

# Oligonucleotides purification table



Description	Reference	Purifications							Specifications						
		SePOP	RP- Cartridge Gold™	RP-HPLC	IEX-HPLC	PAGE	DUAL HPLC (RP+RP)	DUAL HPLC (RP+HPLC)	Ultra Pure Gold™	Position 5'	Position 3'	Post Synthesis Coupling*	Ab <sub>max</sub> (nm)	Em <sub>max</sub> (nm)	MW (g.mol <sup>-1</sup> )
<b>Custom oligonucleotides - Bases</b>															
<b>DNA Bases</b>															
DNA Base 10 nmol scale (15-59 bases)	BA-DN001-001	IN									✓	✓		-	-
DNA Base 40 nmol scale (10-99 bases)	BA-DN001-004	IN									✓	✓		-	-
DNA Base 200 nmol scale (5-139 bases)	BA-DN001-020	IN									✓	✓		-	-
DNA Base 1000 nmol scale (5-139 bases)	BA-DN001-100	IN									✓	✓		-	-
DNA Base 2,5 µmol scale (5-139 bases)	BA-DN001-M02	IN									✓	✓		-	-
DNA Base 5 µmol scale (5-139 bases)	BA-DN001-M05	IN									✓	✓		-	-
DNA Base 10 µmol scale (5-139 bases)	BA-DN001-M10	IN									✓	✓		-	-
DNA Base 20 µmol scale (5-139 bases)	BA-DN001-M20	IN									✓	✓		-	-
DNA Base + Phosphorothioate Link 40 nmol scale	BA-DN010-004	IN									✓	✓		-	-
DNA Base + Phosphorothioate Link 200 nmol scale	BA-DN010-020	IN									✓	✓		-	-
DNA Base + Phosphorothioate Link 1000 nmol scale	BA-DN010-100	IN									✓	✓		-	-
DNA Base + Phosphorothioate Link 2,5 µmol scale	BA-DN010-M02	IN									✓	✓		-	-
DNA Base + Phosphorothioate Link 5 µmol scale	BA-DN010-M05	IN									✓	✓		-	-
DNA Base + Phosphorothioate Link 10 µmol scale	BA-DN010-M10	IN									✓	✓		-	-
DNA Base + Phosphorothioate Link 20 µmol scale (5-139 bases)	BA-DN010-M20	IN									✓	✓		-	-
DNA Base + Methylphosphonate Link 40 nmol scale	BA-DN020-004										✓	✓		-	-
DNA Base + Methylphosphonate Link 200 nmol scale	BA-DN020-020										✓	✓		-	-
DNA Base + Methylphosphonate Link 1000 nmol scale	BA-DN020-100										✓	✓		-	-
<b>RNA Bases</b>															
RNA Base 40 nmol scale (15-29 bases)	BA-RN001-004	IN									✓	✓		-	-
RNA Base 200 nmol scale (8-59 bases)	BA-RN001-020	IN									✓	✓		-	-
RNA Base 1000 nmol scale (8-59 bases)	BA-RN001-100	IN									✓	✓		-	-
RNA Base 2,5 µmol scale (8-59 bases)	BA-RN001-M02	IN									✓	✓		-	-
RNA Base 5 µmol scale (8-59 bases)	BA-RN001-M05	IN									✓	✓		-	-
RNA Base 10 µmol scale (8-59 bases)	BA-RN001-M10	IN									✓	✓		-	-
RNA Base 20 µmol scale (8-59 bases)	BA-RN001-M20	IN									✓	✓		-	-
RNA Base + Phosphorothioate Link 40 nmol scale	BA-RN010-004	IN									✓	✓		-	-
RNA Base + Phosphorothioate Link 200 nmol scale	BA-RN010-020	IN									✓	✓		-	-
RNA Base + Phosphorothioate Link 1000 nmol scale	BA-RN010-100	IN									✓	✓		-	-
RNA Base + Phosphorothioate Link 2,5 µmol scale	BA-RN010-M02	IN									✓	✓		-	-
RNA Base + Phosphorothioate Link 5 µmol scale	BA-RN010-M05	IN									✓	✓		-	-
RNA Base + Phosphorothioate Link 10 µmol scale	BA-RN010-M10	IN									✓	✓		-	-
RNA Base + Phosphorothioate Link 20 µmol scale	BA-RN010-M20	IN									✓	✓		-	-
<b>LNA Bases</b>															
LNA Base 40 nmol scale	BA-LN001-004	IN									✓	✓		-	-
LNA Base 200 nmol scale	BA-LN001-020	IN									✓	✓		-	-
LNA Base 1000 nmol scale	BA-LN001-100	IN									✓	✓		-	-
LNA Base 2,5 µmol scale	BA-LN001-M02	IN									✓	✓		-	-
LNA Base 5 µmol scale	BA-LN001-M05	IN									✓	✓		-	-
LNA Base 10 µmol scale	BA-LN001-M10	IN									✓	✓		-	-
LNA Base 20 µmol scale	BA-LN001-M20	IN									✓	✓		-	-
LNA Base + Phosphorothioate Link 40 nmol scale	BA-LN010-004	IN									✓	✓		-	-
LNA Base + Phosphorothioate Link 200 nmol scale	BA-LN010-020	IN									✓	✓		-	-
LNA Base + Phosphorothioate Link 1000 nmol scale	BA-LN010-100	IN									✓	✓		-	-
LNA Base + Phosphorothioate Link 2500 nmol scale	BA-LN010-M02	IN									✓	✓		-	-
LNA Base + Phosphorothioate Link 5000 nmol scale	BA-LN010-M05	IN									✓	✓		-	-
LNA Base + Phosphorothioate Link 10000 nmol scale	BA-LN010-M10	IN									✓	✓		-	-
LNA Base + Phosphorothioate Link 20000 nmol scale	BA-LN010-M20	IN									✓	✓		-	-
<b>2'-OMe RNA Bases</b>															
2'-OMe RNA Base 40 nmol scale (10-99 bases)	BA-2M001-004	IN									✓	✓		-	-
2'-OMe RNA Base 200 nmol scale (5-139 bases)	BA-2M001-020	IN									✓	✓		-	-
2'-OMe RNA Base 1000 nmol scale (5-139 bases)	BA-2M001-100	IN									✓	✓		-	-
2'-OMe RNA Base 2,5 µmol scale (5-139 bases)	BA-2M001-M02	IN									✓	✓		-	-
2'-OMe RNA Base 5 µmol scale (5-139 bases)	BA-2M001-M05	IN									✓	✓		-	-
2'-OMe RNA Base 10 µmol scale (5-139 bases)	BA-2M001-M10	IN									✓	✓		-	-
2'-OMe RNA Base 20 µmol scale (5-139 bases)	BA-2M001-M20	IN									✓	✓		-	-
2'-OMe RNA Base + Phosphorothioate Link 40 nmol scale	BA-2M010-004	IN									✓	✓		-	-
2'-OMe RNA Base + Phosphorothioate Link 200 nmol scale	BA-2M010-020	IN									✓	✓		-	-
2'-OMe RNA Base + Phosphorothioate Link 1000 nmol scale	BA-2M010-100	IN									✓	✓		-	-
2'-OMe RNA Base + Phosphorothioate Link 2,5 µmol scale	BA-2M010-M02	IN									✓	✓		-	-
2'-OMe RNA Base + Phosphorothioate Link 5 µmol scale	BA-2M010-M05	IN									✓	✓		-	-
2'-OMe RNA Base + Phosphorothioate Link 10 µmol scale	BA-2M010-M10	IN									✓	✓		-	-
