

Item Number	Modifications for custom oligonucleotides Nonnatural bases
MD-NB001-03004	3'-dInosine 40 nmol
MD-NB001-03020	3'-dInosine 200 nmol
MD-NB001-03100	3'-dInosine 1000 nmol
MD-NB001-IN004	dInosine 40 nmol
MD-NB001-IN020	dInosine 200 nmol
MD-NB001-IN100	dInosine 1000 nmol
MD-NB010-03004	5-Nitroindole 40 nmol
MD-NB010-03020	5-Nitroindole 200 nmol
MD-NB010-03100	5-Nitroindole 1000 nmol
MD-NB010-IN004	5-Nitroindole 40 nmol
MD-NB010-IN020	5-Nitroindole 200 nmol
MD-NB010-IN100	5-Nitroindole 1000 nmol
MD-NB020-03004	3-Nitropyrrole 40 nmol
MD-NB020-03020	3-Nitropyrrole 200 nmol
MD-NB020-03100	3-Nitropyrrole 1000 nmol
MD-NB020-IN004	3-Nitropyrrole 40 nmol
MD-NB020-IN020	3-Nitropyrrole 200 nmol
MD-NB020-IN100	3-Nitropyrrole 1000 nmol
MD-NB030-IN004	dP (pyrimidine hybrid) 40 nmol
MD-NB030-IN020	dP (pyrimidine hybrid) 200 nmol
MD-NB030-IN100	dP (pyrimidine hybrid) 1000 nmol
MD-NB035-IN004	dK (purine Analogue) 40 nmol
MD-NB035-IN020	dK (purine Analogue) 200 nmol
MD-NB035-IN100	dK (purine Analogue) 1000 nmol
MD-NB052-IN004	isoCytosine (iso-dC) 40 nmol
MD-NB052-IN020	isoCytosine (iso-dC) 200 nmol
MD-NB052-IN100	isoCytosine (iso-dC) 1000 nmol
MD-NB054-IN004	isoGuanosine (iso-dG) 40 nmol
MD-NB054-IN020	isoGuanosine (iso-dG) 200 nmol
MD-NB054-IN100	isoGuanosine (iso-dG) 1000 nmol
MD-NB068-03004	3' dUracil 40 nmol
MD-NB068-03020	3' dUracil 200 nmol
MD-NB068-03100	3' dUracil 1000 nmol
MD-NB068-IN004	dUracil 40 nmol
MD-NB068-IN020	dUracil 200 nmol
MD-NB068-IN100	dUracil 1000 nmol
MD-NB070-03004	3' dA Blocked 40 nmol
MD-NB070-03020	3' dA Blocked 200 nmol
MD-NB070-03100	3' dA Blocked 1000 nmol
MD-NB072-03004	3' dC Blocked 40 nmol
MD-NB072-03020	3' dC Blocked 200 nmol
MD-NB072-03100	3' dC Blocked 1000 nmol
MD-NB074-03004	3' dG Blocked 40 nmol
MD-NB074-03020	3' dG Blocked 200 nmol
MD-NB074-03100	3' dG Blocked 1000 nmol
MD-NB076-03004	3' dT Blocked 40 nmol
MD-NB076-03020	3' dT Blocked 200 nmol
MD-NB076-03100	3' dT Blocked 1000 nmol
MD-NB080-03004	2', 3' ddC Blocked 40 nmol
MD-NB080-03020	2', 3' ddC Blocked 200 nmol

