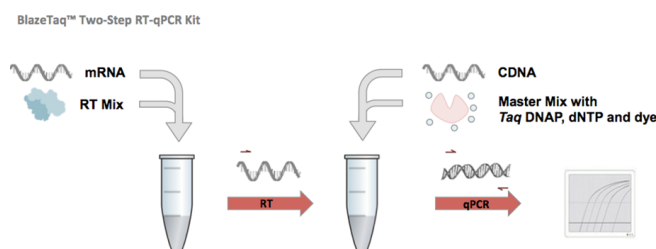


BlazeTaq™ Two-Step RT-qPCR System

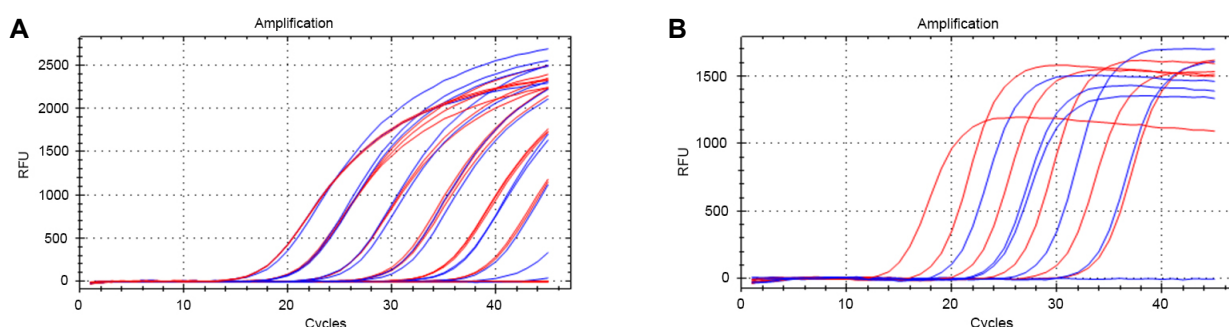
GeneCopoeia's BlazeTaq™ Two-Step SYBR® Green RT-qPCR System includes a SureScript™ First-Strand cDNA Synthesis Kit and a BlazeTaq™ qPCR Master Mix. The 5X qPCR master mix contains an antibody-modified hot start Taq DNA polymerase and optimized buffer system that allows wide dynamic range, high specificity and fast cycling time. Simply add DNA template, primers, and water to start your reaction. The First-Strand cDNA Synthesis Kit uses an optimized reaction buffer, M-MLV reverse transcriptase and murine RNase inhibitor.



Advantages

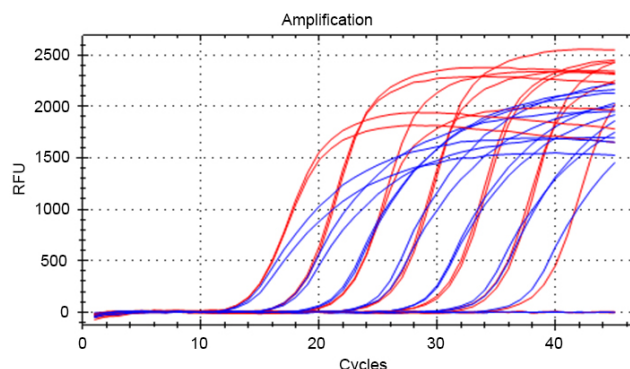
- ⦿ High sensitivity. Detects as little as 10 pg RNA template
- ⦿ High specificity with minimal level of primer-dimer and non-specific product formation
- ⦿ High amplification efficiency over wide GC-content range

BlazeTaq™ Two-Step RT-qPCR Performance



Comparison of the real-time amplification plots of the BlazeTaq™ Two-Step RT-qPCR System (red) with competitor A's (A) and competitor B's (B) products (blue). A 10-fold dilution series of B2M gene from total RNA of HeLa cell extract with amounts ranging from 100 ng to 10 pg.

