen reagent 5'-OMe-dT-CE Phosphoramidite	10-1031-90	100 µmol	193,00	222,00	-	-	
Glen reagent 5'-Pyrene Cap Phosphoramidite	10-1987-02	0.25 g	709,00	815,00	-	-	
Glen reagent 5'-Pyrene Cap Phosphoramidite	10-1987-90	100 µmol	279,00	321,00	-	-	
Glen reagent 5'-Tetrachloro-Fluorescein Phosphoramidite	10-5903-02	0.25 g	879,00	-	-	-	
Glen reagent 5'-Tetrachloro-Fluorescein Phosphoramidite	10-5903-90	100 µmol	407,00	468,00	-	-	
Glen reagent 5'-Tetrachloro-Fluorescein Phosphoramidite	10-5903-95	50 µmol	218,00	251,00	-	-	
Glen reagent 5'-Thiol-Modifier C6	10-1926-02	0.25 g	286,00	329,00	-	-	
Glen reagent 5'-Thiol-Modifier C6	10-1926-90	100 µmol	86,00	98,90	-	-	
Glen reagent 5'-Thiol-Modifier C6, Expedite	10-1926-90E	100 µmol	86,00	-	-	-	
Glen reagent 5'-Trimethoxystilbene Cap Phosphoramidite	10-1986-02	0.25 g	709,00	815,00	-	-	
Glen reagent 5'-Trimethoxystilbene Cap Phosphoramidite	10-1986-90	100 µmol	279,00	321,00	-	-	
Glen reagent 6-Fluorescein Phosphoramidite	10-1964-02	0.25 g	852,00	979,00	-	-	
Glen reagent 6-Fluorescein Phosphoramidite	10-1964-90	100 µmol	422,00	485,00	-	-	
Glen reagent 6-Fluorescein Phosphoramidite	10-1964-95	50 µmol	236,00	271,00	-	-	
Glen reagent 6-thio-dG-CE Phosphoramidite	10-1072-02	0.25 g	1.396,00	1.605,00	-	-	
Glen reagent 6-thio-dG-CE Phosphoramidite	10-1072-90	100 µmol	508,00	584,00	-	-	
Glen reagent 6-thio-dG-CE Phosphoramidite	10-1072-95	50 µmol	254,00	292,00	-	-	
Glen reagent 6-Thio-G-CE Phosphoramidite	10-3072-02	0.25 g	1.718,00	1.975,00	-	-	
Glen reagent 6-Thio-G-CE Phosphoramidite	10-3072-90	100 µmol	716,00	823,00	-	-	
Glen reagent 6-Thio-G-CE Phosphoramidite	10-3072-95	50 µmol	358,00	411,00	-	-	
Glen reagent 7-deaza-8-aza-dA-CE Phosphoramidite	10-1083-02	0.25 g	1.396,00	1.605,00	-	-	
Glen reagent 7-deaza-8-aza-dA-CE Phosphoramidite	10-1083-90	100 µmol	508,00	584,00	-	-	
Glen reagent 7-deaza-8-aza-dA-CE Phosphoramidite	10-1083-95	50 µmol	254,00	292,00	-	-	
Glen reagent 7-Deaza-8-aza-dG-CE Phosphoramidite (PPG)	10-1073-02	0.25 g	1.422,00	1.634,00	-	-	
Glen reagent 7-Deaza-8-aza-dG-CE Phosphoramidite (PPG)	10-1073-90	100 µmol	488,00	561,00	-	-	
Glen reagent 7-Deaza-8-aza-dG-CE Phosphoramidite (PPG)	10-1073-95	50 µmol	257,00	295,00	-	-	
Glen reagent 7-deaza-A-CE Phosphoramidite	10-3001-02	0.25 g	1.574,00	1.809,00	-	-	
Glen reagent 7-deaza-A-CE Phosphoramidite	10-3001-90	100 µmol	644,00	740,00	-	-	
Glen reagent 7-deaza-A-CE Phosphoramidite	10-3001-95	50 µmol	322,00	370,00	-	-	
Glen reagent 7-Deaza-Adenosine a-thiotriphosphate (10mM) (7-DATTPaS)	80-3303-01	0.1 mL	143,00	164,00	-	-	
Glen reagent 7-Deaza-dA-CE Phosphoramidite	10-1001-02	0.25 g	1.396,00	1.605,00	-	-	
Glen reagent 7-Deaza-dA-CE Phosphoramidite	10-1001-90	100 µmol	508,00	584,00	-	-	
Glen reagent 7-Deaza-dA-CE Phosphoramidite	10-1001-95	50 µmol	254,00	292,00	-	-	
Glen reagent 7-deaza-dG-CE Phosphoramidite	10-1021-02	0.25 g	1.396,00	1.605,00	-	-	
Glen reagent 7-deaza-dG-CE Phosphoramidite	10-1021-90	100 µmol	508,00	584,00	-	-	
Glen reagent 7-deaza-dG-CE Phosphoramidite	10-1021-95	50 µmol	254,00	292,00	-	-	
Glen reagent 7-deaza-dX-CE Phosphoramidite	10-1076-02	0.25 g	1.396,00	1.605,00	-	-	
Glen reagent 7-deaza-dX-CE Phosphoramidite	10-1076-90	100 µmol	508,00	584,00	-	-	
Glen reagent 7-deaza-dX-CE Phosphoramidite	10-1076-95	50 µmol	254,00	292,00	-	-	
Glen reagent 7-deaza-guanosine thiotriphosphate (1.0mM) (7-DGTPaS)	80-3321-01	0.1 mL	143,00	164,00	-	-	
Glen reagent 8,5'-Cyclo-dA CE Phosphoramidite	10-1098-95	50 µmol	1.360,00	1.563,00	-	-	
Glen reagent 8-Amino-dA-CE Phosphoramidite	10-1086-02	0.25 g	1.396,00	1.605,00	-	-	
Glen reagent 8-Amino-dA-CE Phosphoramidite	10-1086-90	100 µmol	508,00	584,00	-	-	
Glen reagent 8-Amino-dA-CE Phosphoramidite	10-1086-95	50 µmol	254,00	292,00	-	-	
Glen reagent 8-Amino-dG-CE Phosphoramidite	10-1079-02	0.25 g	1.396,00	1.605,00	-	-	
Glen reagent 8-Amino-dG-CE Phosphoramidite	10-1079-90	100 µmol	508,00	584,00	-	-	
Glen reagent 8-Amino-dG-CE Phosphoramidite	10-1079-95	50 µmol	254,00	292,00	-	-	
Glen reagent 8-Aza-7-deaza-A-CE Phosphoramidite	10-3083-02	0.25 g	2.147,00	2.468,00	-	-	
Glen reagent 8-Aza-7-deaza-A-CE Phosphoramidite	10-3083-90	100 µmol	859,00	987,00	-	-	
Glen reagent 8-Aza-7-deaza-A-CE Phosphoramidite	10-3083-95	50 µmol	429,00	493,00	-	-	
Glen reagent 8-Br-dA-CE Phosphoramidite	10-1007-02	0.25 g	422,00	485,00	-	-	
Glen reagent 8-Br-dA-CE Phosphoramidite	10-1007-90	100 µmol	165,00	190,00	-	-	
Glen reagent 8-Br-dG-CE Phosphoramidite	10-1027-02	0.25 g	365,00	420,00	-	-	
Glen reagent 8-Br-dG-CE Phosphoramidite	10-1027-90	100 µmol	150,00	172,00	-	-	
Glen reagent 8-OMe-dG-CE Phosphoramidite	10-1075-02	0.25 g	1.396,00	1.605,00	-	-	
Glen reagent 8-OMe-dG-CE Phosphoramidite	10-1075-90	100 µmol	508,00	584,00	-	-	
Glen reagent 8-OMe-dG-CE Phosphoramidite	10-1075-95	50 µmol	254,00	292,00	-	-	
Glen reagent 8-oxo-dA-CE Phosphoramidite	10-1008-02	0.25 g	508,00	584,00	-	-	
Glen reagent 8-oxo-dA-CE Phosphoramidite	10-1008-90	100 µmol	193,00	222,00	-	-	
Glen reagent 8-oxo-dG-CE Phosphoramidite	10-1028-02	0.25 g	1.396,00	1.605,00	-	-	
Glen reagent 8-oxo-dG-CE Phosphoramidite	10-1028-90	100 µmol	508,00	584,00	-	-	
Glen reagent 8-oxo-dG-CE Phosphoramidite	10-1028-95	50 µmol	254,00	292,00	-	-	
Glen reagent ABI (240mL capacity)	40-4036-A1	each	7,70	8,85	-	-	

Description	Reference	Scale/Quantity	Price List 2018				Ice Shipping
			EUR	CHF	GBP	USD	
Glen reagent ABI (450mL capacity)	40-4036-A2	each	7,70	8,85	-	-	
Glen reagent Ac-dC-CE Phosphoramidite	10-1015-5S	5 g	83,00	95,40	-	-	
Glen reagent Ac-dC-CPG 1000	20-2015-65	200 x 50 nmol	661,00	760,00	-	-	
Glen reagent Ac-dC-CPG 500	20-2013-10	1 g	53,00	60,90	-	-	
Glen reagent Ac-dC-Me Phosphoramidite	10-1115-02	0.25 g	72,00	82,80	-	-	
Glen reagent Ac-dC-Me Phosphoramidite	10-1115-05	0.5 g	143,00	164,00	-	-	
Glen reagent Ac-dC-Me Phosphoramidite	10-1315-02	0.25 g	35,80	41,10	-	-	
Glen reagent Ac-dC-Me Phosphoramidite	10-1315-05	0.5 g	72,00	82,80	-	-	
Glen reagent Ac-dC-Me Phosphoramidite	10-1315-10	1 g	143,00	164,00	-	-	
Glen reagent Ac-dC-Q-CPG 500	21-2013-01	0.1 g	12,10	13,90	-	-	
Glen reagent Ac-dC-Q-CPG 500	21-2013-02	0.25 g	27,50	31,60	-	-	
Glen reagent Ac-dC-Q-CPG 500	21-2013-10	1 g	105,00	121,00	-	-	
Glen reagent Ac-dC-Q-CPG 500	21-2113-41	4 x 1 µmol	53,00	60,90	-	-	
Glen reagent Ac-dC-Q-CPG 500	21-2113-42	4 x 0.2 µmol	35,20	40,50	-	-	
Glen reagent Ac-dC-Q-CPG 500	21-2213-41	4 x 1 µmol	53,00	60,90	-	-	
Glen reagent Ac-dC-Q-CPG 500	21-2213-42	4 x 0.2 µmol	35,20	40,50	-	-	
Glen reagent Ac-G-CE Phosphoramidite	10-3025-05	0.5 g	53,00	60,90	-	-	
Glen reagent Acridine Phosphoramidite	10-1973-02	0.25 g	966,00	1,110,00	-	-	
Glen reagent Acridine Phosphoramidite	10-1973-90	100 µmol	422,00	485,00	-	-	
Glen reagent Acridine Phosphoramidite	10-1973-95	50 µmol	236,00	271,00	-	-	
Glen reagent Adenosine a-thiotriphosphate (0.5mM) (ATTPaS)	80-3000-01	0.1 mL	143,00	164,00	-	-	
Glen reagent Amino-Modifier C2 dT	10-1037-02	0.25 g	515,00	592,00	-	-	
Glen reagent Amino-Modifier C2 dT	10-1037-05	0.5 g	1,031,00	1,185,00	-	-	
Glen reagent Amino-Modifier C2 dT	10-1037-90	100 µmol	258,00	297,00	-	-	
Glen reagent Amino-Modifier C6 dA	10-1089-02	0.25 g	651,00	748,00	-	-	
Glen reagent Amino-Modifier C6 dA	10-1089-90	100 µmol	293,00	337,00	-	-	
Glen reagent Amino-Modifier C6 dC	10-1019-02	0.25 g	644,00	740,00	-	-	
Glen reagent Amino-Modifier C6 dC	10-1019-90	100 µmol	322,00	370,00	-	-	
Glen reagent Amino-Modifier C6 dT	10-1039-02	0.25 g	476,00	547,00	-	-	
Glen reagent Amino-Modifier C6 dT	10-1039-05	0.5 g	951,00	1,093,00	-	-	
Glen reagent Amino-Modifier C6 dT	10-1039-90	100 µmol	238,00	274,00	-	-	
Glen reagent Amino-Modifier C6 dT, Expedite	10-1039-90E	100 µmol	238,00	274,00	-	-	
Glen reagent Amino-Modifier C6-U Phosphoramidite	10-3039-02	0.25 g	2,111,00	2,426,00	-	-	
Glen reagent Amino-Modifier C6-U Phosphoramidite	10-3039-90	100 µmol	1,031,00	1,185,00	-	-	
