

BLUE MasterMixes for SYBR® and Probe Assays

Avoid aberrant results before they occur!

Get more features at no additional cost

For FAST and REGULAR cycling

For SYBR® and Probe Assays

Easy-to-see reagents

Ready-to-use MasterMixes

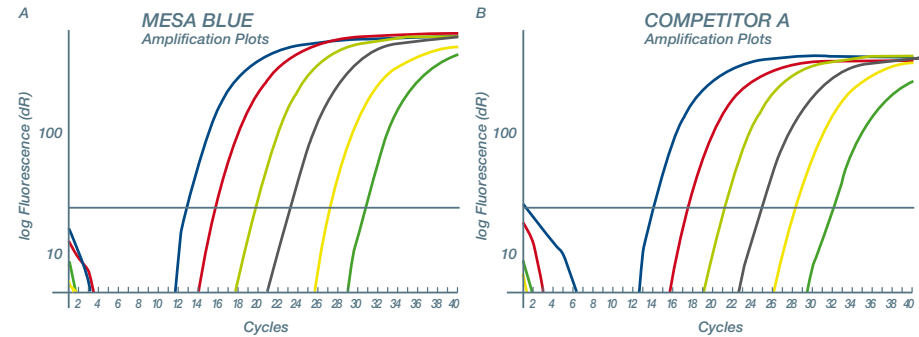


Eurogentec
Experience true partnership

Avoid aberrant results before they occur!

Avoid aberrant results before they occur with Eurogentec's new range of **BLUE MasterMixes** for SYBR® and Probe Assays. The **BLUE MasterMixes** are a unique combination of our highly efficient MeteorTaq together with our new **inert BLUE dye**, which does not affect the qPCR reaction. The **BLUE dye** considerably enhances the contrast between reagent and plasticware, making verification of MasterMix dispensing quick, easy and foolproof.

These ready-to-use, easy-to-see formulated mixes allow you to achieve the most consistent and sensitive results across all different qPCR platforms, in FAST and REGULAR cycling conditions, and avoid aberrant results before they occur.



Increased sensitivity of the MESA BLUE MasterMix Plus No ROX on a MX3000P®. Comparison of the MESA BLUE MasterMix Plus for SYBR® Assay and competitor A (A & B). Experiment was performed using 18S primers and the same cDNA. Eurogentec's new MESA BLUE MasterMix showed superior performance, better sensitivity, higher ΔRn and steeper amplification curves compared to competitor A SYBR® Green reagent.

Get more features at no additional cost

- Visualize dispensing errors at a glance
- Optimized for maximal sensitivity
- High consistency and specificity
- Increased signal intensity on ROX machines
- For FAST and REGULAR protocols

Send us your order

e-mail: order@eurogentec.com
 fax: +32 (0)4 264 07 88

For UK orders

e-mail: order.uk@eurogentec.com
 fax: +44 (0)1794 522 417

Send your sample request

e-mail: info@eurogentec.com
 or contact your local representative

For general information e-mail us at info@eurogentec.com or visit our website www.eurogentec.com

Product	Reference	# RXNs (25 µl)	Sample reference	FAST cycling systems			REGULAR cycling systems					
				ABI Prism® SDS 7900HT ABI StepOne/Plus MasterCycler® ep realplex	ABI Prism® 7500	RotorGene 6000 LC480 SmartCycler® I, II	ABI GeneAmp® 5700 ABI Prism® SDS 7000/7300/7700/7900HT/ MasterCycler® ep realplex StepOne/Plus	ABI Prism® 7500	iCycler iQ® My iQ iQ5	Mx4000® Mx3000P® Mx3005P® with ROX	Mx4000® Mx3000P® Mx3005P® without ROX	RotorGene 2000/3000/6000 MiniOpticon® DNA Engine Opticon® 1,2 Chromo4 Quanta® LC480 SmartCycler® I, II
BLUE MasterMixes for SYBR® Assays												
MESA BLUE qPCR MasterMix Plus for SYBR® Assay, 7.5 ml	RT-SY2X-03+WOUB	600	RT-SY2X-005+WOUB	✓			✓					
MESA BLUE qPCR MasterMix Plus for SYBR® Assay Low ROX, 7.5 ml	RT-SY2X-03+WOUFRB	600	RT-SY2X-005+WOUFRB		✓			✓		✓		
MESA BLUE qPCR MasterMix Plus for SYBR® Assay No ROX, 7.5 ml	RT-SY2X-03+NRWOUB	600	RT-SY2X-005+NRWOUB			✓					✓	✓
MESA BLUE qPCR MasterMix Plus for SYBR® Assay w/fluorescein, 7.5 ml	RT-SY2X-03+WOUFLB	600	RT-SY2X-005+WOUFLB						✓			
BLUE MasterMixes for Probe Assays												
FAST BLUE qPCR MasterMix Plus for Probe Assay, 7.5 ml	RT-QP2X-03+FB	600	RT-QP2X-005+FB	✓			✓					
FAST BLUE qPCR MasterMix Plus for Probe Assay dTTP, 7.5 ml	RT-QP2X-03+WOUNFB	600	RT-QP2X-005+WOUNFB	✓			✓					
FAST BLUE qPCR MasterMix Plus for Probe Assay Low ROX, 7.5 ml	RT-QP2X-03+WOUFRB	600	RT-QP2X-005+WOUFRB		✓			✓		✓		
FAST BLUE qPCR MasterMix Plus for Probe Assay No ROX, 7.5 ml	RT-QP2X-03+NRFB	600	RT-QP2X-005+NRFB			✓			✓		✓	✓