

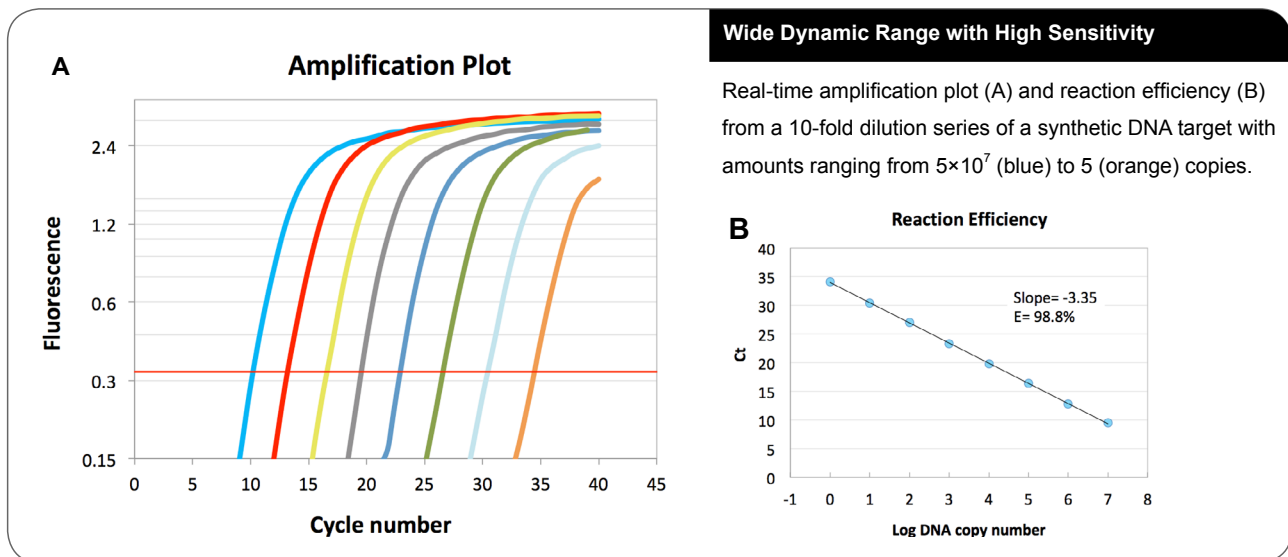


BlazeTaq™ SYBR® Green qPCR Mix

GeneCopoeia's BlazeTaq™ SYBR® Green qPCR Mix is a ready-to-use mix for real-time quantitative PCR (qPCR) on all qPCR instruments. The 2× master mix contains an antibody-modified hot start Taq DNA polymerase and optimized buffer system that provides a wide dynamic range, high specificity and fast cycling time. Simply add DNA template, primers, and sterile water to start your reaction.

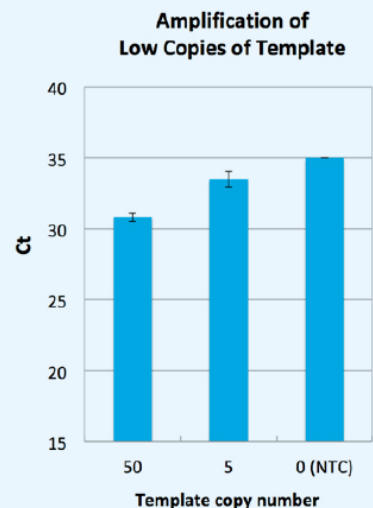
Advantages

- ⊙ High sensitivity. Detects as little as 5 copies of DNA template
- ⊙ High specificity with minimal levels of primer-dimer and non-specific product formation
- ⊙ High amplification efficiency over wide GC-content range

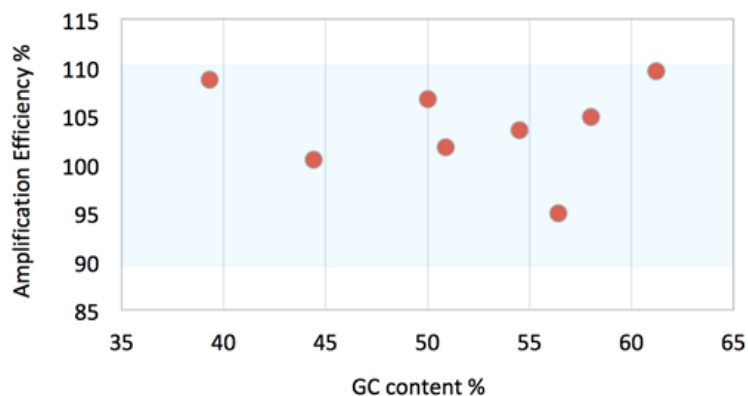


Consistent detection of low copy number DNA templates

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 24 times for each template.



GC content vs Amplification Efficiency



Accurate analysis of templates with wide range of GC content

Amplification efficiency of cDNA templates with various GC content ranging from 90%~110%.

Ordering information for BlazeTaq™ SYBR® Green qPCR Mixes and related products

Catalog #	Product	Description	Price
QP051	BlazeTaq™ SYBR® Green qPCR Mix (200 rxn, with ROX)	High-fidelity, hot-start DNA polymerase, ROX reference dye, optimized reaction buffer and dNTPs	\$199
QP052	BlazeTaq™ SYBR® Green qPCR Mix (600 rxn, with ROX)	High-fidelity, hot-start DNA polymerase, ROX reference dye, optimized reaction buffer and dNTPs	\$465
QP054	BlazeTaq™ SYBR® Green qPCR Mix (1000 rxn, with ROX)	High-fidelity, hot-start DNA polymerase, ROX reference dye, optimized reaction buffer and dNTPs	\$760
QP061	BlazeTaq™ SYBR® Green qPCR Mix (200 rxn, without ROX)	High-fidelity, hot-start DNA polymerase, optimized reaction buffer and dNTPs	\$199
QP062	BlazeTaq™ SYBR® Green qPCR Mix (600 rxn, without ROX)	High-fidelity, hot-start DNA polymerase, optimized reaction buffer and dNTPs	\$465
QP064	BlazeTaq™ SYBR® Green qPCR Mix (1000 rxn, without ROX)	High-fidelity, hot-start DNA polymerase, optimized reaction buffer and dNTPs	\$760
QP008	All-in-One™ First-Strand cDNA Synthesis Kit for gene qPCR array (20 RT reactions)	M-MLV RT (RNase H-), Reaction Buffer, RNase Inhibitor, dNTPs, Oligo(dT)18, Random Primer, dd H2O and spike-in control (for use with gene qPCR arrays)	\$129
QP009	All-in-One™ First-Strand cDNA Synthesis Kit for gene qPCR array (60 synthesis reactions)	M-MLV RT (RNase H-), Reaction Buffer, RNase Inhibitor, dNTPs, Oligo(dT)18, Random Primer, dd H2O and spike-in control (for use with gene qPCR arrays)	\$285
Variable	All-in-One™ qPCR Primer (200 reactions)	Validated gene-specific for human, mouse and rat to ensure specificity and sensitivity	

Contact us:

GeneCopiaea, Inc.

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

Email
Tel
Toll free
Fax
Website

inquiry@genecopiaea.com
+1 (301) 762-0888
+1 (866) 360-9531
+1 (301) 762-3888
www.genecopiaea.com