

Item Number	Modifications for custom oligonucleotides Fluorophores
MD-FL001-03004	3' 6-FAM 40 nmol
MD-FL001-03020	3' 6-FAM 200 nmol
MD-FL001-03100	3' 6-FAM 1000 nmol
MD-FL001-05004	5' 6-FAM 40 nmol
MD-FL001-05020	5' 6-FAM 200 nmol
MD-FL001-05100	5' 6-FAM 1000 nmol
MD-FL001-05M02	5' 6-FAM 2,5 µmol
MD-FL001-05M05	5' 6-FAM 5 µmol
MD-FL001-05M10	5' 6-FAM 10 µmol
MD-FL001-DR004	6-FAM (dR) 40 nmol
MD-FL001-DR020	6-FAM (dR) 200 nmol
MD-FL001-DR100	6-FAM (dR) 1000 nmol
MD-FL001-DT004	6-FAM (dT) 40 nmol
MD-FL001-DT020	6-FAM (dT) 200 nmol
MD-FL001-DT100	6-FAM (dT) 1000 nmol
MD-FL005-03004	3' Fluorescein 40 nmol
MD-FL005-03020	3' Fluorescein 200 nmol
MD-FL005-03100	3' Fluorescein 1000 nmol
MD-FL005-05004	5' Fluorescein 40 nmol
MD-FL005-05020	5' Fluorescein 200 nmol
MD-FL005-05100	5' Fluorescein 1000 nmol
MD-FL010-05004	5' HEX 40 nmol
MD-FL010-05020	5' HEX 200 nmol
MD-FL010-05100	5' HEX 1000 nmol
MD-FL010-DR004	HEX (dR) 40 nmol
MD-FL010-DR020	HEX (dR) 200 nmol
MD-FL010-DR100	HEX (dR) 1000 nmol
MD-FL010-DT004	HEX (dT) 40 nmol
MD-FL010-DT020	HEX (dT) 200 nmol
MD-FL010-DT100	HEX (dT) 1000 nmol
MD-FL012-05004	5' TET 40 nmol
MD-FL012-05020	5' TET 200 nmol
MD-FL012-05100	5' TET 1000 nmol
MD-FL012-DR004	TET (dR) 40 nmol
MD-FL012-DR020	TET (dR) 200 nmol
MD-FL012-DR100	TET (dR) 1000 nmol
MD-FL012-DT004	TET (dT) 40 nmol
MD-FL012-DT020	TET (dT) 200 nmol
MD-FL012-DT100	TET (dT) 1000 nmol
MD-FL014-03004	3' JOE 40 nmol
MD-FL014-03020	3' JOE 200 nmol
MD-FL014-03100	3' JOE 1000 nmol
MD-FL014-05004	5' JOE 40 nmol
MD-FL014-05020	5' JOE 200 nmol
MD-FL014-05100	5' JOE 1000 nmol
MD-FL014-DR004	JOE (dR) 40 nmol
MD-FL014-DR020	JOE (dR) 200 nmol
MD-FL014-DR100	JOE (dR) 1000 nmol
MD-FL014-DT004	JOE (dT) 40 nmol
MD-FL014-DT020	JOE (dT) 200 nmol

MD-FL014-DT100	JOE (dT) 1000 nmol
MD-FL016-03004	3' ROX 40 nmol
MD-FL016-03020	3' ROX 200 nmol
MD-FL016-03100	3' ROX 1000 nmol
MD-FL016-05004	5' ROX 40 nmol
MD-FL016-05020	5' ROX 200 nmol
MD-FL016-05100	5' ROX 1000 nmol
MD-FL016-DR004	ROX (dR) 40 nmol
MD-FL016-DR020	ROX (dR) 200 nmol
MD-FL016-DR100	ROX (dR) 1000 nmol
MD-FL016-DT004	ROX (dT) 40 nmol
MD-FL016-DT020	ROX (dT) 200 nmol
MD-FL016-DT100	ROX (dT) 1000 nmol
MD-FL020-03004	3' Rhodamine-6G 40 nmol
MD-FL020-03020	3' Rhodamine-6G 200 nmol
MD-FL020-03100	3' Rhodamine-6G 1000 nmol
MD-FL020-05004	5' Rhodamine-6G 40 nmol
MD-FL020-05020	5' Rhodamine-6G 200 nmol
MD-FL020-05100	5' Rhodamine-6G 1000 nmol
MD-FL020-DR004	Rhodamine-6G (dR) 40 nmol
MD-FL020-DR020	Rhodamine-6G (dR) 200 nmol
