

*qPCR is a
curve...*



Stay ahead!

FAST MasterMixes for SYBR[®] and Probe Assays

Accelerate your research

For SYBR[®] and Probe Assays

For all qPCR platforms

Available in BLUE format

Ready-to-use MasterMixes



Eurogentec
Experience true partnership

Reduce your run times without compromise

Significantly reduce your run times with Eurogentec's new range of FAST MasterMixes for SYBR® and Probe Assays, without compromising your results. Our new FAST MasterMixes deliver accurate, sensitive and consistent results, using a broad range of template amounts, in FAST as well as in REGULAR cycling conditions, on all qPCR platforms.

Designed for ease of use, these kits are also available with our new **inert BLUE dye**, which does not affect the qPCR reaction. The new **FAST BLUE qPCR kits** for Probe Assays considerably enhance the contrast between reagent and plasticware, making verification of MasterMix dispensing quick, easy and foolproof.

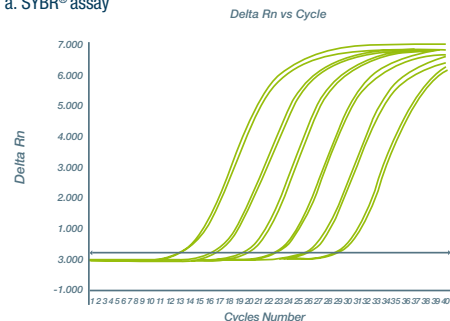
Reduce your run times without compromise

- Tailored for highly sensitive assays
- Optimized for maximum consistency and specificity
- Large dynamic range

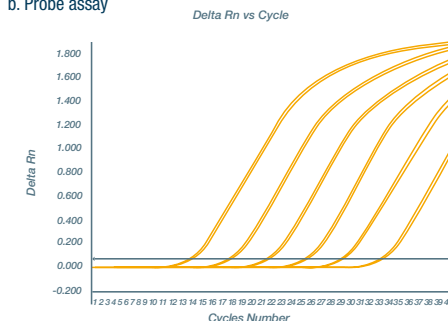
Also available in BLUE format

- New proprietary inert BLUE dye
- Increased signal intensity
- Clear visual dispensing of the reagent

a. SYBR® assay



b. Probe assay



High quality results achieved in 42 minutes with FAST qPCR MasterMixes Low ROX on an ABI Prism® 7500 FAST cycler using ten-times serial dilutions of 18S cDNA. a. MESA FAST qPCR MasterMix Plus Low ROX for SYBR® Assays (E: 99.6 %; R²=0.9998). b. FAST qPCR MasterMix Plus Low ROX for Probe Assays (E: 99.2 %; R²=0.998).

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
FAST MasterMixes for Probe Assays											
FAST qPCR MasterMix Plus for Probe Assay, 7.5 ml	RT-QP2X-03+F	600	RT-QP2X-005+F	✓			✓				
FAST qPCR MasterMix Plus for Probe Assay dTTP, 7.5 ml	RT-QP2X-03+WOUNF	600	RT-QP2X-005+WOUNF	✓			✓				
FAST qPCR MasterMix Plus for Probe Assay Low ROX, 7.5 ml	RT-QP2X-03+WOLURF	600	RT-QP2X-005+WOLURF		✓			✓		✓	
FAST qPCR MasterMix Plus for Probe Assay No ROX, 7.5 ml	RT-QP2X-03+NRFB	600	RT-QP2X-005+NRFB			✓			✓		✓
FAST 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+WOLURFB	600	RT-QP2X-005+WOLURFB		✓			✓		✓	
FAST BLUE qPCR MasterMix Plus for Probe Assay No ROX, 7.5 ml	RT-QP2X-03+NRFB	600	RT-QP2X-005+NRFB			✓			✓		✓
FAST MasterMixes for SYBR® Assays											
MESA FAST qPCR MasterMix Plus for SYBR® Assay, 7.5 ml	RT-SY2X-03+WOUF	600	RT-SY2X-005+WOUF	✓			✓				
MESA FAST qPCR MasterMix Plus for SYBR® Assay Low ROX, 7.5 ml	RT-SY2X-03+WOLURF	600	RT-SY2X-005+WOLURF		✓			✓		✓	
MESA FAST qPCR MasterMix Plus for SYBR® Assay No ROX, 7.5 ml	RT-SY2X-03+NRWOUF	600	RT-SY2X-005+NRWOUF			✓			✓ (*)		✓

(*) Fluorescein must be added to the MasterMix before use • Ref. RT-FLUO-ADD