Glen reagent Amino-Modifier C6-U Phosphoramidite	10-3039-95	50 µmol	515,00	592,00	-	-	
Glen reagent AP-dC-CE Phosphoramidite	10-1097-02	0.25 g	1,682,00	1,933,00	-	-	
Glen reagent AP-dC-CE Phosphoramidite	10-1097-90	100 µmol	658,00	756,00	-	-	
Glen reagent AP-dC-CE Phosphoramidite	10-1097-95	50 µmol	329,00	378,00	-	-	
Glen reagent Ara-C-CE Phosphoramidite	10-4010-02	0.25 g	508,00	584,00	-	-	
Glen reagent Ara-C-CE Phosphoramidite	10-4010-90	100 µmol	193,00	222,00	-	-	
Glen reagent Asymmetric Doubler Phosphoramidite	10-1921-02	0.25 g	429,00	493,00	-	-	
Glen reagent Asymmetric Doubler Phosphoramidite	10-1921-90	100 µmol	258,00	297,00	-	-	
Glen reagent Beckman Oligo 1000 (30mL capacity)	40-4036-B1	each	7,70	8,85	-	-	
Glen reagent BHQ-1-dT	10-5941-02	0.25 g	1,324,00	1,522,00	-	-	
Glen reagent BHQ-1-dT	10-5941-90	100 µmol	751,00	863,00	-	-	
Glen reagent BHQ-1-dT	10-5941-95	50 µmol	379,00	436,00	-	-	
Glen reagent BHQ-2-dT	10-5942-02	0.25 g	1,324,00	1,522,00	-	-	
Glen reagent BHQ-2-dT	10-5942-90	100 µmol	751,00	863,00	-	-	
Glen reagent BHQ-2-dT	10-5942-95	50 µmol	379,00	436,00	-	-	
Glen reagent Biosearch 8000 (100mL capacity)	40-4036-C1	each	11,60	13,30	-	-	
Glen reagent Biosearch 8000 (500mL capacity)	40-4036-C2	each	11,60	13,30	-	-	
Glen reagent Biotin Phosphoramidite	10-1953-02	0.25 g	966,00	1,110,00	-	-	
Glen reagent Biotin Phosphoramidite	10-1953-90	100 µmol	422,00	485,00	-	-	
Glen reagent Biotin Phosphoramidite	10-1953-95	50 µmol	236,00	271,00	-	-	
Glen reagent Biotin-dT	10-1038-02	0.25 g	892,00	1,025,00	-	-	
Glen reagent Biotin-dT	10-1038-90	100 µmol	429,00	493,00	-	-	
Glen reagent Biotin-dT	10-1038-95	50 µmol	238,00	274,00	-	-	
Glen reagent BiotinTEG Phosphoramidite	10-1955-02	0.25 g	526,00	605,00	461,00	631,00	
Glen reagent BiotinTEG Phosphoramidite	10-1955-90	100 µmol	230,00	264,00	202,00	276,00	
Glen reagent BiotinTEG Phosphoramidite	10-1955-95	50 µmol	128,00	147,00	112,00	154,00	
Glen reagent Carboxy-dT	10-1035-02	0.25 g	515,00	592,00	-	-	
Glen reagent Carboxy-dT	10-1035-90	100 µmol	258,00	297,00	-	-	
Glen reagent Chemical Phosphorylation Reagent	10-1900-02	0.25 g	176,00	202,00	-	-	
Glen reagent Chemical Phosphorylation Reagent	10-1900-90	100 µmol	49,50	56,90	-	-	
Glen reagent Chemical Phosphorylation Reagent II	10-1901-02	0.25 g	220,00	253,00	-	-	
Glen reagent Chemical Phosphorylation Reagent II	10-1901-90	100 µmol	66,00	75,90	-	-	
Glen reagent Cholesteryl-TEG Phosphoramidite	10-1975-02	0.25 g	780,00	897,00	-	-	
Glen reagent Cholesteryl-TEG Phosphoramidite	10-1975-90	100 µmol	379,00	-	-	-	
Glen reagent Cholesteryl-TEG Phosphoramidite	10-1975-95	50 µmol	200,00	-	-	-	
Glen reagent Cis-syn Thymine Dimer Phosphoramidite	11-1330-02	0.25 g	14,599,00	16,780,00	-	-	
Glen reagent Cis-syn Thymine Dimer Phosphoramidite	11-1330-90	100 µmol	6,011,00	6,909,00	-	-	
Glen reagent Cis-syn Thymine Dimer Phosphoramidite	11-1330-95	50 µmol	3,006,00	3,455,00	-	-	
Glen reagent Cy3.5* Phosphoramidite	10-5914-02	0.25 g	2,389,00	2,746,00	-	-	
Glen reagent Cy3.5* Phosphoramidite	10-5914-90	100 µmol	1,017,00	1,169,00	-	-	
Glen reagent Cy3.5* Phosphoramidite	10-5914-95	50 µmol	532,00	611,00	-	-	
Glen reagent Cy3* Phosphoramidite	10-5913-02	0.25 g	1,252,00	1,439,00	-	-	
Glen reagent Cy3* Phosphoramidite	10-5913-90	100 µmol	501,00	576,00	-	-	
Glen reagent Cy3* Phosphoramidite	10-5913-95	50 µmol	272,00	313,00	-	-	
Glen reagent Cy3* Phosphoramidite, Cycl.	10-5913-95C	50 µmol	241,00	277,00	211,00	-	
Glen reagent Cy5.5* Phosphoramidite	10-5916-02	0.25 g	1,349,00	1,551,00	-	-	
Glen reagent Cy5.5* Phosphoramidite	10-5916-90	100 µmol	647,00	744,00	-	-	
Glen reagent Cy5.5* Phosphoramidite	10-5916-95	50 µmol	347,00	399,00	-	-	
Glen reagent Cy5* Phosphoramidite	10-5915-02	0.25 g	1,252,00	1,439,00	-	-	
Glen reagent Cy5* Phosphoramidite	10-5915-90	100 µmol	501,00	576,00	-	-	
Glen reagent Cy5* Phosphoramidite	10-5915-95	50 µmol	272,00	313,00	-	-	
Glen reagent Cy5* Phosphoramidite, Pharmacia	10-5915-95P	50 µmol	272,00	313,00	-	-	
Glen reagent Cytidine a-thiotriphosphate (1.0mM) (CTTPaS)	80-3010-01	0.1 mL	143,00	164,00	-	-	
Glen reagent dA PS	26-2600-61	200 x 1 µmol	1,244,00	1,430,00	-	-	
Glen reagent dA+dC+dG+dT-CE Phosphoramidite Mix	10-1023-02	0.25 g	57,00	65,50	-	-	
Glen reagent dA+dG-CE Phosphoramidite	10-1002-02	0.25 g	57,00	65,50	-	-	
Glen reagent dA-5'-CE Phosphoramidite	10-0001-02	0.25 g	91,00	104,60	-	-	
Glen reagent dA-5'-CE Phosphoramidite	10-0001-05	0.5 g	182,00	209,00	-	-	
Glen reagent dA-5'-CE Phosphoramidite	10-0001-10	1 g	363,00	417,00	-	-	

Description	Reference	Scale/Quantity	Price List 2018				Ice Shipping
			EUR	CHF	GBP	USD	
Glen reagent dA-5'-CE Phosphoramidite, Expedite	10-0001-05E	0.5 g	182,00	209,00	-	-	
Glen reagent dA-5'-CPG	20-0002-01	0.1 g	72,00	82,80	-	-	
Glen reagent dA-5'-CPG	20-0002-10	1 g	537,00	617,00	-	-	
Glen reagent dA-5'-CPG	20-0012-13	1 x 10 µmol	322,00	370,00	-	-	
Glen reagent dA-5'-CPG	20-0012-14	1 x 15 µmol	429,00	493,00	-	-	
Glen reagent dA-5'-CPG	20-0012-41	4 x 1 µmol	143,00	164,00	-	-	
Glen reagent dA-5'-CPG	20-0012-42	4 x 0.2 µmol	107,00	123,00	-	-	
Glen reagent Dabacyl-dT	10-1058-02	0.25 g	966,00	1.110,00	-	-	
Glen reagent Dabacyl-dT	10-1058-90	100 µmol	465,00	534,00	-	-	
Glen reagent Dabacyl-dT	10-1058-95	50 µmol	258,00	297,00	-	-	
Glen reagent dA-CPG 1000	20-2001-65	200 x 50 µmol	661,00	760,00	-	-	
Glen reagent dA-H-Phosphonate, TEA Salt	10-1200-02	0.25 g	57,00	65,50	-	-	
Glen reagent dA-H-Phosphonate, TEA Salt	10-1200-05	0.5 g	115,00	132,00	-	-	
Glen reagent dA-Me Phosphoramidite	10-1100-02	0.25 g	72,00	82,80	-	-	
Glen reagent dA-Me Phosphoramidite	10-1100-05	0.5 g	143,00	164,00	-	-	
Glen reagent dA-Q-CPG 500	21-2000-01	0.1 g	12,10	13,90	-	-	
Glen reagent dA-Q-CPG 500	21-2000-02	0.25 g	27,50	31,60	-	-	
Glen reagent dA-Q-CPG 500	21-2000-10	1 g	105,00	121,00	-	-	
Glen reagent dA-Q-CPG 500	21-2010-01	0.1 g	12,10	13,90	-	-	
Glen reagent dA-Q-CPG 500	21-2010-02	0.25 g	27,50	31,60	-	-	
Glen reagent dA-Q-CPG 500	21-2010-10	1 g	105,00	121,00	-	-	
Glen reagent dA-Q-CPG 500	21-2100-41	4 x 1 µmol	53,00	60,90	-	-	
Glen reagent dA-Q-CPG 500	21-2100-42	4 x 0.2 µmol	35,20	40,50	-	-	
Glen reagent dA-Q-CPG 500	21-2200-41	4 x 1 µmol	53,00	60,90	-	-	
Glen reagent dA-Q-CPG 500	21-2200-42	4 x 0.2 µmol	35,20	40,50	-	-	
Glen reagent dC PS	26-2610-61	200 x 1 µmol	1.244,00	1.430,00	-	-	
Glen reagent dC+dT-CE Phosphoramidite	10-1013-02	0.25 g	57,00	65,50	-	-	
Glen reagent dC-5'-CE Phosphoramidite	10-0101-02	0.25 g	91,00	104,60	-	-	
Glen reagent dC-5'-CE Phosphoramidite	10-0101-05	0.5 g	182,00	209,00	-	-	
Glen reagent dC-5'-CE Phosphoramidite	10-0101-10	1 g	363,00	417,00	-	-	
Glen reagent dC-5'-CE Phosphoramidite, Expedite	10-0101-02E	0.25 g	93,00	106,90	-	-	
Glen reagent dC-5'-CE Phosphoramidite, Expedite	10-0101-05E	0.5 g	182,00	209,00	-	-	
Glen reagent dC-5'-CPG	20-0102-01	0.1 g	72,00	82,80	-	-	
Glen reagent dC-5'-CPG	20-0102-10	1 g	537,00	617,00	-	-	
Glen reagent dC-5'-CPG	20-0112-13	1 x 10 µmol	322,00	370,00	-	-	
Glen reagent dC-5'-CPG	20-0112-14	1 x 15 µmol	429,00	493,00	-	-	
Glen reagent dC-5'-CPG	20-0112-41	4 x 1 µmol	143,00	164,00	-	-	
Glen reagent dC-5'-CPG	20-0112-42	4 x 0.2 µmol	107,00	123,00	-	-	
Glen reagent dC-5'-CPG	20-0202-01	0.1 g	72,00	82,80	-	-	
Glen reagent dC-H-Phosphonate, DBU Salt	10-1210-02	0.25 g	57,00	65,50	-	-	
Glen reagent dC-H-Phosphonate, DBU Salt	10-1210-05	0.5 g	115,00	132,00	-	-	
Glen reagent dC-Q-CPG 500	21-2110-41	4 x 1 µmol	53,00	60,90	-	-	
Glen reagent dC-Q-CPG 500	21-2110-42	4 x 0.2 µmol	35,20	40,50	-	-	
Glen reagent dC-Q-CPG 500	21-2210-41	4 x 1 µmol	53,00	60,90	-	-	
Glen reagent dC-Q-CPG 500	21-2210-42	4 x 0.2 µmol	35,20	40,50	-	-	
Glen reagent DesthiobiotinTEG Phosphoramidite	10-1952-02	0.25 g	1.109,00	1.275,00	-	-	
Glen reagent DesthiobiotinTEG Phosphoramidite	10-1952-90	100 µmol	480,00	552,00	-	-	
