

# A Practical Approach to RT-qPCR — Publishing Data That Conform to the MIQE Guidelines

Sean Taylor, Michael Wakem, Greg Dijkman, Marwan Alsarraj, and Marie Nguyen  
Bio-Rad Laboratories, Inc., Hercules, CA 94547

Tech  
Note

Amplification

Bulletin 5859

## Introduction

Real-time quantitative PCR (qPCR) has become a definitive technique for quantitating differences in gene expression levels between samples. Over the past 10 years, the popularity of this method has grown exponentially, with the publication of well over 25,000 papers from diverse fields of science, including agricultural, environmental, industrial, and medical research, making reference to reverse transcription qPCR (RT-qPCR) data.

Apart from the broad applicability of the technique, one of the central factors that has stimulated this impressive growth is the increased demand from journal review panels for the use of RT-qPCR to support phenotypic observations with quantitative, molecular data. Furthermore, gene expression analysis is now being used to support protein expression data from proteomics-based assays. The biotechnology industry has responded to the rapid adoption of this technique with the concomitant development of reagents and instruments to perform RT-qPCR experiments. However, since no strict guidelines have been established, researchers have generally designed their experiments based on information gathered from disparate sources, which has resulted in data of variable quality.

A number of technical deficiencies can affect RT-qPCR assay performance, including improper experimental design, inadequate controls and replicates, lack of well-defined experimental conditions and sample handling techniques, poor quality of the RNA sample, suboptimal choice of primers for reverse transcription and qPCR reactions, lack of validation of reference genes, and inappropriate methods of data analysis. In an effort to assist the scientific community in producing consistent, high-quality data from qPCR experiments, the minimum information for publication of quantitative real-time PCR experiments (MIQE) guidelines have been recently published (Bustin et al. 2009). This has been followed by the development of an XML-based real-time PCR data markup language (RDML) for the consistent reporting of real-time PCR data by the RDML consortium (Lefever et al. 2009). This consortium is active in the development of appropriate and standardized terminology, guidelines on minimum information for biological and biomedical investigations, and a flexible and universal data file structure with tools to create, process, and validate RDML files. All relevant information about the RDML project is available at [www.rdml.org](http://www.rdml.org).

The ultimate goal of RDML and MIQE is to establish a clear framework within which to conduct RT-qPCR experiments and to provide guidelines for reviewers and editors to use in the evaluation of the technical quality of submitted manuscripts against an established yardstick. As a consequence, investigations that use this widely applied methodology will produce data that are more consistent, more comparable, and ultimately more reliable.

Given the highly dynamic nature of mRNA transcription and the potential variables introduced in sample handling and in the downstream processing steps (Garson et al. 2009), a standardized approach to each step of the RT-qPCR workflow is critical for reliable and reproducible results. The MIQE provides this approach with a checklist that contains 85 parameters to ensure quality results that will meet the acceptance criteria of any journal (Bustin et al. 2009).

In this paper we demonstrate how to apply the MIQE guidelines ([www.rdml.org/miqe](http://www.rdml.org/miqe)) to establish a solid experimental approach.

**BIO-RAD**

## 1. Experimental Design

Proper experimental design is the key to any gene expression study. Since mRNA transcription can be sensitive to external stimuli that are unrelated to the processes studied, it is important to work under tightly controlled and well-defined conditions. Taking the time to define experimental procedures, control groups, type and number of replicates, experimental conditions, and sample handling methods within each group is essential to minimize variability (Table 1). Each of these parameters should be carefully recorded prior to conducting gene expression experiments to ensure good biological reproducibility for published data.

**Table 1. RT-qPCR experimental design and sample management.**

This table summarizes the workflow of a typical RT-qPCR experiment from experimental design to defining control groups, replicates, and experimental conditions to the detailed procedures for sample handling. This ultimately ensures that the key steps in RT-qPCR data production lead to high-quality, reproducible, publishable data.

Experiment Design	Control Groups	Replicates	Experiment Conditions	Sample Handling
Disease or treatment groups	Time course study (i.e., t = 0)	Biological (different sample per well)	Growth conditions (media and time or absorbance)	Precise time to harvest cells or tissue
Target genes implicated	Normal vs. disease (i.e., normal)	Technical (same sample per well)	Days of embryonic development	Sample extraction method
Potential reference genes	Untreated vs. drug treated (i.e., untreated)	—	Amount per mass of drug or compound	Preservation method and time
Number of data points to draw statistically significant conclusions	—	—	Sex, phenotype	Thaw and homogenization procedure
—	—	—	Incubation time	Total RNA extraction procedure

## 2. RNA Extraction

If samples must be collected over a period of time or in too large a number to process immediately, they should be stored in appropriate conditions (frozen at  $-80^{\circ}\text{C}$  and/or in RNA storage solution) until use. To minimize handling time during the RNA extraction procedure, it is recommended that samples be processed in relatively small batches of 10 to 20. The RNA extraction procedure should include a DNase I treatment step to remove any contaminating genomic DNA. Many commercially available kits such as the Aurum™ Total RNA Extraction Kits (Bio-Rad Laboratories, Inc.) can be used for virtually any sample type from plant to animal tissue and include an on-column DNase I treatment of the samples.