2'-OMe RNA Base + Phosphorothioate Link 20 µmol scale	BA-2M010-M20	IN									✓	✓		-	-
<b>2'-MOE RNA Bases</b>															
2'-MOE RNA Base 40 nmol scale (10-99 bases)	BA-ME001-004	IN									✓	✓		-	-
2'-MOE RNA Base 200 nmol scale (5-139 bases)	BA-ME001-020	IN									✓	✓		-	-
2'-MOE RNA Base 1000 nmol scale (5-139 bases)	BA-ME001-100	IN									✓	✓		-	-
2'-MOE RNA Base 2,5 µmol scale (5-139 bases)	BA-ME001-M02	IN									✓	✓		-	-
2'-MOE RNA Base 5 µmol scale (5-139 bases)	BA-ME001-M05	IN									✓	✓		-	-
2'-MOE RNA Base 10 µmol scale (5-139 bases)	BA-ME001-M10	IN									✓	✓		-	-
2'-MOE RNA Base 20 µmol scale (5-139 bases)	BA-ME001-M20	IN									✓	✓		-	-
2'-MOE RNA Base + Phosphorothioate Link 40 nmol scale (10-99 bases)	BA-ME010-004	IN									✓	✓		-	-
2'-MOE RNA Base + Phosphorothioate Link 200 nmol scale (5-139 bases)	BA-ME010-020	IN									✓	✓		-	-
2'-MOE RNA Base + Phosphorothioate Link 1000 nmol scale (5-139 bases)	BA-ME010-100	IN									✓	✓		-	-
2'-MOE RNA Base + Phosphorothioate Link 2,5 µmol scale (5-139 bases)	BA-ME010-M02	IN									✓	✓		-	-
2'-MOE RNA Base + Phosphorothioate Link 5 µmol scale (5-139 bases)	BA-ME010-M05	IN									✓	✓		-	-
2'-MOE RNA Base + Phosphorothioate Link 10 µmol scale (5-139 bases)	BA-ME010-M10	IN									✓	✓		-	-
2'-MOE RNA Base + Phosphorothioate Link 20 µmol scale (5-139 bases)	BA-ME010-M20	IN									✓	✓		-	-
<b>PNA Bases</b>															
PNA Base 50 nmol (6-23 bases), Unlabeled 90 % Purity	BA-PN001-005			IN							✓	✓		-	-
PNA Base 100 nmol (6-23 bases), Unlabeled 90 % Purity	BA-PN001-010			IN							✓	✓		-	-
PNA Base 50 nmol (6-23 bases), Unlabeled 95 % Purity	BA-PN010-005			IN							✓	✓		-	-
PNA Base 100 nmol (6-23 bases), Unlabeled 95 % Purity	BA-PN010-010			IN							✓	✓		-	-
<b>NGS Bases</b>															
NGS DNA Base 10 nmol (20-85 bases) + Cartridge purification	BA-NGSCA-001		IN								✓	✓		-	-
NGS DNA Base + Cartridge purification + Phosphorothioate Link	BA-NGSCA-PT-001		IN								✓	✓		-	-
NGS DNA Base 10 nmol (20-85 bases) + HPLC purification	BA-NGSHP-001			IN							✓	✓		-	-
NGS DNA Base + HPLC purification + Phosphorothioate Link	BA-NGSHP-PT-001			IN							✓	✓		-	-
<b>Custom oligonucleotides - Modifications</b>															
<b>Conjugates</b>															
3' Digoxigenin 40 nmol scale	MD-CN001-03004										✓	✓		680.77	
3' Digoxigenin 200 nmol scale	MD-CN001-03020										✓	✓		680.77	
3' Digoxigenin 1000 nmol scale	MD-CN001-03100										✓	✓		680.77	
5'-Digoxigenin 40 nmol scale	MD-CN001-05004										✓	✓		680.77	
5' Digoxigenin 200 nmol scale	MD-CN001-05020										✓	✓		680.77	
5' Digoxigenin 1000 nmol scale	MD-CN001-05100										✓	✓		680.77	
Digoxigenin (dR) 40 nmol scale	MD-CN001-DR004										✓	✓		766.86	
Digoxigenin (dR) 200 nmol scale	MD-CN001-DR020										✓	✓		766.86	
Digoxigenin (dR) 1000 nmol scale	MD-CN001-DR100										✓	✓		766.86	
Digoxigenin (dT) 40 nmol scale	MD-CN001-DT004										✓	✓		930	
Digoxigenin (dT) 200 nmol scale	MD-CN001-DT020										✓	✓		930	
Digoxigenin (dT) 1000 nmol scale	MD-CN001-DT100										✓	✓		930	
3' TEG-Dinitrophenol 40 nmol scale	MD-CN010-03004										✓	✓		509.41	

■ Required Purification\*   
■ Compatible Purification\*   
■ Best Purification\*

Description	Reference	Purifications								Specifications						
		SePOP	RP- Cartridge Gold™	RP-HPLC	IEX-HPLC	PAGE	DUAL HPLC (RP+RP)	DUAL HPLC (RP+HPLC)	Ultra Pure Gold™	Position			Ab <sub>max</sub> (nm)	Em <sub>max</sub> (nm)	MW (g·mol <sup>-1</sup> )	
										5'	IN	3'				
3' TEG-Dinitrophenol 200 nmol scale	MD-CN010-03020														509.41	
3' TEG-Dinitrophenol 1000 nmol scale	MD-CN010-03100														509.41	
5' TEG-Dinitrophenol 40 nmol scale	MD-CN010-05004														509.41	
5' TEG-Dinitrophenol 200 nmol scale	MD-CN010-05020														509.41	
5' TEG-Dinitrophenol 1000 nmol scale	MD-CN010-05100														509.41	
3' Glyceryl 40 nmol scale	MD-CN030-03004	IN													154.06	
3' Glyceryl 200 nmol scale	MD-CN030-03020	IN													154.06	
3' Glyceryl 1000 nmol scale	MD-CN030-03100	IN													154.06	
3' TEG-Cholesteryl 40 nmol scale	MD-CN040-03004	IN													755.97	
3' TEG-Cholesteryl 200 nmol scale	MD-CN040-03020	IN													755.97	
3' TEG-Cholesteryl 1000 nmol scale	MD-CN040-03100	IN													755.97	
5' TEG-Cholesteryl 40 nmol scale	MD-CN040-05004	IN													755.97	
5' TEG-Cholesteryl 200 nmol scale	MD-CN040-05020	IN													755.97	
5' TEG-Cholesteryl 1000 nmol scale	MD-CN040-05100	IN													755.97	
Cholesteryl (dR) 40 nmol scale	MD-CN040-DR004	IN													706.44	
Cholesteryl (dR) 200 nmol scale	MD-CN040-DR020	IN													706.44	
Cholesteryl (dR) 1000 nmol scale	MD-CN040-DR100	IN													706.44	
3' MeO-Cl-Acridine 40 nmol scale	MD-CN050-03004	IN													450.86	
3' MeO-Cl-Acridine 200 nmol scale	MD-CN050-03020	IN													450.86	
3' MeO-Cl-Acridine 1000 nmol scale	MD-CN050-03100	IN													450.86	
5' MeO-Cl-Acridine 40 nmol scale	MD-CN050-05004	IN													450.86	
