

Item Number	Modifications for custom oligonucleotides Quenchers
MD-QU100-03004	3' TAMRA 40 nmol
MD-QU100-03020	3' TAMRA 200 nmol
MD-QU100-03100	3' TAMRA 1000 nmol
MD-QU100-DT004	TAMRA (dT) 40 nmol
MD-QU100-DT020	TAMRA (dT) 200 nmol
MD-QU100-DT100	TAMRA (dT) 1000 nmol
MD-QU200-03004	3' DABCYL 40 nmol
MD-QU200-03020	3' DABCYL 200 nmol
MD-QU200-03100	3' DABCYL 1000 nmol
MD-QU200-05004	5' DABCYL 40 nmol
MD-QU200-05020	5' DABCYL 200 nmol
MD-QU200-05100	5' DABCYL 1000 nmol
MD-QU200-DT004	DABCYL (dT) 40 nmol
MD-QU200-DT020	DABCYL (dT) 200 nmol
MD-QU200-DT100	DABCYL (dT) 1000 nmol
MD-QU300-03004	3' Eclipse® Dark Quencher 40 nmol
MD-QU300-03020	3' Eclipse® Dark Quencher 200 nmol
MD-QU300-03100	3' Eclipse® Dark Quencher 1000 nmol
MD-QU300-05004	5' Eclipse® Dark Quencher 40 nmol
MD-QU300-05020	5' Eclipse® Dark Quencher 200 nmol
MD-QU300-05100	5' Eclipse® Dark Quencher 1000 nmol
MD-QU400-03004	3' Deep Dark Quencher I 40 nmol
MD-QU400-03020	3' Deep Dark Quencher I 200 nmol
MD-QU400-03100	3' Deep Dark Quencher I 1000 nmol
MD-QU400-DR004	Deep Dark Quencher I (dR) 40 nmol
MD-QU400-DR020	Deep Dark Quencher I (dR) 200 nmol
MD-QU400-DR100	Deep Dark Quencher I (dR) 1000 nmol
MD-QU400-DT004	Deep Dark Quencher I (dT) 40 nmol
MD-QU400-DT020	Deep Dark Quencher I (dT) 200 nmol
MD-QU400-DT100	Deep Dark Quencher I (dT) 1000 nmol
MD-QU500-03004	3' Deep Dark Quencher II 40 nmol
MD-QU500-03020	3' Deep Dark Quencher II 200 nmol
MD-QU500-03100	3' Deep Dark Quencher II 1000 nmol
MD-QU500-05004	5' Deep Dark Quencher II 40 nmol
MD-QU500-05020	5' Deep Dark Quencher II 200 nmol
MD-QU500-05100	5' Deep Dark Quencher II 1000 nmol
MD-QU500-DR004	Deep Dark Quencher II (dR) 40 nmol
MD-QU500-DR020	Deep Dark Quencher II (dR) 200 nmol
MD-QU500-DR100	Deep Dark Quencher II (dR) 1000 nmol
MD-QU600-03004	3' BHQ-1™ / BHQ-2™ / BHQ-3™ 40 nmol
MD-QU600-03020	3' BHQ-1™ / BHQ-2™ / BHQ-3™ 200 nmol
MD-QU600-03100	3' BHQ-1™ / BHQ-2™ / BHQ-3™ 1000 nmol
MD-QU600-05004	5' BHQ-1™ / BHQ-2™ 40 nmol
MD-QU600-05020	5' BHQ-1™ / BHQ-2™ 200 nmol
MD-QU600-05100	5' BHQ-1™ / BHQ-2™ 1000 nmol
MD-QU600-DR004	Internal BHQ-2# (dT) 40 nmol
MD-QU600-DR020	Internal BHQ-2# (dT) 200 nmol
MD-QU600-DR100	Internal BHQ-2# (dT) 1000 nmol
MD-QU600-DT004	BHQ-1™/BHQ-2™(dT) 40 nmol
MD-QU600-DT020	BHQ-1™/BHQ-2™ (dT) 200 nmol

MD-QU600-DT100	BHQ-1 TM /BHQ-2 TM (dT) 1000 nmol
MD-QU700-03004	3' QXLTM 490/570/610/670/680 NEW 40 nmol
MD-QU700-03020	3' QXLTM 490/570/610/670/680 NEW 200 nmol
MD-QU700-03100	3' QXLTM 490/570/610/670/680 NEW 1000 nmol