

Item Number	Plexor™ Human control gene primer sets
PB-PP014-005	Plexor™ Labelled Primer ACTIN 5' HEX 5 nmol
PB-PP019-005	Plexor™ Labelled Primer GAPDH 5' HEX 5 nmol
PB-PP034-005	Plexor™ Labelled Primer ACTIN 5' Texas Red® 5 nmol
PB-PP039-005	Plexor™ Labelled Primer GAPDH 5' Texas Red® 5 nmol