5' MeO-Cl-Acridine 200 nmol scale	MD-CN050-05020	IN													450.86	
5' MeO-Cl-Acridine 1000 nmol scale	MD-CN050-05100	IN													450.86	
Internal MeO-Cl-Acridine 40 nmol scale	MD-CN050-IN004	IN													450.86	
Internal MeO-Cl-Acridine 200 nmol scale	MD-CN050-IN020	IN													450.86	
Internal MeO-Cl-Acridine 1000 nmol scale	MD-CN050-IN100	IN													450.86	
5' C2 Psoralen 40 nmol scale	MD-CN062-05004	IN													364.29	
5' C2 Psoralen 200 nmol scale	MD-CN062-05020	IN													364.29	
5' C2 Psoralen 1000 nmol scale	MD-CN062-05100	IN													364.29	
5' C6 Psoralen 40 nmol scale	MD-CN066-05004	IN													420.4	
5' C6 Psoralen 200 nmol scale	MD-CN066-05020	IN													420.4	
5' C6 Psoralen 1000 nmol scale	MD-CN066-05100	IN													420.4	
Carboxy (dT) 40 nmol scale	MD-CN070-DT004	IN													360.22	
Carboxy (dT) 200 nmol scale	MD-CN070-DT020	IN													360.22	
Carboxy (dT) 1000 nmol scale	MD-CN070-DT100	IN													360.22	
5' AP Conjugation 2.5 μmol scale	MD-CN090-05M02	IN														
5' HRP Conjugation 2.5 μmol scale	MD-CN092-05M02	IN														
5' BSA Conjugation 2.5 μmol scale	MD-CN094-05M02	IN														
5' SBP Conjugation 2.5 μmol scale	MD-CN096-05M02	IN														
5' Peptide Conjugation 2.5 μmol scale	MD-CN099-05M02	IN														
3' Oligo Affinity Support (PS) 1000 nmol scale	MD-CN100-03100	IN														
<b>Degenerate bases - Wobblers</b>																
3' Wobble 40 nmol scale	MD-DB001-03004	IN														KB2B12
3' Wobble 200 nmol scale	MD-DB001-03020	IN														
3' Wobble 1000 nmol scale	MD-DB001-03100	IN														
Spiked Oligonucleotide (1-4 spikes) 200 nmol scale	MD-DB002-IN020	IN														
Spiked Oligonucleotide (1-5 spikes) 1000 nmol scale	MD-DB002-IN100	IN														
<b>Fluorescent dyes</b>																
3' 6-FAM 40 nmol scale	MD-FL001-03004	IN														KB2C12
3' 6-FAM 200 nmol scale	MD-FL001-03020	IN														
3' 6-FAM 1000 nmol scale	MD-FL001-03100	IN														
5' 6-FAM 40 nmol scale	MD-FL001-05004	IN														
5' 6-FAM 200 nmol scale	MD-FL001-05020	IN														
5' 6-FAM 1000 nmol scale	MD-FL001-05100	IN														
5' 6-FAM 2.5 μmol scale	MD-FL001-05M02	IN														
5' 6-FAM 5 μmol scale	MD-FL001-05M05	IN														
5' 6-FAM 10 μmol scale	MD-FL001-05M10	IN														
6-FAM (dR) 40 nmol scale	MD-FL001-DR004	IN														
6-FAM (dR) 200 nmol scale	MD-FL001-DR020	IN														
6-FAM (dR) 1000 nmol scale	MD-FL001-DR100	IN														
6-FAM (dT) 40 nmol scale	MD-FL001-DT004	IN														
6-FAM (dT) 200 nmol scale	MD-FL001-DT020	IN														
6-FAM (dT) 1000 nmol scale	MD-FL001-DT100	IN														
6-FAM (dT) 2.5 μmol scale	MD-FL001-DTM02	IN														
6-FAM (dT) 5 μmol scale	MD-FL001-DTM05	IN														
6-FAM (dT) 10 μmol scale	MD-FL001-DTM10	IN														
3' Fluorescein 40 nmol scale	MD-FL005-03004	IN														
3' Fluorescein 200 nmol scale	MD-FL005-03020	IN														
3' Fluorescein 1000 nmol scale	MD-FL005-03100	IN														
5' Fluorescein 40 nmol scale	MD-FL005-05004	IN														
5' Fluorescein 200 nmol scale	MD-FL005-05020	IN														
5' Fluorescein 1000 nmol scale	MD-FL005-05100	IN														
5' HEX 40 nmol scale	MD-FL010-05004	IN														
5' HEX 200 nmol scale	MD-FL010-05020	IN														
5' HEX 1000 nmol scale	MD-FL010-05100	IN														
HEX (dR) 40 nmol scale	MD-FL010-DR004	IN														
HEX (dR) 200 nmol scale	MD-FL010-DR020	IN														
HEX (dR) 1000 nmol scale	MD-FL010-DR100	IN														
HEX (dT) 40 nmol scale	MD-FL010-DT004	IN														
HEX (dT) 200 nmol scale	MD-FL010-DT020	IN														
HEX (dT) 1000 nmol scale	MD-FL010-DT100	IN														
5' TET 40 nmol scale	MD-FL012-05004	IN														
5' TET 200 nmol scale	MD-FL012-05020	IN														
5' TET 1000 nmol scale	MD-FL012-05100	IN														
TET (dR) 40 nmol scale	MD-FL012-DR004	IN														
TET (dR) 200 nmol scale	MD-FL012-DR020	IN														
TET (dR) 1000 nmol scale	MD-FL012-DR100	IN														
TET (dT) 40 nmol scale	MD-FL012-DT004	IN														
TET (dT) 200 nmol scale	MD-FL012-DT020	IN														
TET (dT) 1000 nmol scale	MD-FL012-DT100	IN														
3' JOE 40 nmol scale	MD-FL014-03004	IN														
3' JOE 200 nmol scale	MD-FL014-03020	IN														
3' JOE 1000 nmol scale	MD-FL014-03100	IN														
5' JOE 40 nmol scale	MD-FL014-05004	IN														
5' JOE 200 nmol scale	MD-FL014-05020	IN														
5' JOE 1000 nmol scale	MD-FL014-05100	IN														
JOE (dR) 40 nmol scale	MD-FL014-DR004	IN														
JOE (dR) 200 nmol scale	MD-FL014-DR020	IN														
JOE (dR) 1000 nmol scale	MD-FL014-DR100	IN														
JOE (dT) 40 nmol scale	MD-FL014-DT004	IN														
JOE (dT) 200 nmol scale	MD-FL014-DT020	IN														
JOE (dT) 1000 nmol scale	MD-FL014-DT100	IN														

Description	Reference	Purifications									Specifications					
