

Reference	Product name	# Rxn (25 µl)	Assay type	ROX	Format	Template type DNA, cDNA, gDNA RNA 1-Step RNA 2-Step (combine with our first strand kit: RTCTCK-03)	Carry-over prevention C: dUTP + UNG D: dUTP; no UNG (add UNG: RT-0610-03)	Maximal sensitivity S: (dTTP only, no UNG)	Cycling compatibility F: FAST cycling R: Regular cycling	Blue dye	Corresponding sample
<b>ABI PRISM 5700, 7000, 7300, 7700, 7900, STEPONE/PLUS, MASTERCYCLER® EP REALPLEX (WITH ROX PASSIVE REFERENCE)</b>											
RT-SY2X-03+WOUUB	MESA BLUE qPCR MasterMix Plus for SYBR® Assay 7.5 ml	600	SYBR	Y	MasterMix	DNA	D		F + R	Y	RT-SY2X-005+WOUUB
RT-SY2X-03+WOUF	MESA FAST qPCR MasterMix Plus for SYBR® Assay 7.5 ml	600	SYBR	Y	MasterMix	DNA	D		F + R	N	RT-SY2X-005+WOUF
<b>ABI 7500, MX4000®, MX3000P®, MX3005P® (WITH LOW ROX PASSIVE REFERENCE)</b>											
RT-SY2X-03+WOU LRB	MESA BLUE qPCR MasterMix Plus for SYBR® Assay Low ROX w/o UNG 7.5 ml	600	SYBR	Y	MasterMix	DNA	D		F + R	Y	RT-SY2X-005+WOU LRB
RT-SY2X-03+WOU LRF	MESA FAST qPCR MasterMix Plus for SYBR® Assay Low ROX w/o UNG 7.5 ml	600	SYBR	Y	MasterMix	DNA	D		F + R	N	RT-SY2X-005+WOU LRF
<b>ICYCLER IQ®, MY IQ, DNA ENGINE OPTICON® 1, 2, CHROMO 4, MINIOPTICON, CFX96, ROTORGENE® 2000/3000/6000/Q AND QUANTICA®, MX4000®, MX3000P®, MX3005P®, SMARTCYCLER® 1 &amp; 2, LC480 (WITHOUT ROX PASSIVE REFERENCE)</b>											
RT-SY2X-03+NRWOU B	MESA BLUE qPCR MasterMix Plus for SYBR® Assay No ROX 7.5 ml	600	SYBR	N	MasterMix	DNA	D		F + R	Y	RT-SY2X-005+NRWOU B
RT-SY2X-03+NRWOUN F	MESA FAST qPCR MasterMix Plus dTTP for SYBR® Assay No ROX 7.5 ml	600	SYBR	N	MasterMix	DNA	S		F + R	N	RT-SY2X-005NRWOUN F
RT-SY2X-03+NRWOU F	MESA FAST qPCR MasterMix Plus for SYBR® Assay No ROX 7.5 ml	600	SYBR	N	MasterMix	DNA	D		F + R	N	RT-SY2X-005+NRWOU F
<b>ICYCLER IQ®, MY IQ, IQ™5 (WITHOUT ROX BUT WITH FLUORESC EIN)</b>											
RT-SY2X-03+WOU FLB	MESA BLUE qPCR MasterMix Plus for SYBR® Assay w/fluorescein 7.5 ml	600	SYBR	N	MasterMix	DNA	D		F + R	Y	RT-SY2X-005+WOU FLB