MD-NB080-03100	2', 3' ddC Blocked 1000 nmol
MD-NB085-03004	3'-3' Terminus Linkage (inverted 3' base) 40 nmol
MD-NB085-03020	3'-3' Terminus Linkage (inverted 3' base) 200 nmol
MD-NB085-03100	3'-3' Terminus Linkage (inverted 3' base) 1000 nmol
MD-NB090-IN004	8-Br-dA 40 nmol
MD-NB090-IN020	8-Br-dA 200 nmol
MD-NB090-IN100	8-Br-dA 1000 nmol
MD-NB092-IN004	5-Br-dC 40 nmol
MD-NB092-IN020	5-Br-dC 200 nmol
MD-NB092-IN100	5-Br-dC 1000 nmol
MD-NB094-IN004	8-Br-dG 40 nmol
MD-NB094-IN020	8-Br-dG 200 nmol
MD-NB094-IN100	8-Br-dG 1000 nmol
MD-NB098-03004	3' (5-Br-dU) 40 nmol
MD-NB098-03020	3' (5-Br-dU) 200 nmol
MD-NB098-03100	3' (5-Br-dU) 1000 nmol
MD-NB098-IN004	5-Br-dU 40 nmol
MD-NB098-IN020	5-Br-dU 200 nmol
MD-NB098-IN100	5-Br-dU 1000 nmol
MD-NB108-IN004	5.6-dihydro-dU 40 nmol
MD-NB108-IN020	5.6-dihydro-dU 200 nmol
MD-NB108-IN100	5.6-dihydro-dU 1000 nmol
MD-NB110-IN004	7-deaza-dA 40 nmol
MD-NB110-IN020	7-deaza-dA 200 nmol
MD-NB110-IN100	7-deaza-dA 1000 nmol
MD-NB114-IN004	7-deaza-dG 40 nmol
MD-NB114-IN020	7-deaza-dG 200 nmol
MD-NB114-IN100	7-deaza-dG 1000 nmol
MD-NB120-IN004	8-Oxo-dA 40 nmol
MD-NB120-IN020	8-Oxo-dA 200 nmol
MD-NB120-IN100	8-Oxo-dA 1000 nmol
MD-NB124-IN004	8-Oxo-dG 40 nmol
MD-NB124-IN020	8-Oxo-dG 200 nmol
MD-NB124-IN100	8-Oxo-dG 1000 nmol
MD-NB134-IN004	O6-Me-dG 40 nmol
MD-NB134-IN020	O6-Me-dG 200 nmol
MD-NB134-IN100	O6-Me-dG 1000 nmol
MD-NB142-IN004	C5-propyne dC 40 nmol
MD-NB142-IN020	C5-propyne dC 200 nmol
MD-NB142-IN100	C5-propyne dC 1000 nmol
MD-NB148-IN004	C5-propyne dU 40 nmol
MD-NB148-IN020	C5-propyne dU 200 nmol
MD-NB148-IN100	C5-propyne dU 1000 nmol
MD-NB152-03004	3' (5-Me-dC) 40 nmol
MD-NB152-03020	3' (5-Me-dC) 200 nmol
MD-NB152-03100	3' (5-Me-dC) 1000 nmol
MD-NB152-IN004	5-Me-dC 40 nmol
MD-NB152-IN020	5-Me-dC 200 nmol
MD-NB152-IN100	5-Me-dC 1000 nmol
MD-NB160-IN004	2-Amino-dA NEW 40 nmol
MD-NB160-IN020	2-Amino-dA NEW 200 nmol

MD-NB160-IN100	2-Amino-dA NEW 1000 nmol
MD-NB170-IN004	2-Aminopurine (DNA) NEW
MD-NB170-IN020	2-Aminopurine (DNA) NEW
MD-NB170-IN100	2-Aminopurine (DNA) NEW
MD-NB180-IN004	N4-Et-dC 40 nmol
MD-NB180-IN020	N4-Et-dC 200 nmol
MD-NB180-IN100	N4-Et-dC 1000 nmol
MD-NB190-IN004	AP-dC 40 nmol
MD-NB190-IN020	AP-dC 200 nmol
MD-NB190-IN100	AP-dC 1000 nmol
MD-NB202-IN004	2' Fluoro RNA C 40 nmol
MD-NB202-IN020	2' Fluoro RNA C 200 nmol
MD-NB202-IN100	2' Fluoro RNA C 1000 nmol
MD-NB208-IN004	2' Fluoro RNA U 40 nmol
MD-NB208-IN020	2' Fluoro RNA U 200 nmol
MD-NB208-IN100	2' Fluoro RNA U 1000 nmol
MD-NB210-IN004	2' O-Me RNA Inosine 40 nmol
MD-NB210-IN020	2' O-Me RNA Inosine 200 nmol
MD-NB210-IN100	2' O-Me RNA Inosine 1000 nmol
MD-NB222-IN004	2' O-Me RNA-5-Me-C 200 nmol
MD-NB222-IN020	2' O-Me RNA-5-Me-C 200 nmol
MD-NB222-IN100	2' O-Me RNA-5-Me-C 1000 nmol
MD-NB232-IN004	2' O-Me RNA Propynyl-C 40 nmol
MD-NB232-IN020	2' O-Me RNA Propynyl-C 200 nmol
MD-NB232-IN100	2' O-Me RNA Propynyl-C 1000 nmol
MD-NB238-IN004	2' O-Me RNA Propynyl-U 40 nmol
MD-NB238-IN020	2' O-Me RNA Propynyl-U 200 nmol
MD-NB238-IN100	2' O-Me RNA Propynyl-U 1000 nmol