		SePOP	RP- Cartridge Gold™	RP-HPLC	IEX-HPLC	PAGE	DUAL HPLC (RP+RP)	DUAL HPLC (RP+HPLC)	Ultra Pure Gold™	Position			Ab <sub>max</sub> (nm)	Em <sub>max</sub> (nm)	MW (g.mol <sup>-1</sup> )	
										5'	IN	3'				
5' ROX 200 nmol scale	MD-FL016-05020										✓	✓	✓	575	602	695.75
5' ROX 1000 nmol scale	MD-FL016-05100										✓	✓	✓	575	602	695.75
ROX (dR) 40 nmol scale	MD-FL016-DR004										✓	✓	✓	575	602	812
ROX (dR) 200 nmol scale	MD-FL016-DR020										✓	✓	✓	575	602	812
ROX (dR) 1000 nmol scale	MD-FL016-DR100										✓	✓	✓	575	602	812
ROX (dT) 40 nmol scale	MD-FL016-DT004										✓	✓	✓	575	602	975.01
ROX (dT) 200 nmol scale	MD-FL016-DT020										✓	✓	✓	575	602	975.01
ROX (dT) 1000 nmol scale	MD-FL016-DT100										✓	✓	✓	575	602	975.01
3' Rhodamine-6G 40 nmol scale	MD-FL020-03004											✓	✓	524	550	649.68
3' Rhodamine-6G 200 nmol scale	MD-FL020-03020											✓	✓	524	550	649.68
3' Rhodamine-6G 1000 nmol scale	MD-FL020-03100											✓	✓	524	550	649.68
5' Rhodamine-6G 40 nmol scale	MD-FL020-05004											✓	✓	524	550	649.68
5' Rhodamine-6G 200 nmol scale	MD-FL020-05020											✓	✓	524	550	649.68
5' Rhodamine-6G 1000 nmol scale	MD-FL020-05100											✓	✓	524	550	649.68
Rhodamine-6G (dR) 40 nmol scale	MD-FL020-DR004											✓	✓	524	550	753.8
Rhodamine-6G (dR) 200 nmol scale	MD-FL020-DR020											✓	✓	524	550	753.8
Rhodamine-6G (dR) 1000 nmol scale	MD-FL020-DR100											✓	✓	524	550	753.8
Rhodamine-6G (dT) 40 nmol scale	MD-FL020-DT004											✓	✓	524	550	870.6
Rhodamine-6G (dT) 200 nmol scale	MD-FL020-DT020											✓	✓	524	550	870.6
Rhodamine-6G (dT) 1000 nmol scale	MD-FL020-DT100											✓	✓	524	550	870.6
3' Texas Red® 40 nmol scale	MD-FL030-03004												✓	583	603	881.01
3' Texas Red® 200 nmol scale	MD-FL030-03020												✓	583	603	881.01
3' Texas Red® 1000 nmol scale	MD-FL030-03100												✓	583	603	881.01
5' Texas Red® 40 nmol scale	MD-FL030-05004												✓	583	603	881.01
5' Texas Red® 200 nmol scale	MD-FL030-05020												✓	583	603	881.01
5' Texas Red® 1000 nmol scale	MD-FL030-05100												✓	583	603	881.01
Texas Red® (dR) 40 nmol scale	MD-FL030-DR004												✓	583	603	996.52
Texas Red® (dR) 200 nmol scale	MD-FL030-DR020												✓	583	603	996.52
Texas Red® (dR) 1000 nmol scale	MD-FL030-DR100												✓	583	603	996.52
Texas Red® (dT) 40 nmol scale	MD-FL030-DT004												✓	583	603	1160.26
Texas Red® (dT) 200 nmol scale	MD-FL030-DT020												✓	583	603	1160.26
Texas Red® (dT) 1000 nmol scale	MD-FL030-DT100												✓	583	603	1160.26
5' Yakima Yellow® 40 nmol scale	MD-FL040-05004												✓	530.5	549	718.33
5' Yakima Yellow® 200 nmol scale	MD-FL040-05020												✓	530.5	549	718.33
5' Yakima Yellow® 1000 nmol scale	MD-FL040-05100												✓	530.5	549	718.33
5' Yakima Yellow® - 2.5 µmol scale	MD-FL040-05M02												✓	530.5	549	718.33
5' Yakima Yellow® - 5 µmol scale	MD-FL040-05M05												✓	530.5	549	718.33
5' Yakima Yellow® - 10 µmol scale	MD-FL040-05M10												✓	530.5	549	718.33
5' Yakima Yellow® - 20 µmol scale	MD-FL040-05M20												✓	530.5	549	718.33
5' Dragonfly Orange™ 40 nmol scale	MD-FL045-05004												✓	554	576	755.62
5' Dragonfly Orange™ 200 nmol scale	MD-FL045-05020												✓	554	576	755.62
5' Dragonfly Orange™ 1000 nmol scale	MD-FL045-05100												✓	554	576	755.62
3' Cy® 3 40 nmol scale	MD-FL050-03004												✓	547	563	811.79
3' Cy® 3 200 nmol scale	MD-FL050-03020												✓	547	563	811.79
3' Cy® 3 1000 nmol scale	MD-FL050-03100												✓	547	563	811.79
5' Cy® 3 40 nmol scale	MD-FL050-05004												✓	547	563	811.79
5' Cy® 3 200 nmol scale	MD-FL050-05020												✓	547	563	811.79
5' Cy® 3 1000 nmol scale	MD-FL050-05100												✓	547	563	811.79
Cy® 3 (dT) 40 nmol scale	MD-FL050-DT004												✓	547	563	1070.16
Cy® 3 (dT) 200 nmol scale	MD-FL050-DT020												✓	547	563	1070.16
Cy® 3 (dT) 1000 nmol scale	MD-FL050-DT100												✓	547	563	1070.16
3' Cy® 3.5 40 nmol scale	MD-FL053-03004												✓	591	604	1079.15
3' Cy® 3.5 200 nmol scale	MD-FL053-03020												✓	591	604	1079.15
3' Cy® 3.5 1000 nmol scale	MD-FL053-03100												✓	591	604	1079.15
5' Cy® 3.5 200 nmol scale	MD-FL053-05020												✓	591	604	607.7
5' Cy® 3.5 1000 nmol scale	MD-FL053-05100												✓	591	604	607.7
3' Cy5® 40 nmol scale	MD-FL055-03004												✓	646	662	838.03
3' Cy5® 200 nmol scale	MD-FL055-03020												✓	646	662	838.03
3' Cy5® 1000 nmol scale	MD-FL055-03100												✓	646	662	838.03
5' Cy5® 40 nmol scale	MD-FL055-05004												✓	646	662	838.03
5' Cy5® 200 nmol scale	MD-FL055-05020												✓	646	662	838.03
5' Cy5® 1000 nmol scale	MD-FL055-05100												✓	646	662	838.03
Cy5® (dT) 40 nmol scale	MD-FL055-DT004												✓	646	662	1096.2
Cy5® (dT) 200 nmol scale	MD-FL055-DT020												✓	646	662	1096.2
Cy5® (dT) 1000 nmol scale	MD-FL055-DT100												✓	646	662	1096.2
3' Cy® 5.5 40 nmol scale	MD-FL057-03004												✓	685	706	633.74
3' Cy® 5.5 200 nmol scale	MD-FL057-03020												✓	685	706	633.74
3' Cy® 5.5 1000 nmol scale	MD-FL057-03100												✓	685	706	633.74
