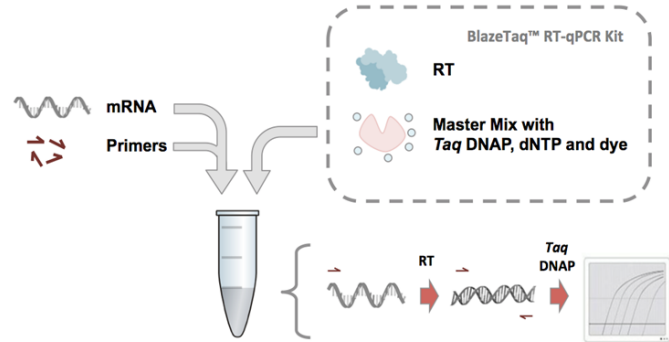


BlazeTaq[™] One-Step SYBR[®] Green RT-qPCR Kit

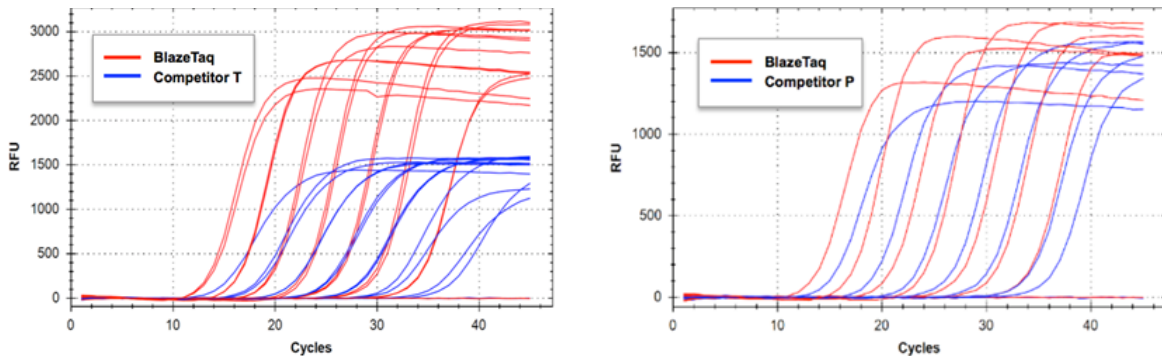
BlazeTaq[™] One-Step SYBR[®] Green RT-qPCR kit is based on quantitative reverse transcription PCR (RT-qPCR) that uses RNA as starting material. The kit is supplied with reverse transcriptase and a 5X master mix with a hot-start antibody-modified Taq DNA polymerase, dNTP and all required buffer components. BlazeTaq[™] One-Step SYBR[®] Green RT-qPCR kit is compatible with diverse real-time PCR instruments and offers best-in-class performance for RNA quantification.



Advantages

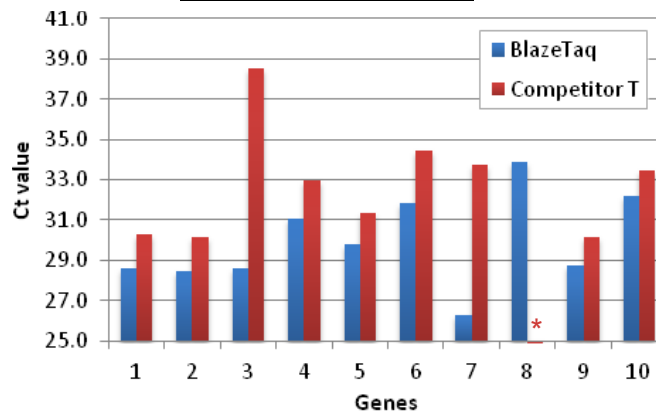
- ⊙ High sensitivity-accurate detection of target RNA from as little as 0.1 pg RNA extract
- ⊙ High specificity, with minimal level of primer-dimer and non-specific product formation
- ⊙ Best-in-class performance with earlier Ct and stronger signals compared to competitor's products