MD-FL020-DR100	Rhodamine-6G (dR) 1000 nmol
MD-FL020-DT004	Rhodamine-6G (dT) 40 nmol
MD-FL020-DT020	Rhodamine-6G (dT) 200 nmol
MD-FL020-DT100	Rhodamine-6G (dT) 1000 nmol
MD-FL030-03004	3' Texas Red® 40 nmol
MD-FL030-03020	3' Texas Red® 200 nmol
MD-FL030-03100	3' Texas Red® 1000 nmol
MD-FL030-05004	5' Texas Red® 40 nmol
MD-FL030-05020	5' Texas Red® 200 nmol
MD-FL030-05100	5' Texas Red® 1000 nmol
MD-FL030-DR004	Texas Red® (dR) 40 nmol
MD-FL030-DR020	Texas Red® (dR) 200 nmol
MD-FL030-DR100	Texas Red® (dR) 1000 nmol
MD-FL030-DT004	Texas Red® (dT) 40 nmol
MD-FL030-DT020	Texas Red® (dT) 200 nmol
MD-FL030-DT100	Texas Red® (dT) 1000 nmol
MD-FL040-05004	5' Yakima Yellow® 40 nmol
MD-FL040-05020	5' Yakima Yellow® 200 nmol
MD-FL040-05100	5' Yakima Yellow® 1000 nmol
MD-FL045-05004	5' Dragonfly Orange™ 40 nmol
MD-FL045-05020	5' Dragonfly Orange™ 200 nmol
MD-FL045-05100	5' Dragonfly Orange™ 1000 nmol
MD-FL050-03004	3' Cy® 3 40 nmol
MD-FL050-03020	3' Cy® 3 200 nmol
MD-FL050-03100	3' Cy® 3 1000 nmol
MD-FL050-05004	5' Cy® 3 40 nmol
MD-FL050-05020	5' Cy® 3 200 nmol
MD-FL050-05100	5' Cy® 3 1000 nmol
MD-FL050-DT004	Cy® 3 (dT) 40 nmol
MD-FL050-DT020	Cy® 3 (dT) 200 nmol

MD-FL050-DT100	Cy@ 3 (dT) 1000 nmol
MD-FL053-03004	3' Cy@ 3.5 40 nmol
MD-FL053-03020	3' Cy@ 3.5 200 nmol
MD-FL053-03100	3' Cy@ 3.5 1000 nmol
MD-FL053-05004	5' Cy@ 3.5 40 nmol
MD-FL053-05020	5' Cy@ 3.5 200 nmol
MD-FL053-05100	5' Cy@ 3.5 1000 nmol
MD-FL053-DT004	Cy@ 3.5 (dT) 40 nmol
MD-FL053-DT020	Cy@ 3.5 (dT) 200 nmol
MD-FL053-DT100	Cy@ 3.5 (dT) 1000 nmol
MD-FL055-03004	3' Cy5@ 40 nmol
MD-FL055-03020	3' Cy5@ 200 nmol
MD-FL055-03100	3' Cy5@ 1000 nmol
MD-FL055-05004	5' Cy5@ 40 nmol
MD-FL055-05020	5' Cy5@ 200 nmol
MD-FL055-05100	5' Cy5@ 1000 nmol
MD-FL055-DT004	Cy5@ (dT) 40 nmol
MD-FL055-DT020	Cy5@ (dT) 200 nmol
MD-FL055-DT100	Cy5@ (dT) 1000 nmol
MD-FL057-03004	3' Cy@ 5.5 40 nmol
MD-FL057-03020	3' Cy@ 5.5 200 nmol
MD-FL057-03100	3' Cy@ 5.5 1000 nmol
MD-FL057-05004	5' Cy@ 5.5 40 nmol
MD-FL057-05020	5' Cy@ 5.5 200 nmol
MD-FL057-05100	5' Cy@ 5.5 1000 nmol
MD-FL057-DT004	Cy@ 5.5 (dT) 40 nmol
MD-FL057-DT020	Cy@ 5.5 (dT) 200 nmol
MD-FL057-DT100	Cy@ 5.5 (dT) 1000 nmol
MD-FL060-35005004	5' Alexa@ 350 40 nmol
MD-FL060-35005020	5' Alexa@ 350 200 nmol
MD-FL060-35005100	5' Alexa@ 350 1000 nmol
MD-FL060-43005004	5' Alexa@ 430 40 nmol
MD-FL060-43005020	5' Alexa@ 430 200 nmol
MD-FL060-43005100	5' Alexa@ 430 1000 nmol
MD-FL060-48005004	5' Alexa@ 488 40 nmol
MD-FL060-48805020	5' Alexa@ 488 200 nmol
MD-FL060-48805100	5' Alexa@ 488 1000 nmol