5' Cy® 5.5 40 nmol scale	MD-FL057-05004												✓	685	706	633.74
5' Cy® 5.5 200 nmol scale	MD-FL057-05020												✓	685	706	633.74
5' Cy® 5.5 1000 nmol scale	MD-FL057-05100												✓	685	706	633.74
5' Alexa® 350 40 nmol scale	MD-FL060-35005004												✓	346	442	474.42
5' Alexa® 350 200 nmol scale	MD-FL060-35005020												✓	346	442	474.42
5' Alexa® 350 1000 nmol scale	MD-FL060-35005100												✓	346	442	474.42
5' Alexa® 430 40 nmol scale	MD-FL060-43005004												✓	434	541	664.8
5' Alexa® 430 200 nmol scale	MD-FL060-43005020												✓	434	541	664.8
5' Alexa® 430 1000 nmol scale	MD-FL060-43005100												✓	434	541	664.8
5' Alexa® 488 40 nmol scale	MD-FL060-48805004												✓	495	519	707.48
5' Alexa® 488 200 nmol scale	MD-FL060-48805020												✓	495	519	707.48
5' Alexa® 488 1000 nmol scale	MD-FL060-48805100												✓	495	519	707.48
5' Alexa® 514 40 nmol scale	MD-FL060-51405004												✓	517	542	
5' Alexa® 514 200 nmol scale	MD-FL060-51405020												✓	517	542	
5' Alexa® 514 1000 nmol scale	MD-FL060-51405100												✓	517	542	
5' Alexa® 532 40 nmol scale	MD-FL060-53205004												✓	532	554	785.07
5' Alexa® 532 200 nmol scale	MD-FL060-53205020												✓	532	554	785.07
5' Alexa® 532 1000 nmol scale	MD-FL060-53205100												✓	532	554	785.07
5' Alexa® 546 40 nmol scale	MD-FL060-54605004												✓	556	573	1120.1
5' Alexa® 546 200 nmol scale	MD-FL060-54605020												✓	556	573	1120.1
5' Alexa® 546 1000 nmol scale	MD-FL060-54605100												✓	556	573	1120.1
5' Alexa® 555 40 nmol scale	MD-FL060-55505004												✓	555	565	1154
5' Alexa® 555 200 nmol scale	MD-FL060-55505020												✓	555	565	1154
5' Alexa® 555 1000 nmol scale	MD-FL060-55505100												✓	555	565	1154
5' Alexa® 568 40 nmol scale	MD-FL060-56805004												✓	578	603	855.87
5' Alexa® 568 200 nmol scale	MD-FL060-56805020												✓	578	603	855.87
5' Alexa® 568 1000 nmol scale	MD-FL060-56805100												✓	578	603	855.87
5' Alexa® 594 40 nmol scale	MD-FL060-59405004												✓	590	617	1080
5' Alexa® 594 200 nmol scale	MD-FL060-59405020												✓	590	617	1080
5' Alexa® 594 1000 nmol scale	MD-FL060-59405100												✓	590	617	1080
5' Alexa® 610 40 nmol scale	MD-FL060-61005004												✓	612	628	1031.33
5' Alexa® 610 200 nmol scale	MD-FL060-61005020												✓	612	628	1031.33
5' Alexa® 610 1000 nmol scale	MD-FL060-61005100												✓	612	628	1031.33
5' Alexa® 633 40 nmol scale	MD-FL060-63305004		</													



Description	Reference	Purifications							Specifications					
		SePOP	RP- Cartridge Gold™	RP-HPLC	IEX-HPLC	PAGE	DUAL HPLC (RP+RP)	DUAL HPLC (RP+HPLC)	Ultra Pure Gold™	Position 5' IN 3'	Post Synthesis Coupling*	Ab <sub>max</sub> (nm)	Em <sub>max</sub> (nm)	MW (g.mol <sup>-1</sup> )
5' ATTO 594 40 nmol scale	MD-FL090-59405004											601	627	968.57
5' ATTO 594 200 nmol scale	MD-FL090-59405020											601	627	968.57
5' ATTO 594 1000 nmol scale	MD-FL090-59405100											601	627	968.57
5' ATTO 610 40 nmol scale	MD-FL090-61005004											615	634	552.67
5' ATTO 610 200 nmol scale	MD-FL090-61005020											615	634	552.67
5' ATTO 610 1000 nmol scale	MD-FL090-61005100											615	634	552.67
5' ATTO 620 40 nmol scale	MD-FL090-62005004											619	643	673.07
5' ATTO 620 200 nmol scale	MD-FL090-62005020											619	643	673.07
5' ATTO 620 1000 nmol scale	MD-FL090-62005100											619	643	673.07
5' ATTO 633 40 nmol scale	MD-FL090-63305004											629	653	713.92
5' ATTO 633 200 nmol scale	MD-FL090-63305020											629	653	713.92
5' ATTO 633 1000 nmol scale	MD-FL090-63305100											629	653	713.92
5' ATTO 637 40 nmol scale	MD-FL090-63705004											635	659	838
5' ATTO 637 200 nmol scale	MD-FL090-63705020											635	659	838
5' ATTO 637 1000 nmol scale	MD-FL090-63705100											635	659	838
5' ATTO 647N 40 nmol scale	MD-FL090-64705004											644	669	807.62
5' ATTO 647N 200 nmol scale	MD-FL090-64705020											644	669	807.62
5' ATTO 647N 1000 nmol scale	MD-FL090-64705100											644	669	807.62
5' ATTO 655 40 nmol scale	MD-FL090-65505004											663	684	690
5' ATTO 655 200 nmol scale	MD-FL090-65505020											663	684	690
5' ATTO 655 1000 nmol scale	MD-FL090-65505100											663	684	690
5' ATTO 680 40 nmol scale	MD-FL090-68005004											680	700	687.07
5' ATTO 680 200 nmol scale	MD-FL090-68005020											680	700	687.07
5' ATTO 680 1000 nmol scale	MD-FL090-68005100											680	700	687.07
5' ATTO 700 40 nmol scale	MD-FL090-70005004											700	719	727.16
5' ATTO 700 200 nmol scale	MD-FL090-70005020											700	719	727.16
5' ATTO 700 1000 nmol scale	MD-FL090-70005100											700	719	727.16
5' ATTO 725 40 nmol scale	MD-FL090-72505004											729	752	
5' ATTO 725 200 nmol scale	MD-FL090-72505020											729	752	
5' ATTO 725 1000 nmol scale	MD-FL090-72505100											729	752	
5' ATTO 740 40 nmol scale	MD-FL090-74005004											740	764	786.2
5' ATTO 740 200 nmol scale	MD-FL090-74005020											740	764	786.2