Glen reagent DesthiobiotinTEG Phosphoramidite	10-1952-95	50 µmol	265,00	305,00	-	-	
Glen reagent DesthiobiotinTEG-CPG	20-2952-01	0.1 g	200,00	230,00	-	-	
Glen reagent DesthiobiotinTEG-CPG	20-2952-10	1 g	1.646,00	1.892,00	-	-	
Glen reagent DesthiobiotinTEG-CPG	20-2952-13	1 x 10 µmol	494,00	568,00	-	-	
Glen reagent DesthiobiotinTEG-CPG	20-2952-14	1 x 15 µmol	744,00	855,00	-	-	
Glen reagent DesthiobiotinTEG-CPG	20-2952-41	4 x 1 µmol	329,00	378,00	-	-	
Glen reagent DesthiobiotinTEG-CPG	20-2952-42	4 x 0.2 µmol	200,00	230,00	-	-	
Glen reagent dG-5'-CPG	20-0202-10	1 g	537,00	617,00	-	-	
Glen reagent dG-5'-CPG	20-0212-13	1 x 10 µmol	322,00	370,00	-	-	
Glen reagent dG-5'-CPG	20-0212-14	1 x 15 µmol	429,00	493,00	-	-	
Glen reagent dG-5'-CPG	20-0212-41	4 x 1 µmol	143,00	164,00	-	-	
Glen reagent dG-5'-CPG	20-0212-42	4 x 0.2 µmol	107,00	123,00	-	-	
Glen reagent dG-CPG 1000	20-2021-65	200 x 50 µmol	661,00	760,00	-	-	
Glen reagent dG-H-Phosphonate, TEA Salt	10-1220-02	0.25 g	57,00	65,50	-	-	
Glen reagent dG-H-Phosphonate, TEA Salt	10-1220-05	0.5 g	115,00	132,00	-	-	
Glen reagent dG-Me Phosphoramidite	10-1120-02	0.25 g	72,00	82,80	-	-	
Glen reagent dG-Me Phosphoramidite	10-1120-05	0.5 g	143,00	164,00	-	-	
Glen reagent dI-CE Phosphoramidite	10-1040-02	0.25 g	132,00	152,00	-	-	
Glen reagent dI-CE Phosphoramidite	10-1040-90	100 µmol	55,00	63,20	-	-	
Glen reagent dI-CPG 1000	20-2041-01	0.1 g	42,90	49,30	-	-	
Glen reagent dI-CPG 1000	20-2191-41	4 x 1 µmol	172,00	198,00	-	-	
Glen reagent dI-CPG 1000	20-2191-42	4 x 0.2 µmol	103,00	118,00	-	-	
Glen reagent dI-CPG 500	20-2040-01	0.1 g	42,90	49,30	-	-	
Glen reagent dI-CPG 500	20-2190-41	4 x 1 µmol	172,00	198,00	-	-	
Glen reagent dI-CPG 500	20-2190-42	4 x 0.2 µmol	103,00	118,00	-	-	
Glen reagent Dithiol Phosphoramidite (DTPA)	10-1937-02	0.25 g	687,00	790,00	-	-	
Glen reagent Dithiol Phosphoramidite (DTPA)	10-1937-90	100 µmol	286,00	329,00	-	-	
Glen reagent dK-CE Phosphoramidite	10-1048-02	0.25 g	852,00	979,00	-	-	
Glen reagent dK-CE Phosphoramidite	10-1048-90	100 µmol	279,00	321,00	-	-	
Glen reagent dmf-5-Me-isodC-CE Phosphoramidite	10-1065-02	0.25 g	309,00	355,00	-	-	
Glen reagent dmf-5-Me-isodC-CE Phosphoramidite	10-1065-90	100 µmol	121,00	139,00	-	-	
Glen reagent dmf-5-Me-isodC-CE Phosphoramidite, Expedite	10-1065-02E	0.25 g	309,00	355,00	-	-	
Glen reagent dmf-dG PS	26-2629-61	200 x 1 µmol	1.244,00	1.430,00	-	-	
Glen reagent dmf-dG PS	26-2629-62	200 x 200 nmol	650,00	747,00	-	-	
Glen reagent dmf-dG PS	26-2629-65	200 x 40 nmol	650,00	747,00	-	-	
Glen reagent dmf-dG-CE Phosphoramidite	10-1029-02	0.25 g	3,90	4,48	-	-	
Glen reagent dmf-dG-CE Phosphoramidite	10-1029-05	0.5 g	7,70	8,85	-	-	
Glen reagent dmf-dG-CE Phosphoramidite	10-1029-10	1 g	15,40	17,70	-	-	
Glen reagent dmf-dG-CE Phosphoramidite	10-1029-1B	1 g	15,40	17,70	-	-	
Glen reagent dmf-dG-CE Phosphoramidite	10-1029-1C	1 g	15,40	17,70	-	-	
Glen reagent dmf-dG-CE Phosphoramidite	10-1029-1P	1 g	15,40	17,70	-	-	
Glen reagent dmf-dG-CE Phosphoramidite	10-1029-1S	10 g	154,00	177,00	-	-	
Glen reagent dmf-dG-CE Phosphoramidite	10-1029-20	2 g	30,80	35,40	-	-	
Glen reagent dmf-dG-CE Phosphoramidite	10-1029-50	5 g	77,00	88,50	-	-	
Glen reagent dmf-dG-CE Phosphoramidite	10-1029-5S	5 g	77,00	88,50	-	-	
Glen reagent dmf-dG-CE Phosphoramidite	10-1029-B2	0.25 g	3,90	4,48	-	-	

Description	Reference	Scale/Quantity	Price List 2018				Ice Shipping
			EUR	CHF	GBP	USD	
Glen reagent dmf-dG-CE Phosphoramidite	10-1029-B5	0.5 g	7,70	8,85	-	-	
Glen reagent dmf-dG-CE Phosphoramidite	10-1029-C2	0.25 g	3,90	4,48	-	-	
Glen reagent dmf-dG-CE Phosphoramidite	10-1029-C5	0.5 g	7,70	8,85	-	-	
Glen reagent dmf-dG-CE Phosphoramidite	10-1029-P4	0.4 g	6,20	7,13	-	-	
Glen reagent dmf-dG-CPG	20-2029-01	0.1 g	6,30	7,24	-	-	
Glen reagent dmf-dG-CPG	20-2029-02	0.25 g	14,10	16,20	-	-	
Glen reagent dmf-dG-CPG	20-2029-10	1 g	53,00	60,90	-	-	
Glen reagent dmf-dG-CPG	20-2129-13	1 x 10 µmol	88,00	101,10	-	-	
Glen reagent dmf-dG-CPG	20-2129-41	4 x 1 µmol	30,80	35,40	-	-	
Glen reagent dmf-dG-CPG	20-2129-42	4 x 0.2 µmol	30,80	35,40	-	-	
Glen reagent dmf-dG-CPG	20-2129-45	4 x 40 nmol	30,80	35,40	-	-	
Glen reagent dmf-dG-CPG	20-2229-14	1 x 15 µmol	132,00	152,00	-	-	
Glen reagent dmf-dG-CPG	20-2229-41	4 x 1 µmol	22,00	25,30	-	-	
Glen reagent dmf-dG-CPG	20-2229-42	4 x 0.2 µmol	15,40	17,70	-	-	
Glen reagent dmf-dG-CPG	20-2229-45	4 x 40 nmol	15,40	17,70	-	-	
Glen reagent dmf-dG-Q-CPG 500	21-2029-01	0.1 g	12,10	13,90	-	-	
Glen reagent dmf-dG-Q-CPG 500	21-2029-02	0.25 g	27,50	31,60	-	-	
Glen reagent dmf-dG-Q-CPG 500	21-2029-10	1 g	105,00	121,00	-	-	
Glen reagent dmf-dG-Q-CPG 500	21-2129-41	4 x 1 µmol	53,00	60,90	-	-	
Glen reagent dmf-dG-Q-CPG 500	21-2129-42	4 x 0.2 µmol	35,20	40,50	-	-	
Glen reagent dmf-dG-Q-CPG 500	21-2229-41	4 x 1 µmol	53,00	60,90	-	-	
Glen reagent dmf-dG-Q-CPG 500	21-2229-42	4 x 0.2 µmol	35,20	40,50	-	-	
Glen reagent dmf-isodG-CE Phosphoramidite	10-1078-02	0.25 g	430,00	494,00	-	-	
Glen reagent dmf-isodG-CE Phosphoramidite	10-1078-90	100 µmol	200,00	230,00	-	-	
Glen reagent DNP-TEG Phosphoramidite	10-1985-02	0.25 g	966,00	1,110,00	-	-	
Glen reagent DNP-TEG Phosphoramidite	10-1985-90	100 µmol	422,00	485,00	-	-	
Glen reagent DNP-TEG Phosphoramidite	10-1985-95	50 µmol	236,00	271,00	-	-	
Glen reagent dP+dK-CE Phosphoramidite	10-1049-02	0.25 g	852,00	979,00	-	-	
Glen reagent dP+dK-CE Phosphoramidite	10-1049-90	100 µmol	279,00	321,00	-	-	
Glen reagent dP-CE Phosphoramidite	10-1047-02	0.25 g	852,00	979,00	-	-	
Glen reagent dP-CE Phosphoramidite	10-1047-90	100 µmol	279,00	321,00	-	-	
Glen reagent dSpacer-CE Phosphoramidite	10-1914-02	0.25 g	328,00	377,00	288,00	394,00	
Glen reagent dSpacer-CE Phosphoramidite	10-1914-90	100 µmol	95,00	109,20	83,30	114,00	
Glen reagent dT-5'-CE Phosphoramidite	10-0301-02	0.25 g	91,00	104,60	-	-	
Glen reagent dT-5'-CE Phosphoramidite	10-0301-05	0.5 g	182,00	209,00	-	-	
Glen reagent dT-5'-CE Phosphoramidite	10-0301-10	1 g	363,00	417,00	-	-	
Glen reagent dT-5'-CE Phosphoramidite, Expedite	10-0301-10E	1 g	1,244,00	1,430,00	-	-	
Glen reagent dT-5'-CPG	20-0302-01	0.1 g	72,00	82,80	-	-	
Glen reagent dT-5'-CPG	20-0302-10	1 g	537,00	617,00	-	-	
Glen reagent dT-5'-CPG	20-0312-13	1 x 10 µmol	322,00	370,00	-	-	
Glen reagent dT-5'-CPG	20-0312-14	1 x 15 µmol	429,00	493,00	-	-	
Glen reagent dT-5'-CPG	20-0312-41	4 x 1 µmol	143,00	164,00	-	-	
Glen reagent dT-5'-CPG	20-0312-42	4 x 0.2 µmol	107,00	123,00	-	-	
Glen reagent dT-5'-Icaa-CPG 500	20-0312-41E	1 µmol	150,00	172,00	-	-	
Glen reagent dT-CPG 1000	20-2031-65	200 x 50 µmol	661,00	760,00	-	-	
Glen reagent dT-H-Phosphonate, TEA Salt	10-1230-02	0.25 g	57,00	65,50	-	-	
Glen reagent dT-H-Phosphonate, TEA Salt	10-1230-05	0.5 g	115,00	132,00	-	-	
Glen reagent dT-Me Phosphonamidite	10-1130-02	0.25 g	72,00	82,80	-	-	
Glen reagent dT-Me Phosphonamidite	10-1130-05	0.5 g	143,00	164,00	-	-	
Glen reagent dT-Me Phosphoramidite	10-1330-02	0.25 g	35,80	41,10	-	-	
Glen reagent dT-Me Phosphoramidite	10-1330-05	0.5 g	72,00	82,80	-	-	
Glen reagent dT-Me Phosphoramidite	10-1330-10	1 g	143,00	164,00	-	-	
Glen reagent dT-PS	26-2630-61	200 x 1 µmol	1,244,00	1,430,00	-	-	
Glen reagent dT-Q-CPG 500	21-2030-01	0.1 g	12,10	13,90	-	-	
Glen reagent dT-Q-CPG 500	21-2030-02	0.25 g	27,50	31,60	-	-	
Glen reagent dT-Q-CPG 500	21-2030-10	1 g	105,00	121,00	-	-	
Glen reagent dT-Q-CPG 500	21-2130-41	4 x 1 µmol	53,00	60,90	-	-	
Glen reagent dT-Q-CPG 500	21-2130-42	4 x 0.2 µmol	35,20	40,50	-	-	
Glen reagent dT-Q-CPG 500	21-2230-41	4 x 1 µmol	53,00	60,90	-	-	
Glen reagent dU-CE Phosphoramidite	10-1050-02	0.25 g	110,00	126,00	-	-	
Glen reagent dU-CE Phosphoramidite	10-1050-90	100 µmol	38,50	44,30	-	-	
Glen reagent dU-CPG 1000	20-2051-01	0.1 g	72,00	82,80	-	-	
Glen reagent dU-CPG 1000	20-2151-41	4 x 1 µmol	286,00	329,00	-	-	
