

Item Number	Standard Scorpions® probes
PB-SP001-020	Standard Scorpions® Primers 200 nmol