MD-FL060-50005004	5' Alexa@ 500 40 nmol
MD-FL060-50005020	5' Alexa@ 500 200 nmol
MD-FL060-50005100	5' Alexa@ 500 1000 nmol
MD-FL060-51405004	5' Alexa@ 514 40 nmol
MD-FL060-51405020	5' Alexa@ 514 200 nmol
MD-FL060-51405100	5' Alexa@ 514 1000 nmol
MD-FL060-53205004	5' Alexa@ 532 40 nmol
MD-FL060-53205020	5' Alexa@ 532 200 nmol
MD-FL060-53205100	5' Alexa@ 532 1000 nmol
MD-FL060-54605004	5' Alexa@ 546 40 nmol
MD-FL060-54605020	5' Alexa@ 546 200 nmol
MD-FL060-54605100	5' Alexa@ 546 1000 nmol
MD-FL060-55505004	5' Alexa@ 555 40 nmol
MD-FL060-55505020	5' Alexa@ 555 200 nmol

MD-FL060-55505100	5' Alexa® 555 1000 nmol
MD-FL060-56805004	5' Alexa® 568 40 nmol
MD-FL060-56805020	5' Alexa® 568 200 nmol
MD-FL060-56805100	5' Alexa® 568 1000 nmol
MD-FL060-59405004	5' Alexa® 594 40 nmol
MD-FL060-59405020	5' Alexa® 594 200 nmol
MD-FL060-59405100	5' Alexa® 594 1000 nmol
MD-FL060-61005004	5' Alexa® 610 40 nmol
MD-FL060-61005020	5' Alexa® 610 200 nmol
MD-FL060-61005100	5' Alexa® 610 1000 nmol
MD-FL060-63305004	5' Alexa® 633 40 nmol
MD-FL060-63305020	5' Alexa® 633 200 nmol
MD-FL060-63305100	5' Alexa® 633 1000 nmol
MD-FL060-64705004	5' Alexa® 647 40 nmol
MD-FL060-64705020	5' Alexa® 647 200 nmol
MD-FL060-64705100	5' Alexa® 647 1000 nmol
MD-FL060-66005004	5' Alexa® 660 40 nmol
MD-FL060-66005020	5' Alexa® 660 200 nmol
MD-FL060-66005100	5' Alexa® 660 1000 nmol
MD-FL060-68005004	5' Alexa® 680 40 nmol
MD-FL060-68005020	5' Alexa® 680 200 nmol
MD-FL060-68005100	5' Alexa® 680 1000 nmol
MD-FL060-70005004	5' Alexa® 700 40 nmol
MD-FL060-70005020	5' Alexa® 700 200 nmol
MD-FL060-70005100	5' Alexa® 700 1000 nmol
MD-FL060-75005004	5' Alexa® 750 40 nmol
MD-FL060-75005020	5' Alexa® 750 200 nmol
MD-FL060-75005100	5' Alexa® 750 1000 nmol
MD-FL065-40505004	5' HiLyte Fluor™ 405 NEW 40 nmol
MD-FL065-40505020	5' HiLyte Fluor™ 405 NEW 200 nmol
MD-FL065-40505100	5' HiLyte Fluor™ 405 NEW 1000 nmol
MD-FL065-48805004	5' HiLyte Fluor™ 488 NEW 40 nmol
MD-FL065-48805020	5' HiLyte Fluor™ 488 NEW 200 nmol
MD-FL065-48805100	5' HiLyte Fluor™ 488 NEW 1000 nmol
MD-FL065-55505004	5' HiLyte Fluor™ 555 NEW 40 nmol
MD-FL065-55505020	5' HiLyte Fluor™ 555 NEW 200 nmol
MD-FL065-55505100	5' HiLyte Fluor™ 555 NEW 1000 nmol
MD-FL065-59405004	5' HiLyte Fluor™ 594 NEW 40 nmol
MD-FL065-59405020	5' HiLyte Fluor™ 594 NEW 200 nmol
MD-FL065-59405100	5' HiLyte Fluor™ 594 NEW 1000 nmol
MD-FL065-64705004	5' HiLyte Fluor™ 647 NEW 40 nmol
MD-FL065-64705020	5' HiLyte Fluor™ 647 NEW 200 nmol
MD-FL065-64705100	5' HiLyte Fluor™ 647 NEW 1000 nmol