Glen reagent dU-CPG 1000	20-2151-42	4 x 0.2 µmol	172,00	198,00	-	-	
Glen reagent dU-CPG 500	20-2050-01	0.1 g	42,90	49,30	-	-	
Glen reagent dU-CPG 500	20-2150-41	4 x 1 µmol	172,00	198,00	-	-	
Glen reagent dU-CPG 500	20-2150-42	4 x 0.2 µmol	103,00	118,00	-	-	
Glen reagent EDTA-C2-dT-CE Phosphoramidite	10-1059-02	0.25 g	1,396,00	1,605,00	-	-	
Glen reagent EDTA-C2-dT-CE Phosphoramidite	10-1059-90	100 µmol	709,00	815,00	-	-	
Glen reagent EDTA-C2-dT-CE Phosphoramidite	10-1059-95	50 µmol	358,00	411,00	-	-	
Glen reagent Epoch Eclipse™ Quencher CPG	20-5925-01	0.1 g	329,00	378,00	-	-	
Glen reagent Epoch Eclipse™ Quencher CPG	20-5925-10	1 g	2,755,00	3,167,00	-	-	
Glen reagent Epoch Eclipse™ Quencher CPG	20-5925-13	1 x 10 µmol	1,424,00	1,637,00	-	-	
Glen reagent Epoch Eclipse™ Quencher CPG	20-5925-14	1 x 15 µmol	2,140,00	2,460,00	-	-	
Glen reagent Epoch Eclipse™ Quencher CPG	20-5925-41	4 x 1 µmol	501,00	576,00	-	-	
Glen reagent Epoch Eclipse™ Quencher CPG	20-5925-42	4 x 0.2 µmol	251,00	289,00	-	-	
Glen reagent Epoch Eclipse™ Quencher Phosphoramidite	10-5925-02	0.25 g	1,465,00	1,684,00	-	-	
Glen reagent Epoch Eclipse™ Quencher Phosphoramidite	10-5925-90	100 µmol	593,00	682,00	-	-	
Glen reagent Epoch Eclipse™ Quencher Phosphoramidite	10-5925-95	50 µmol	309,00	355,00	-	-	
Glen reagent Epoch Gig Harbor Green™ Phosphoramidite	10-5922-02	0.25 g	1,138,00	1,308,00	-	-	
Glen reagent Epoch Gig Harbor Green™ Phosphoramidite	10-5922-90	100 µmol	427,00	491,00	-	-	
Glen reagent Epoch Gig Harbor Green™ Phosphoramidite	10-5922-95	50 µmol	226,00	260,00	-	-	
Glen reagent Epoch Redmond Red™ CPG	20-5920-01	0.1 g	258,00	297,00	-	-	
Glen reagent Epoch Redmond Red™ CPG	20-5920-10	1 g	2,147,00	2,468,00	-	-	
Glen reagent Epoch Redmond Red™ CPG	20-5920-13	1 x 10 µmol	1,073,00	1,233,00	-	-	
Glen reagent Epoch Redmond Red™ CPG	20-5920-14	1 x 15 µmol	1,610,00	1,851,00	-	-	
Glen reagent Epoch Redmond Red™ CPG	20-5920-41	4 x 1 µmol	429,00	493,00	-	-	
Glen reagent Epoch Redmond Red™ CPG	20-5920-42	4 x 0.2 µmol	215,00	247,00	-	-	
Glen reagent Epoch Redmond Red™ Phosphoramidite	10-5920-02	0.25 g	1,292,00	1,485,00	-	-	
Glen reagent Epoch Redmond Red™ Phosphoramidite	10-5920-90	100 µmol	519,00	597,00	-	-	
Glen reagent Epoch Redmond Red™ Phosphoramidite	10-5920-95	50 µmol	272,00	313,00	-	-	
Glen reagent Epoch Yakima Yellow™ CPG	20-5921-01	0.1 g	258,00	297,00	-	-	
Glen reagent Epoch Yakima Yellow™ CPG	20-5921-10	1 g	2,147,00	2,468,00	-	-	

Description	Reference	Scale/Quantity	Price List 2018				Ice Shipping
			EUR	CHF	GBP	USD	
Glen reagent Epoch Yakima Yellow™ CPG	20-5921-13	1 x 10 µmol	1,073,00	1,233,00	-	-	
Glen reagent Epoch Yakima Yellow™ CPG	20-5921-14	1 x 15 µmol	1,610,00	1,851,00	-	-	
Glen reagent Epoch Yakima Yellow™ CPG	20-5921-41	4 x 1 µmol	429,00	493,00	-	-	
Glen reagent Epoch Yakima Yellow™ CPG	20-5921-42	4 x 0.2 µmol	215,00	247,00	-	-	
Glen reagent Epoch Yakima Yellow™ Phosphoramidite	10-5921-02	0.25 g	1,292,00	1,485,00	-	-	
Glen reagent Epoch Yakima Yellow™ Phosphoramidite	10-5921-90	100 µmol	544,00	625,00	-	-	
Glen reagent Epoch Yakima Yellow™ Phosphoramidite	10-5921-95	50 µmol	284,00	326,00	-	-	
Glen reagent Etheno-dA-CE Phosphoramidite	10-1006-02	0.25 g	365,00	420,00	-	-	
Glen reagent Etheno-dA-CE Phosphoramidite	10-1006-90	100 µmol	150,00	172,00	-	-	
Glen reagent Expedite 8905 (150mL capacity)	40-4036-C3	each	7,70	8,85	-	-	
Glen reagent Expedite 8909 and Cyclone (240mL capacity)	40-4036-C4	each	7,70	8,85	-	-	
Glen reagent Fluorescein Phosphoramidite	10-1963-02	0.25 g	852,00	979,00	-	-	
Glen reagent Fluorescein Phosphoramidite	10-1963-90	10 µmol	422,00	485,00	-	-	
Glen reagent Fluorescein Phosphoramidite	10-1963-95	50 µmol	236,00	271,00	-	-	
Glen reagent Fluorescein-dT Phosphoramidite	10-1056-02	0.25 g	751,00	863,00	659,00	902,00	
Glen reagent Fluorescein-dT Phosphoramidite	10-1056-90	100 µmol	362,00	416,00	318,00	435,00	
Glen reagent Fluorescein-dT Phosphoramidite	10-1056-95	50 µmol	200,00	230,00	175,00	240,00	
Glen reagent Fluorescein-dUTP-dabyl (1mM, 25nmoles, 10mM Tris, 1mM EDTA)	88-1056-01	0.1 mL	386,00	444,00	-	-	
Glen reagent Fmoc Amino-Modifier C6 dT	10-1536-02	0.25 g	515,00	592,00	-	-	
Glen reagent Fmoc Amino-Modifier C6 dT	10-1536-90	100 µmol	258,00	297,00	-	-	
Glen reagent Glen-Pak™ DNA purification cartridge (for use with disposable syringes), Pk/1	60-5200-01		9,70	11,15	-	-	
Glen reagent Glen-Pak™ DNA purification cartridge (for use with disposable syringes), Pk/10	60-5200-10		79,00	90,80	-	-	
Glen reagent Glen-Pak™ DNA purification cartridge, Pk/10	60-5100-10		88,00	101,10	-	-	
Glen reagent Glen-Pak™ DNA purification cartridge, Pk/30	60-5100-30		176,00	202,00	-	-	
Glen reagent Glen-Pak™ DNA purification cartridge, Pk/96	60-5100-96		317,00	364,00	-	-	
Glen reagent Glen-Pak™ RNA purification cartridge (for use with disposable syringes), Pk/1	60-6200-01		11,50	13,20	-	-	
Glen reagent Glen-Pak™ RNA purification cartridge (for use with disposable syringes), Pk/10	60-6200-10		94,00	108,00	-	-	
Glen reagent Glen-Pak™ RNA purification cartridge, Pk/10	60-6100-10		146,00	168,00	-	-	
Glen reagent Glen-Pak™ RNA purification cartridge, Pk/30	60-6100-30		297,00	341,00	-	-	
Glen reagent Glen-Pak™ RNA purification cartridge, Pk/96	60-6100-96		633,00	728,00	-	-	
Glen reagent Guanosine a-thiotriphosphate (0.5mM) (GTPaS)	80-3020-01	0.1 mL	143,00	164,00	-	-	
Glen reagent I-CE Phosphoramidite	10-3040-02	0.25 g	680,00	782,00	-	-	
Glen reagent Inosine a-thiotriphosphate (4mM) (ITTPaS)	80-3050-01	0.1 mL	143,00	164,00	-	-	
Glen reagent iPr-Pac-dG-CE Phosphoramidite	10-1621-02	0.25 g	21,50	24,70	-	-	
Glen reagent iPr-Pac-dG-CE Phosphoramidite	10-1621-05	0.5 g	42,90	49,30	-	-	
Glen reagent iPr-Pac-dG-CE Phosphoramidite	10-1621-10	1 g	86,00	98,90	-	-	
Glen reagent iPr-Pac-dG-CE Phosphoramidite, Expedite	10-1621-05E	0.5 g	42,90	-	-	-	
Glen reagent iPr-Pac-dG-CPG	20-2621-01	0.1 g	25,80	29,70	-	-	
Glen reagent iPr-Pac-dG-CPG	20-2621-02	0.25 g	57,00	65,50	-	-	
Glen reagent iPr-Pac-dG-CPG	20-2621-10	1 g	215,00	247,00	-	-	
Glen reagent iPr-Pac-dG-CPG	20-2721-13	1 x 10 µmol	143,00	164,00	-	-	
Glen reagent iPr-Pac-dG-CPG	20-2721-41	4 x 1 µmol	86,00	98,90	-	-	
Glen reagent iPr-Pac-dG-CPG	20-2721-42	4 x 0.2 µmol	57,00	65,50	-	-	
Glen reagent iPr-Pac-dG-CPG	20-2721-45	4 x 40 nmol	57,00	65,50	-	-	
Glen reagent iPr-Pac-dG-CPG	20-2821-14	1 x 15 µmol	215,00	247,00	-	-	
Glen reagent iPr-Pac-dG-CPG	20-2821-41	4 x 1 µmol	86,00	98,90	-	-	
Glen reagent iPr-Pac-dG-CPG	20-2821-42	4 x 0.2 µmol	57,00	65,50	-	-	
Glen reagent iPr-Pac-dG-CPG	20-2821-45	4 x 40 nmol	57,00	65,50	-	-	
Glen reagent iPr-Pac-dG-Me Phosphoramidite	10-1321-02	0.25 g	35,80	41,10	-	-	
Glen reagent iPr-Pac-dG-Me Phosphoramidite	10-1321-05	0.5 g	72,00	82,80	-	-	
Glen reagent iPr-Pac-dG-Me Phosphoramidite	10-1321-10	1 g	143,00	164,00	-	-	
Glen reagent iPr-Pac-G-RNA-CPG	20-3321-10	1 g	508,00	584,00	-	-	
Glen reagent Long Trebler Phosphoramidite	10-1925-02	0.25 g	429,00	493,00	-	-	
Glen reagent Long Trebler Phosphoramidite	10-1925-90	100 µmol	286,00	329,00	-	-	
Glen reagent N3-Cyanoethyl-dT	10-1531-02	0.25 g	859,00	987,00	-	-	
Glen reagent N3-Cyanoethyl-dT	10-1531-90	100 µmol	286,00	329,00	-	-	
Glen reagent N4-Et-dC-CE Phosphoramidite	10-1068-02	0.25 g	966,00	1,110,00	-	-	
Glen reagent N4-Et-dC-CE Phosphoramidite	10-1068-90	100 µmol	322,00	370,00	-	-	
Glen reagent N4-Et-dC-CE Phosphoramidite	10-1068-95	50 µmol	179,00	206,00	-	-	
Glen reagent N4-Et-dC-CE Phosphoramidite	10-1068-SP	1 g	1,670,00	-	-	-	
Glen reagent N6-Me-Adenosine a-thiotriphosphate (4mM) (N6-Me-ATTPaS)	80-3302-01	0.1 mL	143,00	164,00	-	-	
Glen reagent N6-Me-dA-CE Phosphoramidite	10-1003-02	0.25 g	709,00	815,00	-	-	
Glen reagent N6-Me-dA-CE Phosphoramidite	10-1003-90	100 µmol	233,00	268,00	-	-	
Glen reagent Nebularine-CE Phosphoramidite	10-3041-02	0.25 g	1,789,00	2,056,00	-	-	