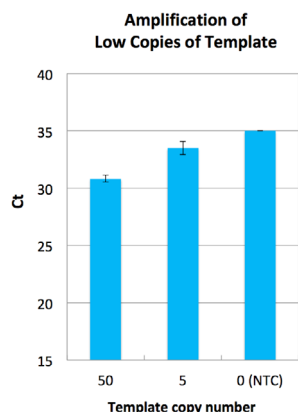
BlazeTaq™ SYBR Green qPCR Mix Performance



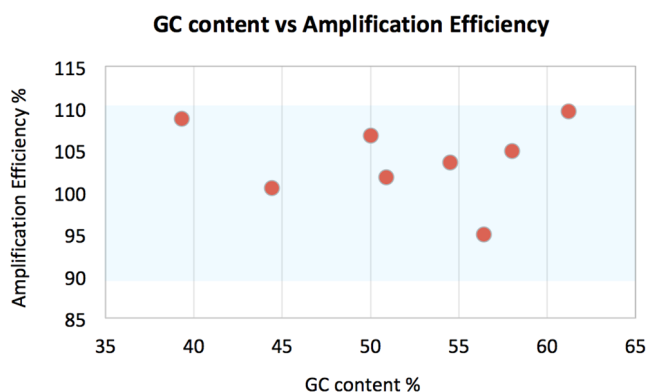
Comparison of the real-time amplification plots of the BlazeTaq™ SYBR Green qPCR Master Mix (red) with competitor A's product (blue). A 10-fold dilution series of input template concentrations from 100 ng to 10 pg. cDNA was generated from HeLa total RNA using GeneCopoeia's SureScript™ First-Strand cDNA Synthesis Kit.

BlazeTaq™ SYBR Green qPCR Mix Performance

A



B



Performance Analysis of the BlazeTaq™ qPCR Master Mix. **A.** Mean Ct values from real-time PCR of 5 and 50 copies of synthetic DNA template amplified with BlazeTaq™ SYBR Green qPCR Mix. The reactions were repeated for 24 times for each template. **B.** Amplification efficiency of cDNA templates with various GC contents ranging from 90%~110%.

Order information for BlazeTaq™ Two-Step RT-qPCR System and related products

Catalog #	Product	Description
SureScript™ First-Strand cDNA Synthesis Kit		
QP056	SureScript™ First-Strand cDNA Synthesis Kit (20 rxn)	M-MLV RT (RNase H-), Reaction Buffer, RNase Inhibitor, dNTPs, Oligo-(dT), Random Primer, ddH2O and spike-in control
QP057	SureScript™ First-Strand cDNA Synthesis Kit (60 rxn)	M-MLV RT (RNase H-), Reaction Buffer, RNase Inhibitor, dNTPs, Oligo-(dT), Random Primer, ddH2O and spike-in control

BlazeTaq™ SYBR Green qPCR Master Mix

QP031	BlazeTaq™ SYBR Green qPCR Mix 2.0 (200 rxn, with ROX)	High-fidelity, hot-start DNA polymerase, ROX reference dye, optimized reaction buffer and dNTPs
QP032	BlazeTaq™ SYBR Green qPCR Mix 2.0 (600 rxn, with ROX)	High-fidelity, hot-start DNA polymerase, ROX reference dye, optimized reaction buffer and dNTPs
QP034	BlazeTaq™ SYBR Green qPCR Mix 2.0 (1000 rxn, with ROX)	High-fidelity, hot-start DNA polymerase, ROX reference dye, optimized reaction buffer and dNTPs
QP041	BlazeTaq™ SYBR Green qPCR Mix 2.0 (200 rxn, without ROX)	High-fidelity, hot-start DNA polymerase, optimized reaction buffer and dNTPs
QP042	BlazeTaq™ SYBR Green qPCR Mix 2.0 (600 rxn, without ROX)	High-fidelity, hot-start DNA polymerase, optimized reaction buffer and dNTPs
QP044	BlazeTaq™ SYBR Green qPCR Mix 2.0 (1000 rxn, without ROX)	High-fidelity, hot-start DNA polymerase, optimized reaction buffer and dNTPs

Validated qPCR Primers

Variable	All-in-One™ qPCR Primer (200 reactions)	Validated gene-specific for human, mouse and rat to ensure specificity and sensitivity
-----------------	---	--

Contact us:

GeneCopoeia, Inc.

9620 Medical Center Dr. Suite 101,
Rockville, MD 20850 USA

Email inquiry@genecopoeia.com
Tel +1 (301) 762-0888
Toll free +1 (866) 360-9531
Fax +1 (301) 762-3888
Website www.genecopoeia.com