

Item Number	WavelengthShifting Molecular Beacons
PB-WB016-020	WS Molecular Beacon ROX / FAM FRET 200 nmol
PB-WB030-020	WS Molecular Beacon Texas Red® / FAM FRET 200 nmol
PB-WB100-020	WS Molecular Beacon TAMRA / FAM FRET 200 nmol