MD-FL065-68005004	5' HiLyte Fluor™ 680 NEW 40 nmol
MD-FL065-68005020	5' HiLyte Fluor™ 680 NEW 200 nmol
MD-FL065-68005100	5' HiLyte Fluor™ 680 NEW 1000 nmol
MD-FL065-75005004	5' HiLyte Fluor™ 750 NEW 40 nmol
MD-FL065-75005020	5' HiLyte Fluor™ 750 NEW 200 nmol
MD-FL065-75005100	5' HiLyte Fluor™ 750 NEW 1000 nmol
MD-FL070-CBL03004	3' Cascade Blue® 40 nmol
MD-FL070-CBL03020	3' Cascade Blue® 200 nmol

MD-FL070-CBL03100	3' Cascade Blue® 1000 nmol
MD-FL070-CBL05004	5' Cascade Blue® 40 nmol
MD-FL070-CBL05020	5' Cascade Blue® 200 nmol
MD-FL070-CBL05100	5' Cascade Blue® 1000 nmol
MD-FL070-CBLDT004	Cascade Blue® (dT-40 nmol)
MD-FL070-CBLDT020	Cascade Blue® (dT-200 nmol)
MD-FL070-CBLDT100	Cascade Blue® (dT-1000 nmol)
MD-FL070-MBL05004	5' Marina Blue® 40 nmol
MD-FL070-MBL05020	5' Marina Blue® 200 nmol
MD-FL070-MBL05100	5' Marina Blue® 1000 nmol
MD-FL070-OGR05004	5' Oregon Green® 488 40 nmol
MD-FL070-OGR05020	5' Oregon Green® 488 200 nmol
MD-FL070-OGR05100	5' Oregon Green® 488 1000 nmol
MD-FL070-OGX05004	5' Oregon Green® 488 X 40 nmol
MD-FL070-OGX05020	5' Oregon Green® 488 X 200 nmol
MD-FL070-OGX05100	5' Oregon Green® 488 X 1000 nmol
MD-FL070-PBL05004	5' Pacific Blue® 40 nmol
MD-FL070-PBL05020	5' Pacific Blue® 200 nmol
MD-FL070-PBL05100	5' Pacific Blue® 1000 nmol
MD-FL080-68105004	5' DYOMICS 681 40 nmol
MD-FL080-68105020	5' DYOMICS 681 200 nmol
MD-FL080-68105100	5' DYOMICS 681 1000 nmol
MD-FL080-78105004	5' DYOMICS 781 40 nmol
MD-FL080-78105020	5' DYOMICS 781 200 nmol
MD-FL080-78105100	5' DYOMICS 781 1000 nmol
MD-FL080-78205004	5' DYOMICS 782 40 nmol
MD-FL080-78205020	5' DYOMICS 782 200 nmol
MD-FL080-78205100	5' DYOMICS 782 1000 nmol
MD-FL080-80005004	5' DYOMICS 800 40 nmol
MD-FL080-80005020	5' DYOMICS 800 200 nmol
MD-FL080-80005100	5' DYOMICS 800 1000 nmol
MD-FL085-53005004	5' BODIPY® 530/550 40 nmol
MD-FL085-53005020	5' BODIPY® 530/550 200 nmol
MD-FL085-53005100	5' BODIPY® 530/550 1000 nmol
MD-FL085-FL05004	5' BODIPY® FL 40 nmol
MD-FL085-FL05020	5' BODIPY® FL 200 nmol
MD-FL085-FL05100	5' BODIPY® FL 1000 nmol
MD-FL085-TMR05004	5' BODIPY® TMR 40 nmol
MD-FL085-TMR05020	5' BODIPY® TMR 200 nmol
MD-FL085-TMR05100	5' BODIPY® TMR 1000 nmol
MD-FL085-TR05004	5' BODIPY® TR 40 nmol
MD-FL085-TR05020	5' BODIPY® TR 200 nmol
MD-FL085-TR05100	5' BODIPY® TR 1000 nmol
MD-FL090-39005004	5' ATTO 390 40 nmol
MD-FL090-39005020	5' ATTO 390 200 nmol
MD-FL090-39005100	5' ATTO 390 1000 nmol
MD-FL090-42505004	5' ATTO 425 40 nmol
MD-FL090-42505020	5' ATTO 425 200 nmol