5' ATTO 740 1000 nmol scale	MD-FL090-74005100											740	764	786.2
5' TAMRA 40 nmol scale	MD-FL100-05004											565	580	623.6
5' TAMRA 200 nmol scale	MD-FL100-05020											565	580	623.6
5' TAMRA 1000 nmol scale	MD-FL100-05100											565	580	623.6
<b>Modifiers</b>														<b>KB2D12</b>
5' Amino Modifier C3 40 nmol scale	MD-MF003-05004	IN												137.08
5' Amino Modifier C3 200 nmol scale	MD-MF003-05020	IN												137.08
5' Amino Modifier C3 1000 nmol scale	MD-MF003-05100	IN												137.08
3' Amino Modifier C6 40 nmol scale	MD-MF006-03004	IN												209.18
3' Amino Modifier C6 200 nmol scale	MD-MF006-03020	IN												209.18
3' Amino Modifier C6 1000 nmol scale	MD-MF006-03100	IN												209.18
3' Amino Modifier C6 2.5 µmol scale	MD-MF006-03M02	IN												209.18
3' Amino Modifier C6 5 µmol scale	MD-MF006-03M05	IN												209.18
3' Amino Modifier C6 10 µmol scale	MD-MF006-03M10	IN												209.18
5' Amino Modifier C6 40 nmol scale	MD-MF006-05004	IN												209.18
5' Amino Modifier C6 200 nmol scale	MD-MF006-05020	IN												209.18
5' Amino Modifier C6 1000 nmol scale	MD-MF006-05100	IN												209.18
5' Amino Modifier C6 2.5 µmol scale	MD-MF006-05M02	IN												209.18
5' Amino Modifier C6 5 µmol scale	MD-MF006-05M05	IN												209.18
5' Amino Modifier C6 10 µmol scale	MD-MF006-05M10	IN												209.18
5' Amino Modifier C6 20 µmol scale	MD-MF006-05M20	IN												209.18
Amino Modifier C6 (dT) 40 nmol scale	MD-MF006-DT004	IN												458.41
Amino Modifier C6 (dT) 200 nmol scale	MD-MF006-DT020	IN												458.41
Amino Modifier C6 (dT) 1000 nmol scale	MD-MF006-DT100	IN												458.41
5' Amino Modifier C12 40 nmol scale	MD-MF012-05004	IN												263.32
5' Amino Modifier C12 200 nmol scale	MD-MF012-05020	IN												263.32
5' Amino Modifier C12 1000 nmol scale	MD-MF012-05100	IN												263.32
Internal Amino Modifier Propargyl-dU 40 nmol scale	MD-MF030-IN004	IN												344.22
Internal Amino Modifier Propargyl-dU 200 nmol scale	MD-MF030-IN020	IN												344.22
Internal Amino Modifier Propargyl-dU 1000 nmol scale	MD-MF030-IN100	IN												344.22
Amine (dR) 40 nmol scale	MD-MF040-DR004	IN												295.27
Amine (dR) 200 nmol scale	MD-MF040-DR020	IN												295.27
Amine (dR) 1000 nmol scale	MD-MF040-DR100	IN												295.27
3' Thiol C3 40 nmol scale	MD-MF053-03004	IN												154.12
3' Thiol C3 200 nmol scale	MD-MF053-03020	IN												154.12
3' Thiol C3 1000 nmol scale	MD-MF053-03100	IN												154.12
3' Thiol C6 40 nmol scale	MD-MF056-03004	IN												196.2
3' Thiol C6 200 nmol scale	MD-MF056-03020	IN												196.2
3' Thiol C6 1000 nmol scale	MD-MF056-03100	IN												196.2
3' Thiol C6 2.5 µmol scale	MD-MF056-03M02	IN												196.2
3' Thiol C6 5 µmol scale	MD-MF056-03M05	IN												196.2
3' Thiol C6 10 µmol scale	MD-MF056-03M10	IN												196.2
5' Thiol C6 40 nmol scale	MD-MF056-05004	IN												196.2
5' Thiol C6 200 nmol scale	MD-MF056-05020	IN												196.2
5' Thiol C6 1000 nmol scale	MD-MF056-05100	IN												196.2
5' Thiol C6 2.5 µmol scale	MD-MF056-05M02	IN												196.2
5' Thiol C6 5 µmol scale	MD-MF056-05M05	IN												196.2
5' Thiol C6 10 µmol scale	MD-MF056-05M10	IN												196.2
5' Thiol C6 20 µmol scale	MD-MF056-05M20	IN												196.2
3' Thiophosphate 40 nmol scale	MD-MF060-03004	IN												96.04
3' Thiophosphate 200 nmol scale	MD-MF060-03020	IN												96.04
3' Thiophosphate 1000 nmol scale	MD-MF060-03100	IN												96.04
5' Thiophosphate 40 nmol scale	MD-MF060-05004	IN												96.04
5' Thiophosphate 200 nmol scale	MD-MF060-05020	IN												96.04
5' Thiophosphate 1000 nmol scale	MD-MF060-05100	IN												96.04
5' Triphosphate NEW 40 nmol scale	MD-MF070-05004	IN												
5' Triphosphate NEW 200 nmol scale	MD-MF070-05020	IN												
5' Triphosphate NEW 1000 nmol scale	MD-MF070-05100	IN												
<b>Non natural bases</b>														<b>KB2E17</b>
3'-dinosine 40 nmol scale	MD-NB001-03004	IN												314.19
3'-dinosine 200 nmol scale	MD-NB001-03020	IN												314.19
3'-dinosine 1000 nmol scale	MD-NB001-03100	IN												314.19
dinosine 40 nmol scale	MD-NB001-IN004	IN												314.19
dinosine 200 nmol scale	MD-NB001-IN020	IN												314.19
dinosine 1000 nmol scale	MD-NB001-IN100	IN												314.19
5-Nitroindole 40 nmol scale	MD-NB010-IN004	IN												340.23
5-Nitroindole 200 nmol scale	MD-NB010-IN020	IN												340.23
5-Nitroindole 1000 nmol scale	MD-NB010-IN100	IN												340.23
dP (pyrimidine hybrid) 40 nmol scale	MD-NB030-IN004	IN												330.23
dP (pyrimidine hybrid) 200 nmol scale	MD-NB030-IN020	IN												330.23
dP (pyrimidine hybrid) 1000 nmol scale	MD-NB030-IN100	IN												330.23
dK (purine Analogue) 40 nmol scale	MD-NB035-IN004	IN												358.25
dK (purine Analogue) 200 nmol scale	MD-NB035-IN020	IN												358.25
dK (purine Analogue) 1000 nmol scale	MD-NB035-IN100	IN												358.25

■ Required Purification\*   
