

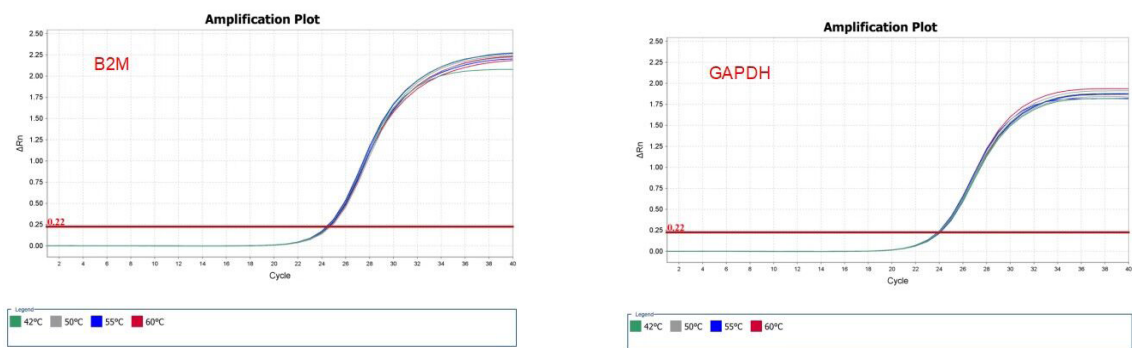
## POPULAR ENZYMES FOR SINGLE CELL RNA-SEQ LIBRARIES

If you are interested in these products, please feel free to contact us.

Tel: 301-762-0888 or 1-888-860-4093 Email: [Inquiry@genecopoeia.com](mailto:Inquiry@genecopoeia.com)

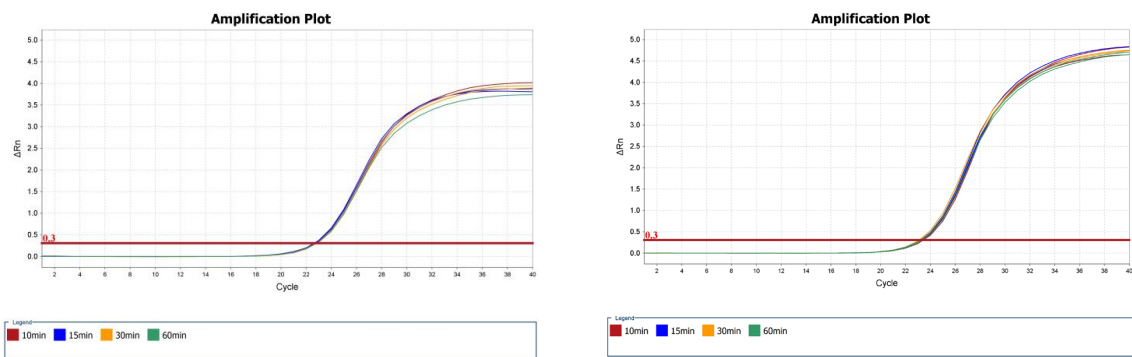


## Broad thermostability



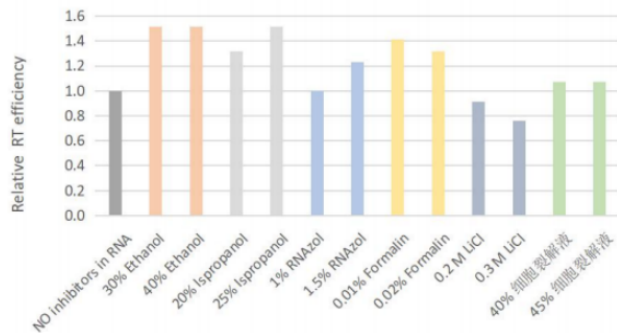
**Figure 4.** Broad thermal stability of AccelerRT® 5G Template Switching RT Enzyme Mix. 1ng Hela total RNA was reverse transcribed using AccelerRT® 5G Template Switching Reverse Transcriptase at different temperatures from 42 °C to 60°C. The GAPDH and B2M cDNAs were amplified with BlazeTaq™ SYBR® Green qPCR mix 2.0(Cat.# QP031).