Glen reagent Nebularine-CE Phosphoramidite	10-3041-90	100 µmol	644,00	740,00	-	-	
Glen reagent Nebularine-CE Phosphoramidite	10-3041-95	50 µmol	322,00	370,00	-	-	
Glen reagent NHS-Carboxy-dT	10-1535-02	0.25 g	787,00	905,00	-	-	
Glen reagent NHS-Carboxy-dT	10-1535-90	100 µmol	301,00	346,00	-	-	
Glen reagent O4-Me-dT-CE Phosphoramidite	10-1032-02	0.25 g	508,00	584,00	-	-	
Glen reagent O4-Me-dT-CE Phosphoramidite	10-1032-90	100 µmol	193,00	222,00	-	-	
Glen reagent O4-Triazolyl-dU-CE Phosphoramidite	10-1051-02	0.25 g	508,00	584,00	-	-	
Glen reagent O4-Triazolyl-dU-CE Phosphoramidite	10-1051-90	100 µmol	193,00	222,00	-	-	
Glen reagent O6-Me-dG-CE Phosphoramidite	10-1070-02	0.25 g	365,00	420,00	-	-	
Glen reagent O6-Me-dG-CE Phosphoramidite	10-1070-90	100 µmol	150,00	172,00	-	-	
Glen reagent O6-Phenyl-dl-CE Phosphoramidite	10-1042-02	0.25 g	508,00	584,00	-	-	
Glen reagent O6-Phenyl-dl-CE Phosphoramidite	10-1042-90	100 µmol	193,00	222,00	-	-	
Glen reagent Oligo-Affinity Support (PS)	26-4001-01	0.1 g	258,00	297,00	-	-	
Glen reagent Oligo-Affinity Support (PS)	26-4001-02	0.25 g	608,00	699,00	-	-	
Glen reagent Oligo-Affinity Support (PS)	26-4001-10	1 g	2,276,00	2,616,00	-	-	
Glen reagent Oligo-Affinity Support (PS)	26-4101-41	4 x 1 µmol	429,00	493,00	-	-	
Glen reagent Oligo-Affinity Support (PS), Expedite	26-4101-41E	4 x 1 µmol	429,00	493,00	-	-	
Glen reagent Oligonucleotide-Peptide Conjugation Kit (OPeC)	50-1000-01	1 g	709,00	815,00	-	-	
Glen reagent Pac-A-RNA-CPG	20-3300-10	1 g	508,00	584,00	-	-	
Glen reagent Pac-dA-CE Phosphoramidite	10-1601-02	0.25 g	21,50	24,70	-	-	
Glen reagent Pac-dA-CE Phosphoramidite	10-1601-05	0.5 g	42,90	49,30	-	-	
Glen reagent Pac-dA-CE Phosphoramidite	10-1601-10	1 g	86,00	98,90	-	-	
Glen reagent Pac-dA-CE Phosphoramidite, Expedite	10-1601-05E	0.5 g	42,90	-	-	-	
Glen reagent Pac-dA-CPG	20-2601-01	0.1 g	25,80	29,70	-	-	

Description	Reference	Scale/Quantity	Price List 2018				Ice Shipping
			EUR	CHF	GBP	USD	
Glen reagent Pac-dA-CPG	20-2601-02	0.25 g	57,00	65,50	-	-	
Glen reagent Pac-dA-CPG	20-2601-10	1 g	215,00	247,00	-	-	
Glen reagent Pac-dA-CPG	20-2701-13	1 x 10 µmol	143,00	164,00	-	-	
Glen reagent Pac-dA-CPG	20-2701-41	4 x 1 µmol	86,00	98,90	-	-	
Glen reagent Pac-dA-CPG	20-2701-42	4 x 0.2 µmol	57,00	65,50	-	-	
Glen reagent Pac-dA-CPG	20-2701-45	4 x 40 nmol	57,00	65,50	-	-	
Glen reagent Pac-dA-CPG	20-2801-14	1 x 15 µmol	215,00	247,00	-	-	
Glen reagent Pac-dA-CPG	20-2801-41	4 x 1 µmol	86,00	98,90	-	-	
Glen reagent Pac-dA-CPG	20-2801-42	4 x 0.2 µmol	57,00	65,50	-	-	
Glen reagent Pac-dA-CPG	20-2801-45	4 x 40 nmol	57,00	65,50	-	-	
Glen reagent Pac-dA-Me Phosphoramidite	10-1301-02	0.25 g	35,80	41,10	-	-	
Glen reagent Pac-dA-Me Phosphoramidite	10-1301-05	0.5 g	72,00	82,80	-	-	
Glen reagent Pac-dA-Me Phosphoramidite	10-1301-10	1 g	143,00	164,00	-	-	
Glen reagent PC Amino-Modifier Phosphoramidite	10-4906-02	0.25 g	565,00	649,00	-	-	
Glen reagent PC Amino-Modifier Phosphoramidite	10-4906-90	100 µmol	193,00	222,00	-	-	
Glen reagent PC Biotin Phosphoramidite	10-4950-02	0.25 g	966,00	1.110,00	-	-	
Glen reagent PC Biotin Phosphoramidite	10-4950-90	100 µmol	358,00	411,00	-	-	
Glen reagent PC Biotin Phosphoramidite	10-4950-95	50 µmol	193,00	222,00	-	-	
Glen reagent PC Linker Phosphoramidite	10-4920-02	0.25 g	1.138,00	1.308,00	-	-	
Glen reagent PC Linker Phosphoramidite	10-4920-90	100 µmol	365,00	420,00	-	-	
Glen reagent PC Spacer Phosphoramidite	10-4913-02	0.25 g	565,00	649,00	-	-	
Glen reagent PC Spacer Phosphoramidite	10-4913-90	100 µmol	193,00	222,00	-	-	
Glen reagent pdC-CE Phosphoramidite	10-1014-02	0.25 g	351,00	403,00	-	-	
Glen reagent pdC-CE Phosphoramidite	10-1014-05	0.5 g	701,00	806,00	-	-	
Glen reagent pdC-CE Phosphoramidite	10-1014-90	100 µmol	122,00	140,00	-	-	
Glen reagent pdU-CE Phosphoramidite	10-1054-02	0.25 g	279,00	321,00	-	-	
Glen reagent pdU-CE Phosphoramidite	10-1054-05	0.5 g	558,00	641,00	-	-	
Glen reagent Perylene-dU	10-1591-02	0.25 g	1.031,00	1.185,00	-	-	
Glen reagent Perylene-dU	10-1591-90	100 µmol	429,00	493,00	-	-	
Glen reagent Perylene-dU	10-1591-95	50 µmol	215,00	247,00	-	-	
Glen reagent PseudoUridine-CE Phosphoramidite	10-3055-02	0.25 g	1.424,00	1.637,00	-	-	
Glen reagent PseudoUridine-CE Phosphoramidite	10-3055-90	100 µmol	501,00	576,00	-	-	
Glen reagent PseudoUridine-CE Phosphoramidite	10-3055-95	50 µmol	251,00	289,00	-	-	
Glen reagent Psoralen C2 Phosphoramidite	10-1982-02	0.25 g	709,00	815,00	-	-	
Glen reagent Psoralen C2 Phosphoramidite	10-1982-90	100 µmol	279,00	321,00	-	-	
Glen reagent Psoralen C6 Phosphoramidite	10-1983-02	0.25 g	709,00	815,00	-	-	
Glen reagent Psoralen C6 Phosphoramidite	10-1983-90	100 µmol	279,00	321,00	-	-	
Glen reagent Purine riboside α-thiotriphosphate (20mM) (PR-TTPαS)	80-3301-01	0.1 mL	143,00	164,00	-	-	
Glen reagent Puromycin-CPG	20-4040-01	0.1 g	172,00	198,00	-	-	
Glen reagent Puromycin-CPG	20-4040-10	1 g	1.424,00	1.637,00	-	-	
Glen reagent Puromycin-CPG	20-4140-13	1 x 10 µmol	515,00	592,00	-	-	
Glen reagent Puromycin-CPG	20-4140-14	1 x 15 µmol	773,00	889,00	-	-	
Glen reagent Puromycin-CPG	20-4140-41	4 x 1 µmol	286,00	329,00	-	-	
Glen reagent Puromycin-CPG	20-4140-42	4 x 0.2 µmol	172,00	198,00	-	-	
Glen reagent Pyrene-dU	10-1590-02	0.25 g	787,00	905,00	-	-	
Glen reagent Pyrene-dU	10-1590-90	100 µmol	301,00	346,00	-	-	
Glen reagent Pyrene-dU	10-1590-95	50 µmol	150,00	172,00	-	-	
Glen reagent Pyridin-2-one-CE Phosphoramidite	10-3012-02	0.25 g	1.718,00	1.975,00	-	-	
Glen reagent Pyridin-2-one-CE Phosphoramidite	10-3012-90	100 µmol	601,00	691,00	-	-	
Glen reagent Pyridin-2-one-CE Phosphoramidite	10-3012-95	50 µmol	301,00	346,00	-	-	
Glen reagent Pyrrolidine CE Phosphoramidite	10-1915-02	0.25 g	1.553,00	1.785,00	-	-	
Glen reagent Pyrrolidine CE Phosphoramidite	10-1915-90	100 µmol	544,00	625,00	-	-	
Glen reagent Pyrrolidine CE Phosphoramidite	10-1915-95	50 µmol	272,00	313,00	-	-	
Glen reagent Pyrrolo-CTP (10mM)	81-3017-01	0.1 mL	386,00	444,00	-	-	✓
Glen reagent Pyrrolo-dC-CE Phosphoramidite	10-1017-02	0.25 g	966,00	1.110,00	-	-	
Glen reagent Pyrrolo-dC-CE Phosphoramidite	10-1017-90	100 µmol	315,00	362,00	-	-	
Glen reagent Pyrrolo-dC-CE Phosphoramidite	10-1017-95	50 µmol	157,00	180,00	-	-	
Glen reagent Pyrrolo-dCTP (10mM)	81-1017-01	0.1 mL	215,00	247,00	-	-	
Glen reagent Quenching buffer for Glen-Pak RNA purification	60-4120-80	1 l	220,00	253,00	-	-	
Glen reagent Quenching buffer for Glen-Pak RNA purification	60-4120-82	250 mL	88,00	101,10	-	-	
Glen reagent Replacement Filters-Expedite, Pk/20	20-0021-0F		17,60	20,20	-	-	
Glen reagent Replacement Frits - TWIST 10um/15um, Pk/20	20-0040-0F		26,40	30,30	-	-	
Glen reagent rSpacer CE Phosphoramidite	10-3914-02	0.25 g	709,00	815,00	-	-	
Glen reagent rSpacer CE Phosphoramidite	10-3914-90	100 µmol	258,00	297,00	-	-	
Glen reagent rSpacer CE Phosphoramidite	10-3914-95	50 µmol	129,00	148,00	-	-	
Glen reagent SIMA (HEX) Phosphoramidite	10-5905-90	100 µmol	251,00	289,00	-	-	
Glen reagent Solid Chemical Phosphorylation Reagent II	10-1902-02	0.25 g	286,00	329,00	-	-	
Glen reagent Solid Chemical Phosphorylation Reagent II	10-1902-90	100 µmol	86,00	98,90	-	-	
Glen reagent Spacer C12 CE Phosphoramidite	10-1928-02	0.25 g	344,00	395,00	-	-	
Glen reagent Spacer C12 CE Phosphoramidite	10-1928-90	100 µmol	136,00	156,00	-	-	
Glen reagent Spacer Phosphoramidite 18	10-1918-02	0.25 g	317,00	364,00	-	-	
Glen reagent Spacer Phosphoramidite 18	10-1918-90	100 µmol	126,00	145,00	-	-	
Glen reagent Spacer Phosphoramidite 9	10-1909-02	0.25 g	344,00	395,00	-	-	
Glen reagent Spacer Phosphoramidite 9	10-1909-90	100 µmol	107,00	123,00	-	-	
Glen reagent Spacer Phosphoramidite C3	10-1913-02	0.25 g	344,00	395,00	-	-	
Glen reagent Spacer Phosphoramidite C3	10-1913-90	100 µmol	107,00	123,00	-	-	
Glen reagent Spacer Phosphoramidite C3, Expedite	10-1913-90E	100 µmol	107,00	123,00	-	-	