■ Compatible Purification\*   
■ Best Purification\*   
■ IN Included Purification  
- Not Applicable

\* Post-synthesis modifications may yield 50 % less than the Minimum Guaranteed Yields



Description	Reference	Purifications								Specifications				
		SePOP	RP- Cartridge Gold™	RP-HPLC	IEX-HPLC	PAGE	DUAL HPLC (RP+RP)	DUAL HPLC (RP+IEX)	Ultra Pure Gold™	Position 5' IN 3'	Post Synthesis Coupling*	Ab <sub>max</sub> (nm)	Em <sub>max</sub> (nm)	MW (g.mol <sup>-1</sup> )
5-Me-iso-dC 40 nmol scale	MD-NB052-IN004	IN								✓	✓			303.21
5-Me-iso-dC 200 nmol scale	MD-NB052-IN020	IN								✓	✓			303.21
5-Me-iso-dC 1000 nmol scale	MD-NB052-IN100	IN								✓	✓			303.21
isoGuanosine (iso-dG) 40 nmol scale	MD-NB054-IN004	IN								✓	✓			329.21
isoGuanosine (iso-dG) 200 nmol scale	MD-NB054-IN020	IN								✓	✓			329.21
isoGuanosine (iso-dG) 1000 nmol scale	MD-NB054-IN100	IN								✓	✓			329.21
3' dUracil 40 nmol scale	MD-NB068-03004	IN									✓			290.17
3' dUracil 200 nmol scale	MD-NB068-03020	IN									✓			290.17
3' dUracil 1000 nmol scale	MD-NB068-03100	IN									✓			290.17
dUracil 40 nmol scale	MD-NB068-IN004	IN								✓	✓			290.17
dUracil 200 nmol scale	MD-NB068-IN020	IN								✓	✓			290.17
dUracil 1000 nmol scale	MD-NB068-IN100	IN								✓	✓			290.17
3' dA Blocked 40 nmol scale	MD-NB070-03004	IN									✓			313.21
3' dA Blocked 200 nmol scale	MD-NB070-03020	IN									✓			313.21
3' dA Blocked 1000 nmol scale	MD-NB070-03100	IN									✓			313.21
3' dC Blocked 40 nmol scale	MD-NB072-03004	IN									✓			289.18
3' dC Blocked 200 nmol scale	MD-NB072-03020	IN									✓			289.18
3' dC Blocked 1000 nmol scale	MD-NB072-03100	IN									✓			289.18
3' dG Blocked 40 nmol scale	MD-NB074-03004	IN									✓			329.21
3' dG Blocked 200 nmol scale	MD-NB074-03020	IN									✓			329.21
3' dG Blocked 1000 nmol scale	MD-NB074-03100	IN									✓			329.21
3' dT Blocked 40 nmol scale	MD-NB076-03004	IN									✓			304.21
3' dT Blocked 200 nmol scale	MD-NB076-03020	IN									✓			304.21
3' dT Blocked 1000 nmol scale	MD-NB076-03100	IN									✓			304.21
2', 3' ddC Blocked 40 nmol scale	MD-NB080-03004	IN								✓	✓			273.18
2', 3' ddC Blocked 200 nmol scale	MD-NB080-03020	IN								✓	✓			273.18
2', 3' ddC Blocked 1000 nmol scale	MD-NB080-03100	IN								✓	✓			273.18
3'-3' Terminus Linkage dT (inverted 3' base) 40 nmol scale	MD-NB085-03004	IN									✓			304.2
3'-3' Terminus Linkage dT (inverted 3' base) 200 nmol scale	MD-NB085-03020	IN									✓			304.2
3'-3' Terminus Linkage dT (inverted 3' base) 1000 nmol scale	MD-NB085-03100	IN									✓			304.2
3'-3' Terminus Linkage dA (inverted 3' base) 40 nmol scale	MD-NB085-04004	IN									✓			304.2
3'-3' Terminus Linkage dA (inverted 3' base) 200 nmol scale	MD-NB085-04020	IN									✓			304.2
3'-3' Terminus Linkage dA (inverted 3' base) 1000 nmol scale	MD-NB085-04100	IN									✓			304.2
3'-3' Terminus Linkage dC (inverted 3' base) 40 nmol scale	MD-NB085-05004	IN									✓			304.2
3'-3' Terminus Linkage dC (inverted 3' base) 200 nmol scale	MD-NB085-05020	IN									✓			304.2
3'-3' Terminus Linkage dC (inverted 3' base) 1000 nmol scale	MD-NB085-05100	IN									✓			304.2
3'-3' Terminus Linkage dG (inverted 3' base) 40 nmol scale	MD-NB085-06004	IN									✓			304.2
3'-3' Terminus Linkage dG (inverted 3' base) 200 nmol scale	MD-NB085-06020	IN									✓			304.2
3'-3' Terminus Linkage dG (inverted 3' base) 1000 nmol scale	MD-NB085-06100	IN									✓			304.2
8-Br-dA 40 nmol scale	MD-NB090-IN004	IN								✓	✓			392.11
8-Br-dA 200 nmol scale	MD-NB090-IN020	IN								✓	✓			392.11
8-Br-dA 1000 nmol scale	MD-NB090-IN100	IN								✓	✓			392.11
5-Br-dC 40 nmol scale	MD-NB092-IN004	IN								✓	✓			368.08
5-Br-dC 200 nmol scale	MD-NB092-IN020	IN								✓	✓			368.08
5-Br-dC 1000 nmol scale	MD-NB092-IN100	IN								✓	✓			368.08
8-Br-dG 40 nmol scale	MD-NB094-IN004	IN								✓	✓			408.1
8-Br-dG 200 nmol scale	MD-NB094-IN020	IN								✓	✓			408.1
8-Br-dG 1000 nmol scale	MD-NB094-IN100	IN								✓	✓			408.1
3' (5-Br-dU) 40 nmol scale	MD-NB098-03004	IN									✓			369.07
3' (5-Br-dU) 200 nmol scale	MD-NB098-03020	IN									✓			369.07
3' (5-Br-dU) 1000 nmol scale	MD-NB098-03100	IN									✓			369.07
5-Br-dU 40 nmol scale	MD-NB098-IN004	IN								✓	✓			369.07
5-Br-dU 200 nmol scale	MD-NB098-IN020	IN								✓	✓			369.07
5-Br-dU 1000 nmol scale	MD-NB098-IN100	IN								✓	✓			369.07
5,6-dihydro-dU 40 nmol scale	MD-NB108-IN004	IN								✓	✓			292.19
5,6-dihydro-dU 200 nmol scale	MD-NB108-IN020	IN								✓	✓			292.19
5,6-dihydro-dU 1000 nmol scale	MD-NB108-IN100	IN								✓	✓			292.19
7-deaza-dA 40 nmol scale	MD-NB110-IN004	IN									✓			312.2
7-deaza-dA 200 nmol scale	MD-NB110-IN020	IN									✓			312.2
7-deaza-dA 1000 nmol scale	MD-NB110-IN100	IN									✓			312.2
7-deaza-dG 40 nmol scale	MD-NB114-IN004	IN									✓			328.22
7-deaza-dG 200 nmol scale	MD-NB114-IN020	IN									✓			328.22
7-deaza-dG 1000 nmol scale	MD-NB114-IN100	IN									✓			328.22
8-Oxo-dA 40 nmol scale	MD-NB120-IN004	IN									✓			329.21
