

Item Number	FRET Scorpions® probes
PB-SP210-020	FRET Scorpions® 5' HEX + 3' DABCYL 200 nmol