

Item Number	Modifications for custom oligonucleotides Conjugates
MD-CN001-03004	3' Digoxigenin 40 nmol
MD-CN001-03020	3' Digoxigenin 200 nmol
MD-CN001-03100	3' Digoxigenin 1000 nmol
MD-CN001-05004	5'-Digoxigenin 40 nmol
MD-CN001-05020	5' Digoxigenin 200 nmol
MD-CN001-05100	5' Digoxigenin 1000 nmol
MD-CN001-DR004	Digoxigenin (dR) 40 nmol
MD-CN001-DR020	Digoxigenin (dR) 200 nmol
MD-CN001-DR100	Digoxigenin (dR) 1000 nmol
MD-CN001-DT004	Digoxigenin (dT) 40 nmol
MD-CN001-DT020	Digoxigenin (dT) 200 nmol
MD-CN001-DT100	Digoxigenin (dT) 1000 nmol
MD-CN010-03004	3' TEG-Dinitrophenol 40 nmol
MD-CN010-03020	3' TEG-Dinitrophenol 200 nmol
MD-CN010-03100	3' TEG-Dinitrophenol 1000 nmol
MD-CN010-05004	5' TEG-Dinitrophenol 40 nmol
MD-CN010-05020	5' TEG-Dinitrophenol 200 nmol
MD-CN010-05100	5' TEG-Dinitrophenol 1000 nmol
MD-CN010-DR004	Dinitrophenol (dR) 40 nmol
MD-CN010-DR020	Dinitrophenol (dR) 200 nmol
MD-CN010-DR100	Dinitrophenol (dR) 1000 nmol
MD-CN010-DT004	Dinitrophenol (dT) 40 nmol
MD-CN010-DT020	Dinitrophenol (dT) 200 nmol
MD-CN010-DT100	Dinitrophenol (dT) 1000 nmol
MD-CN020-05004	5' Acrylamide (Acrydite®) 40 nmol
MD-CN020-05020	5' Acrylamide (Acrydite®) 200 nmol
MD-CN020-05100	5' Acrylamide (Acrydite®) 1000 nmol
MD-CN030-03004	3' Glyceryl 40 nmol
MD-CN030-03020	3' Glyceryl 200 nmol
MD-CN030-03100	3' Glyceryl 1000 nmol
MD-CN040-03004	3' TEG-Cholesteryl 40 nmol
MD-CN040-03020	3' TEG-Cholesteryl 200 nmol
MD-CN040-03100	3' TEG-Cholesteryl 1000 nmol
MD-CN040-05004	5' TEG-Cholesteryl 40 nmol
MD-CN040-05020	5' TEG-Cholesteryl 200 nmol
MD-CN040-05100	5' TEG-Cholesteryl 1000 nmol
MD-CN040-DR004	Cholesteryl (dR) 40 nmol
MD-CN040-DR020	Cholesteryl (dR) 200 nmol
MD-CN040-DR100	Cholesteryl (dR) 1000 nmol
MD-CN050-03004	3' MeO-CI-Acridine 40 nmol
MD-CN050-03020	3' MeO-CI-Acridine 200 nmol
MD-CN050-03100	3' MeO-CI-Acridine 1000 nmol
MD-CN050-05004	5' MeO-CI-Acridine 40 nmol
MD-CN050-05020	5' MeO-CI-Acridine 200 nmol
MD-CN050-05100	5' MeO-CI-Acridine 1000 nmol
MD-CN050-IN004	Internal MeO-CI-Acridine 40 nmol
MD-CN050-IN020	Internal MeO-CI-Acridine 200 nmol
MD-CN050-IN100	Internal MeO-CI-Acridine 1000 nmol
MD-CN062-05004	5' C2 Psoralen 40 nmol
MD-CN062-05020	5' C2 Psoralen 200 nmol

MD-CN062-05100	5' C2 Psoralen 1000 nmol
MD-CN066-05004	5' C6 Psoralen 40 nmol
MD-CN066-05020	5' C6 Psoralen 200 nmol
MD-CN066-05100	5' C6 Psoralen 1000 nmol
MD-CN070-DT004	Carboxy (dT) 40 nmol
MD-CN070-DT020	Carboxy (dT) 200 nmol
MD-CN070-DT100	Carboxy (dT) 1000 nmol
MD-CN090-05100	5' AP Conjugation 1000 nmol
MD-CN090-05M02	5' AP Conjugation 2,5 µmol
MD-CN092-05100	5' HRP Conjugation 1000 nmol
MD-CN092-05M02	5' HRP Conjugation 2,5 µmol
MD-CN094-05100	5' BSA Conjugation 1000 nmol
MD-CN094-05M02	5' BSA Conjugation 2,5 µmol
MD-CN096-05100	5' SBP Conjugation 1000 nmol
MD-CN096-05M02	5' SBP Conjugation 2,5 µmol
MD-CN099-05M02	5' Peptide Conjugation 2,5 µmol
MD-CN100-03100	3' Oligo Affinity Support (PS) 1000 nmol
MD-CN110-03100	3' Oligo Affinity Support (CPG) 1000 nmol