## High efficiency of reverse transcription



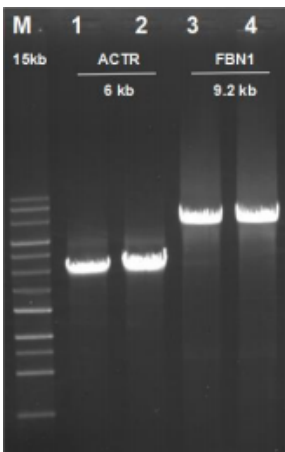
**Figure 5.** RT efficiency of AccelerRT® 5G Template Switching RT Enzyme Mix was demonstrated using 1ng Hela total RNA reverse transcribed for 10min, 15min, 30min and 60min. The GAPDH and B2M cDNAs were amplified using BlazeTaq™ SYBR® Green qPCR mix 2.0(Cat.# QP031).

## Consistent performance in the presence of a variety of inhibitors



**Figure 6.** Various inhibitors were added to total HeLa RNA, then was used in a reverse transcription reaction with AccelerRT® 5G Template Switching RT Enzyme Mix, The synthesized cDNA was used as a template in subsequent qPCR using BlazeTaq™ SYBR® Green qPCR mix 2.0(Cat.# QP031).

## Effective amplification of cDNA Synthesis



**Figure 7.** Total RNA from Hela cells was used in a reverse transcription reaction with AccelerRT® 5G Template Switching RT Enzyme Mix. The synthesized cDNA was used as a template in subsequent PCR using the UltraHiPF® DNA Polymerase Kit (Cat.# PC018).

## Discovery and detection of fusion genes



**Figure 8.** Total RNA from H2228 cells was used in a reverse transcription reaction with AccelerRT® 5G Template Switching RT Enzyme Mix.

The synthesized cDNA was used as a template in subsequent PCR using the 5' -end portion of the TSO primer and the gene-specific primer ALK gene(A). The amplification products in Figure A was verified by amplification with the EML4 forward primer and the ALK reverse primer(B).

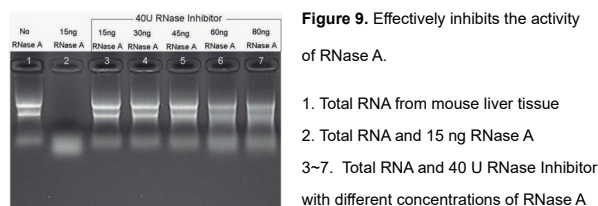
## 2. RNaseLock™ RNase Inhibitor

### Applications:

Effectively inhibits the activity of RNase A, RNase B and RNase C in eukaryotes.

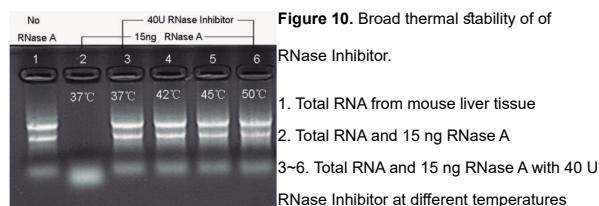
In experiments with potential RNase contamination:

- cDNA synthesis
- RT-PCR
- In vitro transcription
- Isolation and purification of mRNA



**Figure 9.** Effectively inhibits the activity of RNase A.

1. Total RNA from mouse liver tissue
2. Total RNA and 15 ng RNase A
- 3~7. Total RNA and 40 U RNase Inhibitor with different concentrations of RNase A



**Figure 10.** Broad thermal stability of of RNase Inhibitor.

1. Total RNA from mouse liver tissue
2. Total RNA and 15 ng RNase A
- 3~6. Total RNA and 15 ng RNase A with 40 U RNase Inhibitor at different temperatures

## 3. T4 DNA ligase

### Applications:

- Cloning of fragments cutting by restriction endonuclease
- Connect linkers or adapters to the blunt ends of DNA fragments
- Site-directed mutagenesis

## Contact us:

### GeneCopoeia, Inc.

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

Email [inquiry@genecopoeia.com](mailto:inquiry@genecopoeia.com)  
Tel +1 (301) 762-0888  
Fax +1 (301) 762-3888  
Website [www.genecopoeia.com](http://www.genecopoeia.com)