

Item Number	Modifications for custom oligonucleotides Modifiers
MD-MF003-05004	5' Amino Modifier C3 40 nmol
MD-MF003-05020	5' Amino Modifier C3 200 nmol
MD-MF003-05100	5' Amino Modifier C3 1000 nmol
MD-MF006-03004	3' Amino Modifier C6 40 nmol
MD-MF006-03020	3' Amino Modifier C6 200 nmol
MD-MF006-03100	3' Amino Modifier C6 1000 nmol
MD-MF006-05004	5' Amino Modifier C6 40 nmol
MD-MF006-05020	5' Amino Modifier C6 200 nmol
MD-MF006-05100	5' Amino Modifier C6 1000 nmol
MD-MF006-05M02	5' Amino Modifier C6 2,5 µmol
MD-MF006-05M05	5' Amino Modifier C6 5 µmol
MD-MF006-05M10	5' Amino Modifier C6 10 µmol
MD-MF006-DT004	Amino Modifier C6 (dT) 40 nmol
MD-MF006-DT020	Amino Modifier C6 (dT) 200 nmol
MD-MF006-DT100	Amino Modifier C6 (dT) 1000 nmol
MD-MF012-05004	5' Amino Modifier C12 40 nmol
MD-MF012-05020	5' Amino Modifier C12 200 nmol
MD-MF012-05100	5' Amino Modifier C12 1000 nmol
MD-MF020-IN004	Internal Amino Modifier dU 40 nmol
MD-MF020-IN020	Internal Amino Modifier dU 200 nmol
MD-MF020-IN100	Internal Amino Modifier dU 1000 nmol
MD-MF030-IN004	Internal Amino Modifier Propargyl-dU 40 nmol
MD-MF030-IN020	Internal Amino Modifier Propargyl-dU 200 nmol
MD-MF030-IN100	Internal Amino Modifier Propargyl-dU 1000 nmol
MD-MF040-DR004	Amine (dR) 40 nmol
MD-MF040-DR020	Amine (dR) 200 nmol
MD-MF040-DR100	Amine (dR) 1000 nmol
MD-MF053-03004	3' Thiol C3 40 nmol
MD-MF053-03020	3' Thiol C3 200 nmol
MD-MF053-03100	3' Thiol C3 1000 nmol
MD-MF056-03004	3' Thiol C6 40 nmol
MD-MF056-03020	3' Thiol C6 200 nmol
MD-MF056-03100	3' Thiol C6 1000 nmol
MD-MF056-03M02	3' Thiol C6 2,5 µmol
MD-MF056-03M05	3' Thiol C6 5 µmol
MD-MF056-03M10	3' Thiol C6 10 µmol
MD-MF056-05004	5' Thiol C6 40 nmol
MD-MF056-05020	5' Thiol C6 200 nmol
MD-MF056-05100	5' Thiol C6 1000 nmol
MD-MF056-05M02	5' Thiol C6 2,5 µmol
MD-MF056-05M05	5' Thiol C6 5 µmol
MD-MF056-05M10	5' Thiol C6 10 µmol
MD-MF060-03004	3' Thiophosphate 40 nmol
MD-MF060-03020	3' Thiophosphate 200 nmol
MD-MF060-03100	3' Thiophosphate 1000 nmol
MD-MF060-05004	5' Thiophosphate 40 nmol
MD-MF060-05020	5' Thiophosphate 200 nmol
MD-MF060-05100	5' Thiophosphate 1000 nmol
MD-MF070-05004	5' Triphosphate NEW 40 nmol
MD-MF070-05020	5' Triphosphate NEW 200 nmol

MD-MF070-05100

5' Triphosphate NEW 1000 nmol