

Item Number	2' OMethyl RNA Molecular Beacons
PB-OB201-020	O Beacon 5' FAM + 3' DABCYL 200 nmol
PB-OB210-020	O Beacon 5' HEX + 3' DABCYL 200 nmol
PB-OB212-020	O Beacon 5' TET + 3' DABCYL 200 nmol