Glen reagent Sulfurizing Reagent	40-4036-10	1 g	38,50	44,30	-	-	
Glen reagent Sulfurizing Reagent	40-4036-20	2 g	77,00	88,50	-	-	
Glen reagent Symmetric Doubler Phosphoramidite	10-1920-02	0.25 g	344,00	395,00	-	-	
Glen reagent Symmetric Doubler Phosphoramidite	10-1920-90	100 µmol	215,00	247,00	-	-	
Glen reagent TAMRA NHS Ester	50-5910-66	60 µL	345,00	397,00	-	-	
Glen reagent TAMRA-dT	10-1057-02	0.25 g	1.396,00	1.605,00	-	-	
Glen reagent TAMRA-dT	10-1057-90	100 µmol	709,00	815,00	-	-	
Glen reagent TAMRA-dT	10-1057-95	50 µmol	358,00	411,00	-	-	
Glen reagent tC ⁻ -CE Phosphoramidite	10-1517-02	250 mg	2.090,00	2.402,00	-	-	✓
Glen reagent tC ⁻ -CE Phosphoramidite	10-1517-90	100 µmol	701,00	806,00	-	-	✓
Glen reagent tC ⁻ -CE Phosphoramidite	10-1517-95	50 µmol	358,00	411,00	-	-	✓
Glen reagent Thiol-Modifier C6 S-S	10-1936-02	0.25 g	476,00	547,00	-	-	
Glen reagent Thiol-Modifier C6 S-S	10-1936-90	100 µmol	198,00	228,00	-	-	
Glen reagent Thiol-Modifier C6 S-S, Cycl.	10-1936-90C	100 µmol	203,00	233,00	178,00	-	
Glen reagent Thymidine α-thiotriphosphate (15mM) (TTPαS)	80-1030-01	0.1 mL	143,00	164,00	-	-	
Glen reagent Thymidine Glycol CE Phosphoramidite	10-1096-02	0.25 g	1.396,00	1.605,00	-	-	

Description	Reference	Scale/Quantity	Price List 2018				Ice Shipping
			EUR	CHF	GBP	USD	
Glen reagent Thymidine Glycol CE Phosphoramidite	10-1096-90	100 µmol	515,00	592,00	-	-	
Glen reagent Thymidine Glycol CE Phosphoramidite	10-1096-95	50 µmol	258,00	297,00	-	-	
Glen reagent TMP-F-dU-CE Phosphoramidite	10-1016-02	0.25 g	709,00	815,00	-	-	
Glen reagent TMP-F-dU-CE Phosphoramidite	10-1016-90	100 µmol	279,00	321,00	-	-	
Glen reagent Trebler Phosphoramidite	10-1922-02	0.25 g	429,00	493,00	-	-	
Glen reagent Trebler Phosphoramidite	10-1922-90	100 µmol	258,00	297,00	-	-	
Glen reagent Trimer Phosphoramidite Mix 2	13-1992-90	100 µmol	1.474,00	1.694,00	-	-	
Glen reagent Trimer Phosphoramidite Mix 2	13-1992-95	50 µmol	737,00	847,00	-	-	
Glen reagent UniCap Phosphoramidite	10-4410-02	0.25 g	72,00	82,80	-	-	
Glen reagent UniCap Phosphoramidite	10-4410-05	0.5 g	143,00	164,00	-	-	
Glen reagent UniCap Phosphoramidite	10-4410-10	1 g	286,00	329,00	-	-	
Glen reagent UniCap Phosphoramidite	10-4410-20	2 g	573,00	659,00	-	-	
Glen reagent Universal Support II	20-5010-01	0.1 g	12,10	13,90	-	-	
Glen reagent Universal Support II	20-5010-02	0.25 g	27,50	31,60	-	-	
Glen reagent Universal Support II	20-5110-13	1 x 10 µmol	88,00	101,10	-	-	
Glen reagent Universal Support II	20-5110-41	4 x 1 µmol	53,00	60,90	-	-	
Glen reagent Universal Support II	20-5110-42	4 x 0.2 µmol	35,20	40,50	-	-	
Glen reagent Universal Support II	20-5110-45	4 x 40 nmol	35,20	40,50	-	-	
Glen reagent Universal Support II	20-5110-92	96 x 0.2 µmol	385,00	443,00	-	-	
Glen reagent Universal Support II	20-5110-95	96 x 40 nmol	385,00	443,00	-	-	
Glen reagent Universal Support II	20-5210-14	1 x 15 µmol	132,00	152,00	-	-	
Glen reagent Universal Support II	20-5210-41	4 x 1 µmol	53,00	60,90	-	-	
Glen reagent Universal Support II	20-5210-42	4 x 0.2 µmol	35,20	40,50	-	-	
Glen reagent Universal Support II	20-5210-45	4 x 40 nmol	35,20	40,50	-	-	
Glen reagent Universal Support II PS	26-5110-52	10 x 200nmol	88,00	101,10	-	-	
Glen reagent Universal Support II PS	26-5110-55	10x40 nmol	88,00	101,10	-	-	
Glen reagent Universal Support III	20-5010-10	1 g	105,00	121,00	-	-	
Glen reagent Universal Support III PS	27-5010-01	0.1 g	16,50	19,00	-	-	
Glen reagent Universal Support III PS	27-5010-02	0.25 g	38,50	44,30	-	-	
Glen reagent Universal Support III PS	27-5010-10	1 g	105,00	121,00	-	-	
Glen reagent Universal Support III PS	27-5110-92	96 x 0.2 µmol	385,00	443,00	-	-	
Glen reagent Universal Support III PS	27-5110-95	96 x 40 nmol	385,00	443,00	-	-	
Glen reagent Universal Support III PS	27-5110-P1	10 x 1.3 µmol	132,00	152,00	-	-	
Glen reagent Universal Support III PS	27-5110-P2	10 x 0.2 µmol	88,00	101,10	-	-	
Glen reagent Uridine α-thiotriphosphate (0.5mM) (UTTPaS)	80-3040-01	0.1 mL	143,00	164,00	-	-	
Glen reagent Zebularine-CE Phosphoramidite	10-3011-02	0.25 g	930,00	1.069,00	-	-	
Glen reagent Zebularine-CE Phosphoramidite	10-3011-90	100 µmol	358,00	411,00	-	-	
Glen reagent Zebularine-CE Phosphoramidite	10-3011-95	50 µmol	179,00	206,00	-	-	
Glen UnySupport 500, Columns	20-5240-41	4 x 1 µmol	53,00	60,90	-	-	
Methylene Blue C3 Phosphoramidite	10-5960-90	100 µmoles	760,00	-	-	-	
Methylene Blue C3 Phosphoramidite	10-5960-95	50 µmol	389,00	-	-	-	
N2-Amino-Modifier C6 dG	10-1529-02	0.25 g	1.581,00	1.817,00	-	-	✓
N2-Amino-Modifier C6 dG	10-1529-90	100 µM	499,00	-	-	-	
N4-Ac-N4-Et-dC-CE Phosphoramidite	10-1513-90	100 µmol	323,00	371,00	-	-	✓
N6-Ac-N6-Me-dA-CE Phosphoramidite	10-1503-02	0.25 g	515,00	-	-	-	
N6-Ac-N6-Me-dA-CE Phosphoramidite	10-1503-90	100 µmol	169,00	-	-	-	
N6-Methyl-A-CE Phosphoramidite	10-3005-95	50 µmoles	412,00	-	-	-	
N-POM Caged-dT-CE Phosphoramidite	10-1534-90	100 µmoles	520,00	598,00	-	-	
Protected Biotin Serinol Phosphoramidite	10-1993-02	0.25 g	966,00	1.110,00	-	-	
Protected Biotin Serinol Phosphoramidite	10-1993-90	100 µmol	422,00	485,00	-	-	
Protected Biotin Serinol Phosphoramidite	10-1993-95	50 µmol	236,00	271,00	-	-	
Protected BiotinLC Serinol Phosphoramidite	10-1995-02	0.25 g	966,00	1.110,00	-	-	
Protected BiotinLC Serinol Phosphoramidite	10-1995-90	100 µmol	522,00	600,00	-	-	
Protected BiotinLC Serinol Phosphoramidite	10-1995-95	50 µmol	293,00	337,00	-	-	
Psoralen Azide	50-2009-92	25 µmoles	161,00	-	-	-	
S-Bz-Thiol-Modifier C6-dT	10-1538-95	50 µmoles	241,00	-	-	-	
S-Bz-Thiol-Modifier C6-dT	10-1538-90	100 µmol	464,00	-	-	-	
SIMA (HEX) Phosphoramidite	10-5905-02	0.25 g	573,00	659,00	-	-	✓
Spermine Phosphoramidite	10-1939-02	0.25 g	760,00	-	-	-	
Spermine Phosphoramidite	10-1939-90	100 µmol	391,00	-	-	-	
Spermine Phosphoramidite	10-1939-95	50 µmol	210,00	-	-	-	
sTP (10mM)	81-3522-02	0.025 mL	501,00	576,00	-	-	
tC-CE Phosphoramidite	10-1516-95	50 µmoles	313,00	-	-	-	
tNitro-CE Phosphoramidite	10-1518-95	50 µmoles	328,00	-	-	-	
TEMPO Azide, 25 µmole	50-2007-92		143,00	-	-	-	
THPTA	50-1004-90	100 µmole	260,00	-	-	-	
Universal Support III HybridCPGTM	28-5010-10	1.0 g	139,00	-	-	-	
Universal Support III PS	26-5210-41	4x1,0µm	53,00	60,90	-	-	
Universal Support III PS	26-5110-42	4x0,2µmoles	34,40	-	-	-	
Universal Support III PS	26-5210-42	4 x 0,2µm	36,00	-	-	-	
α-Tocopherol-TEG Phosphoramidite	10-1977-02	0.25 µg	826,00	949,00	-	-	✓
α-Tocopherol-TEG Phosphoramidite	10-1977-90	100 µmol	431,00	495,00	-	-	✓
α-Tocopherol-TEG Phosphoramidite	10-1977-95	50 µmol	230,00	264,00	-	-	✓
GlenRNA Synthesis							EA1E1Z
2'-OMe-ibu-G-CE Phosphoramidite, Expedite	10-3120-02E	0,25 grams	32,00	36,80	-	-	
2'-OMe-G-Thiophosphoramidite	10-3172-02	0,25 grams	1.041,00	-	-	-	
2'-OMe-U-Thiophosphoramidite	10-3173-02	0,25 grams	1.041,00	-	-	-	
2'-F-Ac-C-CE Phosphoramidite	10-3415-05	0.5 g	143,00	164,00	-	-	
2'-F-Ac-C-CE Phosphoramidite, Expedite	10-3415-02E	0,25 grams	72,00	-	-	-	
2'-F-A-CE Phosphoramidite, Expedite	10-3400-02E	0,25 grams	145,00	-	-	-	
2'-F-G-CE Phosphoramidite, Expedite	10-3420-02E	0,25 grams	145,00	-	-	-	
2'-F-U-CE Phosphoramidite, Expedite	10-3430-02E	0,25 grams	72,00	-	-	-	
2'-OMe-5-Br-U-CE Phosphoramidite in bulk quantities > 10 g	10-3190-BULK	1 g	2.690,00	3.092,00	-	-	✓
2'-OMe-A-Thiophosphoramidite	10-3170-02	0,25 grams	1.041,00	-	-	-	
2'-OMe-C-CE Phosphoramidite, Expedite	10-3110-02E	0,25 grams	50,00	-	-	-	
2'-OMe-G-CE Phosphoramidite, Expedite	10-3121-02E	0,25 grams	50,00	-	-	-	
Ac-A-RNA-CPG	20-3304-01	0.1 g	57,00	65,50	-	-	
Ac-A-RNA-CPG	20-3304-10	0.1 g	508,00	584,00	-	-	
Ac-G-RNA-CPG	20-3324-01	0.1 g	57,00	65,50	-	-	
Ac-G-RNA-CPG	20-3324-10	1 g	508,00	584,00	-	-	
Glen 2'-oMe-ibu-G-CE Phosphoramidite	10-3120-02	0,25gr	61,00	70,10	-	-	
Glen 3'-Propargyl-5-Me-dC CPG	20-2982-41	4 X 1 µmol column	380,00	437,00	-	-	
