

UNIVERSAL EXOGENOUS qPCR POSITIVE CONTROL



TURN UNCERTAINTY INTO ACCURACY
USE A VALIDATED qPCR POSITIVE CONTROL!

TURN UNCERTAINTY INTO ACCURACY

Real-Time PCR assays are prone to inhibition by various substances found in many samples (clinical, soil, plant...). Carryover of reagents used for the isolation of nucleic acids can also inhibit amplification reactions. Other causes of false-negative results include target nucleic acid degradation, sample processing errors and thermocycler malfunction. Eurogentec's Universal Exogenous qPCR positive Control provides an accurate way to assess the integrity of all the steps in a nucleic acid amplification assay.

- > Distinguish true target negatives from false negatives
- > Monitor loss or degradation of your target during extraction
- > Assess the quality of your nucleic acids extracts

Specific

- > Does not interfere with target amplification
- > Avoid amplification of endogenous genes

Sensitive

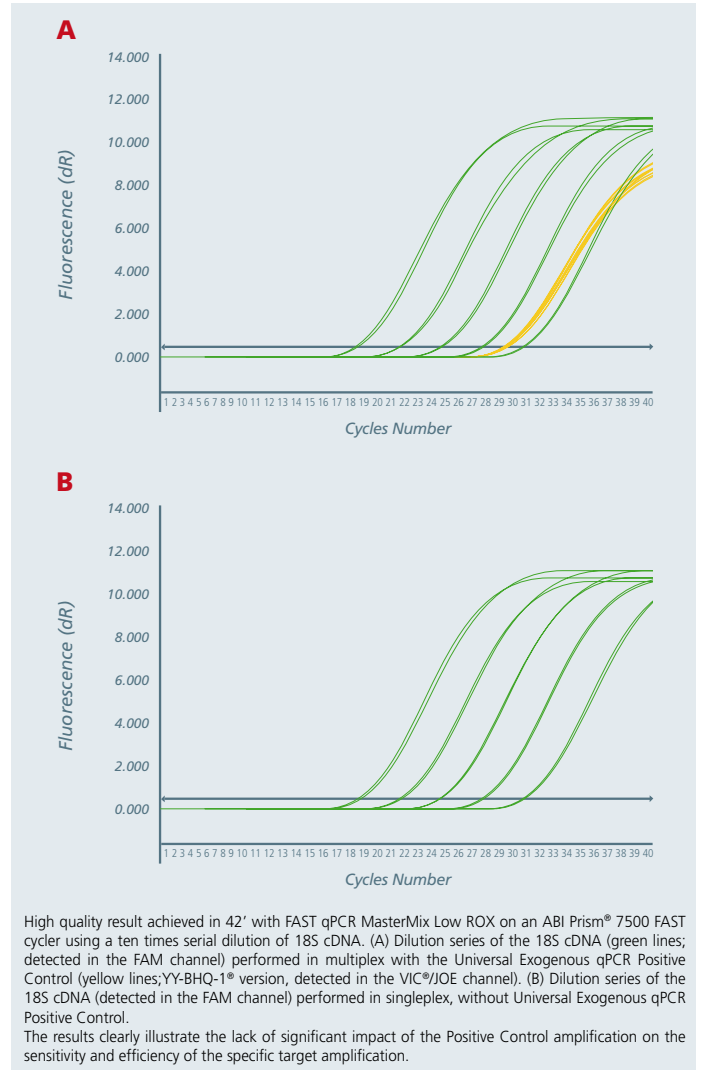
- > Compatible with low copy templates
- > Available with dark quencher (BHQ®) for maximal signal-to-noise ratio

Designed for multiplex assays

- > Detected on the Yakima Yellow®/VIC®/JOE channel available on all thermocyclers
- > Validated in duplex assays with the target in the FAM channel

Amplification		Results Interpretation
Target	IPC	
+	+	Target Positive
-	+	Target Negative
-	-	False Negative ! Reject Result !

Result interpretation of a diagnostic test performed in the presence of the Universal Exogenous qPCR Positive Control. A negative target and IPC amplification suggests the presence of inhibitors, a wrong PCR set-up, a defective mix or an incorrect thermal cycling protocol.



Product	Reference	#RXN (50 µl)	Price *		
			EUR	CHF	GBP
Universal Exogenous qPCR Positive Control, YY-TAMRA	RT-IPCY-T02	200	220	345	175
	RT-IPCY-T10	1000	990	1555	785
Universal Exogenous qPCR Positive Control, YY-BHQ-1®	RT-IPCY-B02	200	220	345	175
	RT-IPCY-B10	1000	990	1555	785

* Mentioned prices are VAT excluded

BHQ-1® is a registered trademark of Biosearch Technologies, Inc., VIC® is a registered trademark of Applied Biosystems, Yakima Yellow® is a registered trademark of Epoch Biosciences, Inc., ABI Prism® is a registered trademark of The Perkin-Elmer Corp.

How to order

- By email at order@eurogentec.com, for UK: order.uk@eurogentec.com
- By fax to: +32 (0) 4 264 07 88
- To learn more about these products and exceptional savings, visit www.eurogentec.com/promotions.htm