High sensitivity



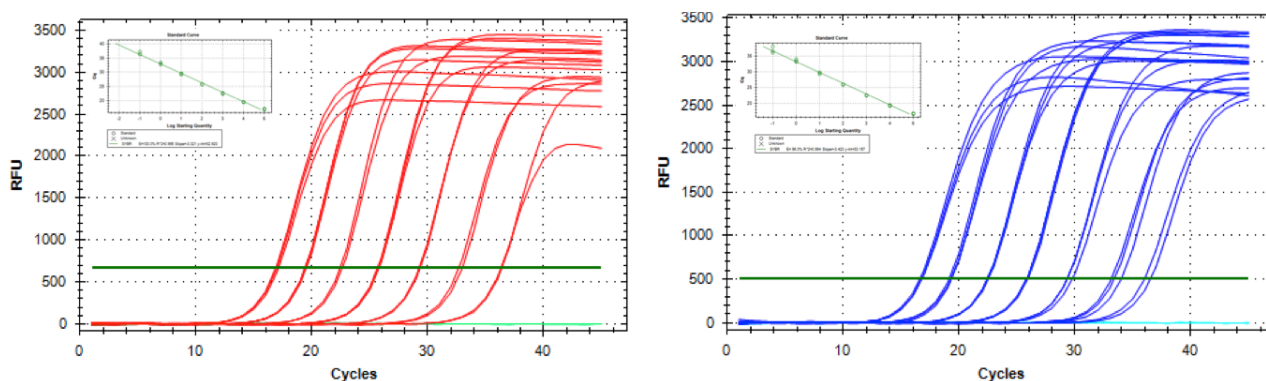
Amplification curves of GAPDH in RNA extracts from HeLa cells ranging from 100 ng to 0.1 pg in 10-fold dilution series with BlazeTaq[™] (red) and competitor's kit (blue) showed increased fluorescent signals and earlier Ct values.

Earlier Ct values



Earlier Ct values achieved from BlazeTaq[™] One-Step RT-qPCR kit with genes tested and compared with competitor's kit.
* Not detected.

Great Amplification Efficiency



Analysis of the level of GAPDH (left) and Actin (right) in RNA extract ranging from 100 ng to 0.1 pg from HeLa cells with amplification efficiency greater than 95%.

Order Information

Catalog #	Product	Description
QP071	BlazeTaq™ One-Step SYBR Green RT-qPCR Kit (200 rxn, with ROX)	High-fidelity, hot-start DNA polymerase, ROX reference dye, optimized reaction buffer and dNTPs
QP072	BlazeTaq™ One-Step SYBR Green RT-qPCR Kit (600 rxn, with ROX)	High-fidelity, hot-start DNA polymerase, ROX reference dye, optimized reaction buffer and dNTPs
QP074	BlazeTaq™ One-Step SYBR Green RT-qPCR Kit (1000 rxn, with ROX)	High-fidelity, hot-start DNA polymerase, ROX reference dye, optimized reaction buffer and dNTPs
QP081	BlazeTaq™ One-Step SYBR Green RT-qPCR Kit (200 rxn, w/o ROX)	High-fidelity, hot-start DNA polymerase, optimized reaction buffer and dNTPs
QP082	BlazeTaq™ One-Step SYBR Green RT-qPCR Kit (600 rxn, w/o ROX)	High-fidelity, hot-start DNA polymerase, optimized reaction buffer and dNTPs
QP084	BlazeTaq™ One-Step SYBR Green RT-qPCR Kit (1000 rxn, w/o ROX)	High-fidelity, hot-start DNA polymerase, optimized reaction buffer and dNTPs

Contact us:

GeneCopoeia, Inc.

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

Email

Tel

Toll free

Fax

Website

inquiry@genecopoeia.com

+1 (301) 762-0888

+1 (866) 360-9531

+1 (301) 762-3888

www.genecopoeia.com