

<b>Item Number</b>	<b>Plexor™ Labelled Primers</b>
PB-PP001-004	Plexor™ Labelled Primer 5' FAM 40 nmol
PB-PP001-020	Plexor™ Labelled Primer 5' FAM 200 nmol
PB-PP010-004	Plexor™ Labelled Primer 5' HEX 40 nmol
PB-PP010-020	Plexor™ Labelled Primer 5' HEX 200 nmol
PB-PP012-004	Plexor™ Labelled Primer 5' TET 40 nmol
PB-PP012-020	Plexor™ Labelled Primer 5' TET 200 nmol
PB-PP016-004	Plexor™ Labelled Primer 5' ROX 40 nmol
PB-PP016-020	Plexor™ Labelled Primer 5' ROX 200 nmol
PB-PP030-004	Plexor™ Labelled Primer 5' Texas Red® 40 nmol
PB-PP030-020	Plexor™ Labelled Primer 5' Texas Red® 200 nmol
PB-PP055-004	Plexor™ Labelled Primer 5' Cy5® 40 nmol
PB-PP055-020	Plexor™ Labelled Primer 5' Cy5® 200 nmol
PB-PP100-004	Plexor™ Labelled Primer 5' TAMRA 40 nmol
PB-PP100-020	Plexor™ Labelled Primer 5' TAMRA 200 nmol