

RNA extraction & QC:

Promega publications

- “Working with RNA” - <http://promega.wordpress.com/2012/01/24/working-with-rna/>
- “RNase Contamination Happens; Recombinant RNasin® Inhibitor Can Safeguard Your Samples. <http://www.promega.com/resources/articles/pubhub/rnase-contamination-happens-recombinant-rnasin-inhibitor-can-safeguard-your-samples/>
- “RNA Purification Kit Comparison: Yield, Quality and Real-Time RT-PCR Performance” - <http://www.promega.com/resources/articles/pubhub/promega-notes-2008/rna-purification-kit-comparison-yield-quality-and-real-time-rt-pcr-performance/>

RT-qPCR Resources (including those referenced in the webinar):

General

Gene Quantification - <http://www.gene-quantification.info/>

A-Z of quantitative PCR (Editor: S.A. Bustin), International University Line (IUL), La Jolla, CA, USA

Design, validation, controls; normalization -

- MIQE Guidelines - Minimum information for publication of quantitative real-time PCR experiments - <http://www.rdml.org/miqe.php> (Bustin et al., Clinical Chemistry, 2009)
- [Sean Taylor, “A MIQE Case Study — Effect of RNA Sample Quality and Reference Gene Stability on Gene Expression Data” 2011, Bio-Rad tech note 6245](#)
- Garbarino-Pico, E., Rollag, M.D., Strayer, C.A., Niu, S., Besharse, J.C., and Green, C.B. (2007) Immediate early response of the circadian polyA ribonuclease nocturnin to two extracellular stimuli. RNA 13: 745-755. [PMID:17400819](https://pubmed.ncbi.nlm.nih.gov/17400819/)

Pre-designed qPCR Primers:

- *Primer Bank* - <http://pqa.mgh.harvard.edu/primerbank/>
- *RTPrimerDB* - <http://www.rtpimerdb.org/>

Primer design software:

- *Primer3* - http://www-genome.wi.mit.edu/cgi-bin/primer/primer3_www.cgi
- *Primer-BLAST* - <http://www.ncbi.nlm.nih.gov/tools/primer-blast/>

Real-Time qPCR Analysis Software:

geNORM - <http://medgen.ugent.be/~jvdesomp/genorm/>
Reference gene normalization software

QPCR - <https://rtqcr.genome.tugraz.at/rtqcr/>
Web-based Real-Time PCR data management and analysis

Real-Time PCR Terms:

ΔR – baseline-corrected fluorescence - R *minus* average R over baseline region

Passive reference – free dye (unassociated with amplification product) used for normalization
(controls for experimental/optical variation)

R_n – normalized fluorescence (R divided by passive reference)

ΔR_n – baseline-corrected, normalized fluorescence.

Experimental Target – target gene/transcript/channel/dye of interest

Control Target – normalizer, endogenous control, reference gene/transcript/channel/dye,
housekeeping gene

Experimental Sample – unknown, treated, e.g., tumor, mutant, +2 hours, etc.

Control Sample – calibrator, reference sample, e.g., untreated control, normal, time zero, wild
type, etc.