

Item Number	Modifications for custom oligonucleotides Biotins
MD-BT001-03004	3' Biotin TEG 40 nmol
MD-BT001-03020	3' Biotin TEG 200 nmol
MD-BT001-03100	3' Biotin TEG 1000 nmol
MD-BT001-05004	5' Biotin TEG 40 nmol
MD-BT001-05020	5' Biotin TEG 200 nmol
MD-BT001-05100	5' Biotin TEG 1000 nmol
MD-BT001-05M02	5' Biotin TEG 2,5 µmol
MD-BT001-05M05	5' Biotin TEG 5 µmol
MD-BT001-05M10	5' Biotin TEG 10 µmol
MD-BT001-DR004	Biotin (dR) 40 nmol
MD-BT001-DR020	Biotin (dR) 200 nmol
MD-BT001-DR100	Biotin (dR) 1000 nmol
MD-BT001-DRM02	Biotin (dR) 2,5 µmol
MD-BT001-DRM05	Biotin (dR) 5 µmol
MD-BT001-DRM10	Biotin (dR) 10 µmol
MD-BT001-DT004	Biotin (dT) 40 nmol
MD-BT001-DT020	Biotin (dT) 200 nmol
MD-BT001-DT100	Biotin (dT) 1000 nmol
MD-BT010-05004	5' PC Biotin 40 nmol
MD-BT010-05020	5' PC Biotin 200 nmol
MD-BT010-05100	5' PC Biotin 1000 nmol
MD-BT010-IN004	5' PC Biotin 40 nmol
MD-BT010-IN020	5' PC Biotin 200 nmol
MD-BT010-IN100	5' PC Biotin 1000 nmol
MD-BT020-05004	5' Biotin New 40 nmol
MD-BT020-05020	5' Biotin New 200 nmol
MD-BT020-05100	5' Biotin New 1000 nmol