8-Oxo-dA 200 nmol scale	MD-NB120-IN020	IN									✓			329.21
8-Oxo-dA 1000 nmol scale	MD-NB120-IN100	IN									✓			329.21
8-Oxo-dG 40 nmol scale	MD-NB124-IN004	IN									✓			345.21
8-Oxo-dG 200 nmol scale	MD-NB124-IN020	IN									✓			345.21
8-Oxo-dG 1000 nmol scale	MD-NB124-IN100	IN									✓			345.21
06-Me-dG 40 nmol scale	MD-NB134-IN004	IN									✓			343.24
06-Me-dG 200 nmol scale	MD-NB134-IN020	IN									✓			343.24
06-Me-dG 1000 nmol scale	MD-NB134-IN100	IN									✓			343.24
C5-propyne dC 40 nmol scale	MD-NB142-IN004	IN									✓			327.23
C5-propyne dC 200 nmol scale	MD-NB142-IN020	IN									✓			327.23
C5-propyne dC 1000 nmol scale	MD-NB142-IN100	IN									✓			327.23
C5-propyne dU 40 nmol scale	MD-NB148-IN004	IN									✓			328.22
C5-propyne dU 200 nmol scale	MD-NB148-IN020	IN									✓			328.22
C5-propyne dU 1000 nmol scale	MD-NB148-IN100	IN									✓			328.22
3' (5-Me-dC) 40 nmol scale	MD-NB152-03004	IN									✓			303.21
3' (5-Me-dC) 200 nmol scale	MD-NB152-03020	IN									✓			303.21
3' (5-Me-dC) 1000 nmol scale	MD-NB152-03100	IN									✓			303.21
5-Me-dC 40 nmol scale	MD-NB152-IN004	IN									✓			303.21
5-Me-dC 200 nmol scale	MD-NB152-IN020	IN									✓			303.21
5-Me-dC 1000 nmol scale	MD-NB152-IN100	IN									✓			303.21
2-Amino-dA NEW 40 nmol scale	MD-NB160-IN004	IN									✓			328.22
2-Amino-dA NEW 200 nmol scale	MD-NB160-IN020	IN									✓			328.22
2-Amino-dA NEW 1000 nmol scale	MD-NB160-IN100	IN									✓			328.22
2-Aminopurine (DNA) 40 nmol scale	MD-NB170-IN004	IN									✓			313.21
2-Aminopurine (DNA) 200 nmol scale	MD-NB170-IN020	IN									✓			313.21
2-Aminopurine (DNA) 1000 nmol scale	MD-NB170-IN100	IN									✓			313.21
N4-Et-dC 40 nmol scale	MD-NB180-IN004	IN									✓			317.42
N4-Et-dC 200 nmol scale	MD-NB180-IN020	IN									✓			317.42
N4-Et-dC 1000 nmol scale	MD-NB180-IN100	IN									✓			317.42
AP-dC 40 nmol scale	MD-NB190-IN004	IN									✓			
AP-dC 200 nmol scale	MD-NB190-IN020	IN									✓			
AP-dC 1000 nmol scale	MD-NB190-IN100	IN									✓			
2' Fluoro RNA A 40 nmol scale	MD-NB201-IN004	IN									✓			307.17
2' Fluoro RNA A 200 nmol scale	MD-NB201-IN020	IN									✓			307.17
2' Fluoro RNA A 1000 nmol scale	MD-NB201-IN100	IN									✓			307.17
2' Fluoro RNA C 40 nmol scale	MD-NB202-IN004	IN									✓			307.17
2' Fluoro RNA C 200 nmol scale	MD-NB202-IN020	IN									✓			307.17
2' Fluoro RNA C 1000 nmol scale	MD-NB202-IN100	IN									✓			307.17
2' Fluoro RNA G 40 nmol scale	MD-NB203-IN004	IN									✓			308.16
2' Fluoro RNA G 200 nmol scale	MD-NB203-IN020	IN									✓			308.16
2' Fluoro RNA G 1000 nmol scale	MD-NB203-IN100	IN									✓			308.16
2' Fluoro RNA U 40 nmol scale	MD-NB208-IN004	IN									✓			308.16

■ Required Purification\* 
 ■ Compatible Purification\* 
 ■ Best Purification\* 
 IN Included Purification  
 - Not Applicable

\* Post-synthesis modifications may yield 50 % less than the Minimum Guaranteed Yields





Description	Reference	Purifications							Specifications						
		SePOP	RP- Cartridge Gold™	RP-HPLC	IEX-HPLC	PAGE	DUAL HPLC (RP+RP)	DUAL HPLC (RP+HEX)	Ultra Pure Gold™	Position			Ab <sub>max</sub> (nm)	Em <sub>max</sub> (nm)	MW (g.mol <sup>-1</sup> )
									5'	IN	3'	Post Synthesis Coupling*			
5' Biotin TEG 1000 nmol scale	MD-BT001-05100														569.61
5' Biotin TEG 2,5 µmol scale	MD-BT001-05M02														569.61
5' Biotin TEG 5 µmol scale	MD-BT001-05M05														569.61
5' Biotin TEG 10 µmol scale	MD-BT001-05M10														569.61
5' Biotin TEG 20 µmol scale	MD-BT001-05M20														569.61
Biotin (dR) 40 nmol scale	MD-BT001-DR004	IN													521.57
Biotin (dR) 200 nmol scale	MD-BT001-DR020	IN													521.57
Biotin (dR) 1000 nmol scale	MD-BT001-DR100	IN													521.57
Biotin (dR) 2,5 µmol scale	MD-BT001-DRM02	IN													521.57
Biotin (dR) 5 µmol scale	MD-BT001-DRM05	IN													521.57
Biotin (dR) 10 µmol scale	MD-BT001-DRM10	IN													521.57
Biotin (dT) 40 nmol scale	MD-BT001-DT004	IN													684.7
Biotin (dT) 200 nmol scale	MD-BT001-DT020	IN													684.7
Biotin (dT) 1000 nmol scale	MD-BT001-DT100	IN													684.7
5' PC Biotin 40 nmol scale	MD-BT010-05004														597.62
5' PC Biotin 200 nmol scale	MD-BT010-05020														597.62
5' PC Biotin 1000 nmol scale	MD-BT010-05100														597.62
5' Biotin 40 nmol scale	MD-BT020-05004	IN													405.45
5' Biotin 200 nmol scale	MD-BT020-05020	IN													405.45
5' Biotin 1000 nmol scale	MD-BT020-05100	IN													405.45
3' Biotin C3 40 nmol scale	MD-BT030-03004	IN													
3' Biotin C3 200 nmol scale	MD-BT030-03020	IN													
3' Biotin C3 1000 nmol scale	MD-BT030-03100	IN													
<b>NGS Modifications</b>															
NGS 5' Biotin-TEG	MD-BT002-05001														569.61
NGS 5' Phosphate	MD-PH002-05001														79.98

■ Required Purification\*   
■ Compatible Purification\*   
■ Best Purification\*   
IN Included Purification  
■ - Not Applicable

\* Post-synthesis modifications may yield 50 % less than the Minimum Guaranteed Yields