MD-FL090-42505100	5' ATTO 425 1000 nmol
MD-FL090-46505004	5' ATTO 465 40 nmol
MD-FL090-46505020	5' ATTO 465 200 nmol

MD-FL090-46505100	5' ATTO 465 1000 nmol
MD-FL090-48805004	5' ATTO 488 40 nmol
MD-FL090-48805020	5' ATTO 488 200 nmol
MD-FL090-48805100	5' ATTO 488 1000 nmol
MD-FL090-49505004	5' ATTO 495 40 nmol
MD-FL090-49505020	5' ATTO 495 200 nmol
MD-FL090-49505100	5' ATTO 495 1000 nmol
MD-FL090-52005004	5' ATTO 520 40 nmol
MD-FL090-52005020	5' ATTO 520 200 nmol
MD-FL090-52005100	5' ATTO 520 1000 nmol
MD-FL090-53205004	5' ATTO 532 40 nmol
MD-FL090-53205020	5' ATTO 532 200 nmol
MD-FL090-53205100	5' ATTO 532 1000 nmol
MD-FL090-55005004	5' ATTO 550 40 nmol
MD-FL090-55005020	5' ATTO 550 200 nmol
MD-FL090-55005100	5' ATTO 550 1000 nmol
MD-FL090-56505004	5' ATTO 565 40 nmol
MD-FL090-56505020	5' ATTO 565 200 nmol
MD-FL090-56505100	5' ATTO 565 1000 nmol
MD-FL090-59005004	5' ATTO 590 40 nmol
MD-FL090-59005020	5' ATTO 590 200 nmol
MD-FL090-59005100	5' ATTO 590 1000 nmol
MD-FL090-59405004	5' ATTO 594 40 nmol
MD-FL090-59405020	5' ATTO 594 200 nmol
MD-FL090-59405100	5' ATTO 594 1000 nmol
MD-FL090-61005004	5' ATTO 610 40 nmol
MD-FL090-61005020	5' ATTO 610 200 nmol
MD-FL090-61005100	5' ATTO 610 1000 nmol
MD-FL090-62005004	5' ATTO 620 40 nmol
MD-FL090-62005020	5' ATTO 620 200 nmol
MD-FL090-62005100	5' ATTO 620 1000 nmol
MD-FL090-63305004	5' ATTO 633 40 nmol
MD-FL090-63305020	5' ATTO 633 200 nmol
MD-FL090-63305100	5' ATTO 633 1000 nmol
MD-FL090-63505004	5' ATTO 635 40 nmol
MD-FL090-63505020	5' ATTO 635 200 nmol
MD-FL090-63505100	5' ATTO 635 1000 nmol
MD-FL090-63705004	5' ATTO 637 40 nmol
MD-FL090-63705020	5' ATTO 637 200 nmol
MD-FL090-63705100	5' ATTO 637 1000 nmol
MD-FL090-64705004	5' ATTO 647N 40 nmol
MD-FL090-64705020	5' ATTO 647N 200 nmol
MD-FL090-64705100	5' ATTO 647N 1000 nmol
MD-FL090-65505004	5' ATTO 655 200 nmol
MD-FL090-65505020	5' ATTO 655 200 nmol
MD-FL090-65505100	5' ATTO 655 1000 nmol
MD-FL090-68005004	5' ATTO 680 200 nmol
MD-FL090-68005020	5' ATTO 680 200 nmol
MD-FL090-68005100	5' ATTO 680 1000 nmol
MD-FL090-70005004	5' ATTO 700 40 nmol
MD-FL090-70005020	5' ATTO 700 200 nmol

MD-FL090-70005100	5' ATTO 700 1000 nmol
MD-FL090-72505004	5' ATTO 725 40 nmol
MD-FL090-72505020	5' ATTO 725 200 nmol
MD-FL090-72505100	5' ATTO 725 1000 nmol
MD-FL090-74005004	5' ATTO 740 40 nmol
MD-FL090-74005020	5' ATTO 740 200 nmol
MD-FL090-74005100	5' ATTO 740 1000 nmol
MD-FL100-05004	5' TAMRA 40 nmol
MD-FL100-05020	5' TAMRA 200 nmol
MD-FL100-05100	5' TAMRA 1000 nmol