Glen 3'-Propargyl-5-Me-dC CPG	20-2982-42	4 X 0,2 µmol column	190,00	218,00	-	-	

Description	Reference	Scale/Quantity	Price List 2018				Ice Shipping
			EUR	CHF	GBP	USD	
			Glen reagent 2,6-Diaminopurine-TOM-CE Phosphoramidite	10-3085-02	0.25 g	1,396,00	
Glen reagent 2-Aminopurine-TBDMs-CE Phosphoramidite	10-3070-02	0.25 g	1,396,00	1,605,00	-	-	
Glen reagent 2-Aminopurine-TBDMs-CE Phosphoramidite	10-3070-90	100 µmol	608,00	699,00	-	-	
Glen reagent 2-Aminopurine-TBDMs-CE Phosphoramidite	10-3070-95	50 µmol	304,00	349,00	-	-	
Glen reagent 2'-F-Ac-C-CE Phosphoramidite	10-3415-02	0.25 g	72,00	82,80	-	-	
Glen reagent 2'-F-Ac-C-CE Phosphoramidite	10-3415-90	100 µmol	241,00	277,00	-	-	
Glen reagent 2'-Fluoro-Adenosine a-thiotriphosphate (10mM) (2'-F-ATTPaS)	80-1102-01	0.1 mL	143,00	164,00	-	-	
Glen reagent 2'-F-U-CE Phosphoramidite	10-3430-02	0.25 g	72,00	82,80	-	-	
Glen reagent 2'-F-U-CE Phosphoramidite	10-3430-90	100 µmol	241,00	277,00	-	-	
Glen reagent 2'-OMe-2,6-Diaminopurine-CE Phosphoramidite	10-3124-02	0.25 g	1,396,00	1,605,00	-	-	
Glen reagent 2'-OMe-2,6-Diaminopurine-CE Phosphoramidite	10-3124-90	100 µmol	508,00	584,00	-	-	
Glen reagent 2'-OMe-2,6-Diaminopurine-CE Phosphoramidite	10-3124-95	50 µmol	254,00	292,00	-	-	
Glen reagent 2'-OMe-2-Aminopurine-CE Phosphoramidite	10-3123-02	0.25 g	1,396,00	1,605,00	-	-	
Glen reagent 2'-OMe-2-Aminopurine-CE Phosphoramidite	10-3123-90	100 µmol	508,00	584,00	-	-	
Glen reagent 2'-OMe-2-Aminopurine-CE Phosphoramidite	10-3123-95	50 µmol	254,00	292,00	-	-	
Glen reagent 2'-OMe-3-deaza-5-aza-C-CE Phosphoramidite	10-3116-02	0.25 g	1,396,00	1,605,00	-	-	
Glen reagent 2'-OMe-3-deaza-5-aza-C-CE Phosphoramidite	10-3116-90	100 µmol	508,00	584,00	-	-	
Glen reagent 2'-OMe-3-deaza-5-aza-C-CE Phosphoramidite	10-3116-95	50 µmol	254,00	292,00	-	-	
Glen reagent 2'-OMe-5-Br-U-CE Phosphoramidite	10-3190-02	0.25 g	966,00	1,110,00	-	-	
Glen reagent 2'-OMe-5-Br-U-CE Phosphoramidite	10-3190-90	100 µmol	344,00	395,00	-	-	
Glen reagent 2'-OMe-5-F-U-CE Phosphoramidite	10-3132-02	0.25 g	1,396,00	1,605,00	-	-	
Glen reagent 2'-OMe-5-F-U-CE Phosphoramidite	10-3132-90	100 µmol	508,00	584,00	-	-	
Glen reagent 2'-OMe-5-F-U-CE Phosphoramidite	10-3132-95	50 µmol	254,00	292,00	-	-	
Glen reagent 2'-OMe-5-Me-C-CE Phosphoramidite	10-3160-02	0.25 g	966,00	1,110,00	-	-	
Glen reagent 2'-OMe-5-Me-C-CE Phosphoramidite	10-3160-90	100 µmol	344,00	395,00	-	-	
Glen reagent 2'-OMe-5-Me-U-CE Phosphoramidite	10-3131-02	0.25 g	515,00	592,00	-	-	
Glen reagent 2'-OMe-5-Me-U-CE Phosphoramidite	10-3131-90	100 µmol	215,00	247,00	-	-	
Glen reagent 2'-OMe-Ac-C-CE Phosphoramidite	10-3115-02	0.25 g	61,00	70,10	-	-	
Glen reagent 2'-OMe-Ac-C-CE Phosphoramidite	10-3115-05	0.5 g	121,00	139,00	-	-	
Glen reagent 2'-OMe-Ac-C-CE Phosphoramidite	10-3115-10	1 g	242,00	278,00	-	-	
Glen reagent 2'-OMe-Ac-C-CE Phosphoramidite	10-3115-90	100 µmol	24,20	27,80	-	-	
Glen reagent 2'-OMe-Ac-C-CE Phosphoramidite, Expedite	10-3115-02E	0.25 g	61,00	70,10	-	-	
Glen reagent 2'-OMe-Ac-C-CE Phosphoramidite, Expedite	10-3115-05E	0.5 g	121,00	139,00	-	-	
Glen reagent 2'-OMe-Ac-C-RNA-CPG	20-3615-01	0.1 g	57,00	65,50	-	-	
Glen reagent 2'-OMe-Ac-C-RNA-CPG	20-3615-02	0.25 g	136,00	156,00	-	-	
Glen reagent 2'-OMe-Ac-C-RNA-CPG	20-3615-10	1 g	508,00	584,00	-	-	
Glen reagent 2'-OMe-Ac-C-RNA-CPG	20-3715-13	1 x 10 µmol	322,00	370,00	-	-	
Glen reagent 2'-OMe-Ac-C-RNA-CPG	20-3715-14	1 x 15 µmol	429,00	493,00	-	-	
Glen reagent 2'-OMe-Ac-C-RNA-CPG	20-3715-41	4 x 1 µmol	143,00	164,00	-	-	
Glen reagent 2'-OMe-Ac-C-RNA-CPG	20-3715-42	4 x 0.2 µmol	107,00	123,00	-	-	
Glen reagent 2'-OMe-A-CE Phosphoramidite	10-3100-02E	0.25 g	61,00	70,10	-	-	
Glen reagent 2'-OMe-A-CE Phosphoramidite	10-3100-05	0.5 g	121,00	139,00	-	-	
Glen reagent 2'-OMe-A-CE Phosphoramidite	10-3100-10	1 g	242,00	278,00	-	-	
Glen reagent 2'-OMe-A-CE Phosphoramidite	10-3100-90	100 µmol	24,20	27,80	-	-	
Glen reagent 2'-OMe-A-CE Phosphoramidite, Expedite	10-3100-02	0.25 g	61,00	70,10	-	-	
Glen reagent 2'-OMe-A-CE Phosphoramidite, Expedite	10-3100-05E	0.5 g	121,00	139,00	-	-	
Glen reagent 2'-OMe-Adenosine a-thiotriphosphate (20mM) (2'-O-Me-ATTPaS)	80-1101-01	0.1 mL	143,00	164,00	-	-	
Glen reagent 2'-OMe-A-RNA-CPG	20-3600-01	0.1 g	57,00	65,50	-	-	
Glen reagent 2'-OMe-A-RNA-CPG	20-3600-02	0.25 g	136,00	156,00	-	-	
Glen reagent 2'-OMe-A-RNA-CPG	20-3600-10	1 g	508,00	584,00	-	-	
Glen reagent 2'-OMe-A-RNA-CPG	20-3700-13	1 x 10 µmol	322,00	370,00	-	-	
Glen reagent 2'-OMe-A-RNA-CPG	20-3700-14	1 x 15 µmol	429,00	493,00	-	-	
Glen reagent 2'-OMe-A-RNA-CPG	20-3700-41	4 x 1 µmol	143,00	164,00	-	-	
Glen reagent 2'-OMe-A-RNA-CPG	20-3700-42	4 x 0.2 µmol	107,00	123,00	-	-	
Glen reagent 2'-OMe-A-RNA-CPG, Expedite	20-3700-42E	4 x 0.2 µmol	107,00	-	-	-	
Glen reagent 2'-OMe-C-CE Phosphoramidite	10-3110-02	0.25 g	61,00	70,10	-	-	
Glen reagent 2'-OMe-C-CE Phosphoramidite	10-3110-05	0.5 g	121,00	139,00	-	-	
Glen reagent 2'-OMe-C-CE Phosphoramidite	10-3110-10	1 g	242,00	278,00	-	-	
Glen reagent 2'-OMe-C-CE Phosphoramidite	10-3110-90	100 µmol	24,20	27,80	-	-	
Glen reagent 2'-OMe-C-RNA-CPG	20-3610-01	0.1 g	57,00	65,50	-	-	
Glen reagent 2'-OMe-C-RNA-CPG	20-3610-02	0.25 g	136,00	156,00	-	-	
Glen reagent 2'-OMe-C-RNA-CPG	20-3610-10	1 g	508,00	584,00	-	-	
Glen reagent 2'-OMe-C-RNA-CPG	20-3710-13	1 x 10 µmol	322,00	370,00	-	-	
Glen reagent 2'-OMe-C-RNA-CPG	20-3710-14	1 x 15 µmol	429,00	493,00	-	-	
Glen reagent 2'-OMe-C-RNA-CPG	20-3710-41	4 x 1 µmol	143,00	164,00	-	-	
Glen reagent 2'-OMe-C-RNA-CPG	20-3710-42	4 x 0.2 µmol	107,00	123,00	-	-	
Glen reagent 2'-OMe-G-CE Phosphoramidite	10-3121-02	0.25 g	61,00	70,10	-	-	
Glen reagent 2'-OMe-G-CE Phosphoramidite	10-3121-05	0.5 g	121,00	139,00	-	-	
Glen reagent 2'-OMe-G-CE Phosphoramidite	10-3121-10	1 g	242,00	278,00	-	-	
Glen reagent 2'-OMe-G-CE Phosphoramidite	10-3121-90	100 µmol	24,20	27,80	-	-	
Glen reagent 2'-OMe-G-CE Phosphoramidite, Expedite	10-3121-05E	0.5 g	121,00	139,00	-	-	
Glen reagent 2'-OMe-G-RNA-CPG	20-3621-01	0.1 g	57,00	65,50	-	-	
Glen reagent 2'-OMe-G-RNA-CPG	20-3621-02	0.25 g	136,00	156,00	-	-	
Glen reagent 2'-OMe-G-RNA-CPG	20-3621-10	1 g	508,00	584,00	-	-	
Glen reagent 2'-OMe-G-RNA-CPG	20-3721-13	1 x 10 µmol	322,00	370,00	-	-	
Glen reagent 2'-OMe-G-RNA-CPG	20-3721-14	1 x 15 µmol	429,00	493,00	-	-	
Glen reagent 2'-OMe-G-RNA-CPG	20-3721-41	4 x 1 µmol	143,00	164,00	-	-	
Glen reagent 2'-OMe-G-RNA-CPG	20-3721-42	4 x 0.2 µmol	107,00	123,00	-	-	
Glen reagent 2'-OMe-I-CE Phosphoramidite	10-3140-02	0.25 g	515,00	592,00	-	-	
Glen reagent 2'-OMe-I-CE Phosphoramidite	10-3140-90	100 µmol	215,00	247,00	-	-	
Glen reagent 2'-OMe-iPr-Pac-G-CE Phosphoramidite	10-3621-02	0.25 g	90,00	103,40	-	-	
Glen reagent 2'-OMe-iPr-Pac-G-CE Phosphoramidite	10-3621-05	0.5 g	179,00	206,00	-	-	
Glen reagent 2'-OMe-iPr-Pac-G-CE Phosphoramidite	10-3621-10	1 g	358,00	411,00	-	-	
Glen reagent 2'-OMe-iPr-Pac-G-CE Phosphoramidite	10-3621-90	100 µmol	35,80	41,10	-	-	
Glen reagent 2'-OMe-Pac-A-CE Phosphoramidite	10-3601-02	0.25 g	90,00	103,40	-	-	
Glen reagent 2'-OMe-Pac-A-CE Phosphoramidite	10-3601-05	0.5 g	179,00	206,00	-	-	
Glen reagent 2'-OMe-Pac-A-CE Phosphoramidite	10-3601-90	100 µmol	35,80	41,10	-	-	
Glen reagent 2'-OMe-TMP-5-F-U-CE Phosphoramidite	10-3111-02	0.25 g	1,396,00	1,605,00	-	-	
Glen reagent 2'-OMe-TMP-5-F-U-CE Phosphoramidite	10-3111-90	100 µmol	508,00	584,00	-	-	
Glen reagent 2'-OMe-TMP-5-F-U-CE Phosphoramidite	10-3111-95	50 µmol	254,00	292,00	-	-	
Glen reagent 2'-OMe-U-CE Phosphoramidite	10-3130-02	0.25 g	61,00	70,10	-	-	
Glen reagent 2'-OMe-U-CE Phosphoramidite	10-3130-05	0.5 g	121,00	139,00	-	-	
Glen reagent 2'-OMe-U-CE Phosphoramidite	10-3130-10	1 g	242,00	278,00	-	-	

Description	Reference	Scale/Quantity	Price List 2018				Ice Shipping
			EUR	CHF	GBP	USD	
Glen reagent 2'-OMe-U-CE Phosphoramidite	10-3130-90	100 µmol	24,20	27,80	-	-	
