

## RNA Expression Analysis Webinar References

### *Promega PubHub articles*

Hendricksen, A. et al. [\*\*RNase Contamination Happens; Recombinant RNasin® Inhibitor Can Safeguard Your Samples.\*\*](#)

Promega Corporation. [\*\*Getting the Best Results from Your Reverse Transcription Quantitative PCR Assay.\*\*](#)

[Ammerschläger, M. et al. \*\*Lining Up the Scripts: Reverse Transcriptase Comparison Study.\*\*](#)

### *PubMed*

Tuma, R.S. et al. [\*\*Characterization of SYBR Gold nucleic acid gel stain: a dye optimized for use with 300-nm ultraviolet transilluminators.\*\*](#) Anal Biochem. 1999, 268(2), 278-88.

Kirsanov, K.I. et al. [\*\*SYBR Gold and SYBR Green II are not mutagenic in the Ames test.\*\*](#) Mutation Research/Genetic Toxicology and Environmental Mutagenesis. 2010, 1–2 (17), 1–4.

### *Previous webinars*

[\*\*Maximize Your Reverse Transcription-qPCR \(RT-qPCR\) Assays\*\*](#) Speaker: Dr. Carl Strayer

### *Promega Products*

[ReliaPrep™ RNA Cell Miniprep System](#)

[Maxwell® 16 LEV simplyRNA Kits](#)

[RNasin® Ribonuclease Inhibitors](#)

[GloMax®-Multi+ Microplate Multimode Reader](#)

[QuantiFluor™ Single-Tube Fluorometers](#)

[QuantiFluor™ RNA System](#)

[Ethidium Bromide Solution, Molecular Grade](#)

[GoTaq® Real-Time PCR Systems](#)

[Plexor® qPCR and qRT-PCR Systems](#)

[GoScript™ Reverse Transcription System](#)

[GoTaq® DNA Polymerase products](#)

[GoTaq® Real-Time PCR Systems](#)