

<b>Item Number</b>	<b>Duplex Scorpions probes</b>
PB-SP212-020	Scorpions® 5' TET + 3' DABCYL 200 nmol