Glen reagent 2'-OMe-U-CE Phosphoramidite, Expedite	10-3130-02E	0.25 g	61,00	70,10	-	-	
Glen reagent 2'-OMe-U-RNA-CPG	20-3630-02	0.25 g	136,00	156,00	-	-	
Glen reagent 2'-OMe-U-RNA-CPG	20-3630-10	1 g	508,00	584,00	-	-	
Glen reagent 2'-OMe-U-RNA-CPG	20-3730-13	1 x 10 µmol	322,00	370,00	-	-	
Glen reagent 2'-OMe-U-RNA-CPG	20-3730-14	1 x 15 µmol	429,00	493,00	-	-	
Glen reagent 2'-OMe-U-RNA-CPG	20-3730-41	4 x 1 µmol	143,00	164,00	-	-	
Glen reagent 2'-OMe-U-RNA-CPG	20-3730-42	4 x 0.2 µmol	107,00	123,00	-	-	
Glen reagent 2'-Se-Me-U-CE Phosphoramidite	10-3730-02	0.25 g	1.396,00	1.605,00	-	-	
Glen reagent 2'-Se-Me-U-CE Phosphoramidite	10-3730-90	100 µmol	501,00	576,00	-	-	
Glen reagent 2'-Se-Me-U-CE Phosphoramidite	10-3730-95	50 µmol	251,00	289,00	-	-	
Glen reagent 4-Thio-U-TOM-CE Phosphoramidite	10-3052-02	0.25 g	1.396,00	1.605,00	-	-	
Glen reagent 4-Thio-U-TOM-CE Phosphoramidite	10-3052-90	100 µmol	608,00	699,00	-	-	
Glen reagent 4-Thio-U-TOM-CE Phosphoramidite	10-3052-95	50 µmol	304,00	349,00	-	-	
Glen reagent 5-Br-U-CE Phosphoramidite	10-3090-02	0.25 g	680,00	782,00	-	-	
Glen reagent 5-Br-U-CE Phosphoramidite	10-3090-90	100 µmol	279,00	321,00	-	-	
Glen reagent 5-I-U-CE Phosphoramidite	10-3091-02	0.25 g	680,00	782,00	-	-	
Glen reagent 5-I-U-CE Phosphoramidite	10-3091-90	100 µmol	279,00	321,00	-	-	
Glen reagent 5-Me-C-TOM-CE Phosphoramidite	10-3064-02	0.25 g	680,00	782,00	-	-	
Glen reagent 5-Me-C-TOM-CE Phosphoramidite	10-3064-90	100 µmol	272,00	313,00	-	-	
Glen reagent 5-Me-C-TOM-CE Phosphoramidite	10-3064-95	50 µmol	136,00	156,00	-	-	
Glen reagent 5-Me-U-CE Phosphoramidite	10-3050-02	0.25 g	680,00	782,00	-	-	
Glen reagent 5-Me-U-CE Phosphoramidite	10-3050-90	100 µmol	272,00	313,00	-	-	
Glen reagent Ac-C-CE Phosphoramidite	10-3015-02	0.25 g	35,20	40,50	-	-	
Glen reagent Ac-C-CE Phosphoramidite	10-3015-05	0.5 g	53,00	60,90	-	-	
Glen reagent Ac-C-CE Phosphoramidite	10-3015-10	1 g	99,00	113,80	-	-	
Glen reagent Ac-C-CE Phosphoramidite, Expedite	10-3015-05E	0.5 g	53,00	60,90	-	-	
Glen reagent Ac-C-RNA-CPG	20-3315-01	0.1 g	57,00	65,50	-	-	
Glen reagent Ac-C-RNA-CPG	20-3315-02	0.25 g	136,00	156,00	-	-	
Glen reagent Ac-C-RNA-CPG	20-3315-10	1 g	508,00	584,00	-	-	
Glen reagent Ac-C-RNA-CPG	20-3415-13	1 x 10 µmol	322,00	370,00	-	-	
Glen reagent Ac-C-RNA-CPG	20-3415-14	1 x 15 µmol	429,00	493,00	-	-	
Glen reagent Ac-C-RNA-CPG	20-3415-41	4 x 1 µmol	143,00	164,00	-	-	
Glen reagent Ac-C-RNA-CPG	20-3415-42	4 x 0.2 µmol	107,00	123,00	-	-	
Glen reagent A-RNA-TOM-CPG	20-3304-02	0.25 g	136,00	156,00	-	-	
Glen reagent A-RNA-TOM-CPG	20-3404-13	1 x 10 µmol	322,00	370,00	-	-	
Glen reagent A-RNA-TOM-CPG	20-3404-14	1 x 15 µmol	429,00	493,00	-	-	
Glen reagent A-RNA-TOM-CPG	20-3404-41	4 x 1 µmol	143,00	164,00	-	-	
Glen reagent A-RNA-TOM-CPG	20-3404-42	4 x 0.2 µmol	107,00	123,00	-	-	
Glen reagent A-TOM-CE Phosphoramidite	10-3004-02	0.25 g	107,00	123,00	-	-	
Glen reagent A-TOM-CE Phosphoramidite	10-3004-05	0.5 g	165,00	190,00	-	-	
Glen reagent A-TOM-CE Phosphoramidite	10-3004-10	1 g	303,00	348,00	-	-	
Glen reagent Bz-A-CE Phosphoramidite	10-3003-02	0.25 g	35,20	40,50	-	-	
Glen reagent Bz-A-CE Phosphoramidite	10-3003-05	0.5 g	53,00	60,90	-	-	
Glen reagent Bz-A-CE Phosphoramidite	10-3003-10	1 g	99,00	113,80	-	-	
Glen reagent Bz-A-RNA-CPG	20-3303-01	0.1 g	57,00	65,50	-	-	
Glen reagent Bz-A-RNA-CPG	20-3303-02	0.25 g	136,00	156,00	-	-	
Glen reagent Bz-A-RNA-CPG	20-3303-10	1 g	508,00	584,00	-	-	
Glen reagent Bz-A-RNA-CPG	20-3403-13	1 x 10 µmol	322,00	370,00	-	-	
Glen reagent Bz-A-RNA-CPG	20-3403-14	1 x 15 µmol	429,00	493,00	-	-	
Glen reagent Bz-A-RNA-CPG	20-3403-41	4 x 1 µmol	143,00	164,00	-	-	
Glen reagent Bz-A-RNA-CPG	20-3403-42	4 x 0.2 µmol	107,00	123,00	-	-	
Glen reagent C-TOM-CE Phosphoramidite	10-3014-02	0.25 g	107,00	123,00	-	-	
Glen reagent C-TOM-CE Phosphoramidite	10-3014-05	0.5 g	165,00	190,00	-	-	
Glen reagent C-TOM-CE Phosphoramidite	10-3014-10	1 g	303,00	348,00	-	-	
Glen reagent dmf-G-RNA-CPG	20-3329-01	0.1 g	57,00	65,50	-	-	
Glen reagent dmf-G-RNA-CPG	20-3329-02	0.25 g	136,00	156,00	-	-	
Glen reagent dmf-G-RNA-CPG	20-3329-10	1 g	508,00	584,00	-	-	
Glen reagent dmf-G-RNA-CPG	20-3429-13	1 x 10 µmol	322,00	370,00	-	-	
Glen reagent dmf-G-RNA-CPG	20-3429-14	1 x 15 µmol	429,00	493,00	-	-	
Glen reagent dmf-G-RNA-CPG	20-3429-41	4 x 1 µmol	143,00	164,00	-	-	
Glen reagent dmf-G-RNA-CPG	20-3429-42	4 x 0.2 µmol	107,00	123,00	-	-	
Glen reagent G-RNA-TOM-CPG	20-3324-02	0.25 g	136,00	156,00	-	-	
Glen reagent G-RNA-TOM-CPG	20-3424-13	1 x 10 µmol	322,00	370,00	-	-	
Glen reagent G-RNA-TOM-CPG	20-3424-14	1 x 15 µmol	429,00	493,00	-	-	
Glen reagent G-RNA-TOM-CPG	20-3424-41	4 x 1 µmol	143,00	164,00	-	-	
Glen reagent G-RNA-TOM-CPG	20-3424-42	4 x 0.2 µmol	107,00	123,00	-	-	
Glen reagent G-TOM-CE Phosphoramidite	10-3024-02	0.25 g	107,00	123,00	-	-	
Glen reagent G-TOM-CE Phosphoramidite	10-3024-05	0.5 g	165,00	190,00	-	-	
Glen reagent G-TOM-CE Phosphoramidite	10-3024-10	1 g	303,00	348,00	-	-	
Glen reagent iPr-Pac-G-CE Phosphoramidite	10-3021-02	0.25 g	107,00	123,00	-	-	
Glen reagent iPr-Pac-G-CE Phosphoramidite	10-3021-05	0.5 g	215,00	247,00	-	-	
Glen reagent iPr-Pac-G-CE Phosphoramidite	10-3021-05E	0.5 g	215,00	247,00	-	-	
Glen reagent iPr-Pac-G-CE Phosphoramidite	10-3021-10	1 g	394,00	453,00	-	-	
Glen reagent iPr-Pac-G-RNA-CPG	20-3321-01	0.1 g	57,00	65,50	-	-	
Glen reagent iPr-Pac-G-RNA-CPG	20-3321-02	0.25 g	136,00	156,00	-	-	
Glen reagent iPr-Pac-G-RNA-CPG	20-3421-13	1 x 10 µmol	322,00	370,00	-	-	
Glen reagent iPr-Pac-G-RNA-CPG	20-3421-14	1 x 15 µmol	429,00	493,00	-	-	
Glen reagent iPr-Pac-G-RNA-CPG	20-3421-41	4 x 1 µmol	143,00	164,00	-	-	
Glen reagent iPr-Pac-G-RNA-CPG	20-3421-42	4 x 0.2 µmol	107,00	123,00	-	-	
Glen reagent Pac-A-CE Phosphoramidite	10-3000-02	0.25 g	107,00	123,00	-	-	
Glen reagent Pac-A-CE Phosphoramidite	10-3000-05	0.5 g	215,00	247,00	-	-	
Glen reagent Pac-A-CE Phosphoramidite	10-3000-10	1 g	394,00	453,00	-	-	
Glen reagent Pac-A-CE Phosphoramidite, Expedite	10-3000-05E	0.5 g	215,00	247,00	-	-	
Glen reagent Pac-A-RNA-CPG	20-3300-01	0.1 g	57,00	65,50	-	-	
Glen reagent Pac-A-RNA-CPG	20-3300-02	0.25 g	136,00	156,00	-	-	
Glen reagent Pac-A-RNA-CPG	20-3400-13	1 x 10 µmol	322,00	370,00	-	-	
Glen reagent Pac-A-RNA-CPG	20-3400-14	1 x 15 µmol	429,00	493,00	-	-	
Glen reagent Pac-A-RNA-CPG	20-3400-41	4 x 1 µmol	143,00	164,00	-	-	
Glen reagent Pac-A-RNA-CPG	20-3400-42	4 x 0.2 µmol	107,00	123,00	-	-	
Glen reagent Pyrrolo-C-TOM-CE Phosphoramidite	10-3017-02	0.25 g	1.396,00	1.605,00	-	-	
Glen reagent Pyrrolo-C-TOM-CE Phosphoramidite	10-3017-90	100 µmol	608,00	699,00	-	-	
Glen reagent Pyrrolo-C-TOM-CE Phosphoramidite	10-3017-95	50 µmol	304,00	349,00	-	-	

Description	Reference	Scale/Quantity	Price List 2018				Ice Shipping
			EUR	CHF	GBP	USD	
Glen reagent U-CE Phosphoramidite	10-3030-02	0.25 g	35,20	40,50	-	-	
Glen reagent U-CE Phosphoramidite	10-3030-05	0.5 g	53,00	60,90	-	-	
Glen reagent U-CE Phosphoramidite	10-3030-10	1 g	99,00	113,80	-	-	
Glen reagent U-CE Phosphoramidite, Expedite	10-3030-05E	0.5 g	53,00	60,90	-	-	
Glen reagent U-RNA-CPG	20-3330-01	0.1 g	57,00	65,50	-	-	
Glen reagent U-RNA-CPG	20-3330-02	0.25 g	136,00	156,00	-	-	
Glen reagent U-RNA-CPG	20-3330-10	1 g	508,00	584,00	-	-	
Glen reagent U-RNA-CPG	20-3430-13	1 x 10 µmol	322,00	370,00	-	-	
Glen reagent U-RNA-CPG	20-3430-14	1 x 15 µmol	429,00	493,00	-	-	
Glen reagent U-RNA-CPG	20-3430-41	4 x 1 µmol	143,00	164,00	-	-	
Glen reagent U-RNA-CPG	20-3430-42	4 x 0.2 µmol	107,00	123,00	-	-	
Glen reagent U-TOM-CE Phosphoramidite	10-3034-02	0.25 g	107,00	123,00	-	-	
Glen reagent U-TOM-CE Phosphoramidite	10-3034-05	0.5 g	165,00	190,00	-	-	
Glen reagent U-TOM-CE Phosphoramidite	10-3034-10	1 g	303,00	348,00	-	-	
Pseudo-Uridine-a-thiotriphosphate (1.0mM)	80-3341-01	0.1 mL	143,00	164,00	-	-	

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

A= Applied Biosystems 3900