## 3. RNA Quality Control

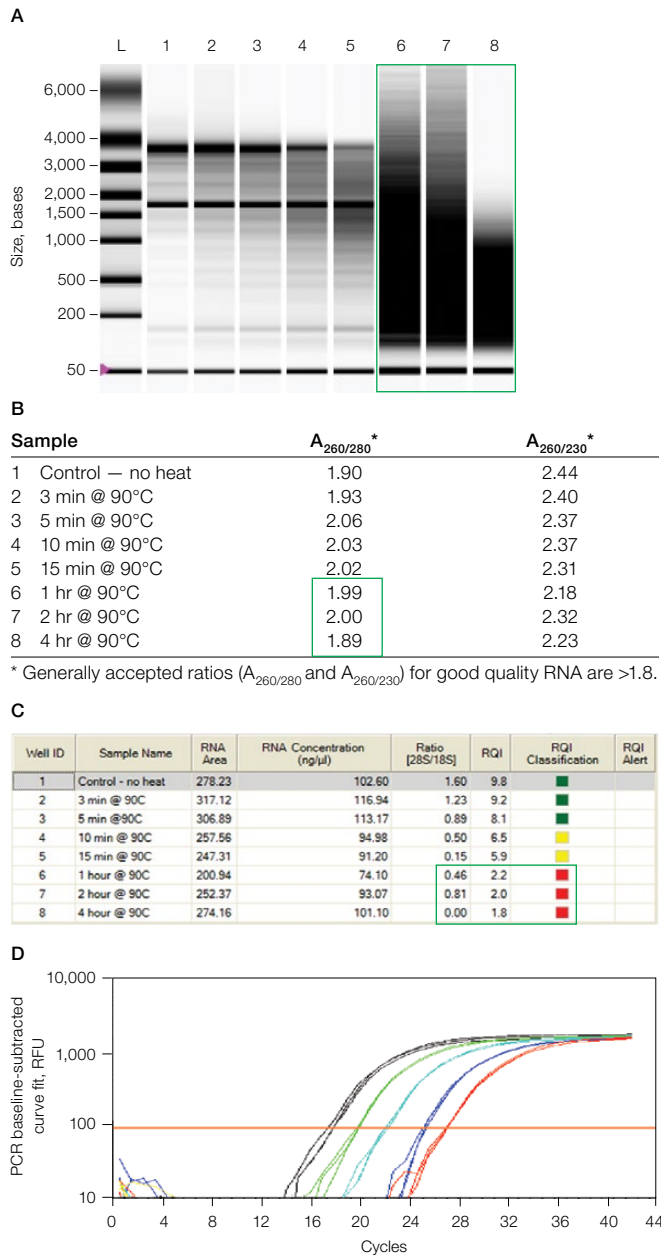
Ensuring that only RNA of high purity (no contaminants) and high integrity (not degraded) is used is one of the most critical points in the RT-qPCR experimental workflow. Impurities in

the RNA sample may lead to inhibition of the RT and PCR reaction, yielding biased data. Using partially degraded RNA can lead to varying and incorrect quantification results (Fleige and Pfaffl 2006, Gingrich et al. 2008). Since sample purity and integrity are not related, both should be assessed to ascertain that the RNA sample meets minimal acceptance criteria for the downstream workflow.

Purity of the sample with respect to protein contamination can be assessed spectrophotometrically by measuring the  $A_{260/280}$  ratio in a pH neutral buffer. An  $A_{260/280}$  ratio of 1.8 to 2.0 indicates good quality RNA that is devoid of protein and phenol contamination (Figure 1B). However, no RNA integrity information can be obtained from a spectrophotometric reading. RNA integrity can be assessed using several methods. The traditional method is visual inspection after electrophoresis on a formaldehyde agarose gel in the presence of a fluorescent dye such as ethidium bromide. Observation of two sharp bands for the large and the small subunit ribosomal RNAs (rRNA) with the intensity of the larger band being about twice that of the smaller band is indicative of intact RNA. While this method is inexpensive, the interpretation of the data is mostly subjective and requires approximately 200 ng of total RNA, which may be difficult if the sample is only available in limited quantities. This method can be refined by quantifying the intensity of the rRNA bands using an imager with densitometry scanning. While the values of the 28S/18S ratio can vary between different tissues or cell types, a ratio between 1 and 2 is indicative of an intact RNA sample.

A major improvement in RNA integrity analysis came with the introduction of microfluidics-based electrophoresis systems (Imbeaud et al. 2005) such as the Experion™ Automated Electrophoresis System (Bio-Rad) that combines integrity and concentration quantification of RNA in a single step from nanogram or picogram quantities (Figures 1A and 1C). In addition to generating a virtual gel, an electropherogram, and calculating the 28S/18S ratio, the Experion System's software automatically calculates and reports an RNA quality indicator (RQI) value (Figure 1C) that reflects the integrity of the RNA sample based on several criteria (Denisov et al. 2008) of the electropherogram. The RQI value ranging from 1 (degraded) to 10 (intact) provides an objective assessment of the integrity of the RNA sample and it can be used to screen out degraded samples.

By ensuring consistency in both purity and quality across all RNA samples, variability between biological replicates can be reduced (Figure 1D) (Imbeaud et al. 2005). When a batch of RNA samples has successfully met the standards of quality control, we recommend their immediate use in RT-qPCR experiments, or their conversion into much more stable cDNA by reverse transcription to preserve RNA integrity after the quality check.



**Fig. 1. Analysis of RNA purity and integrity.** This figure shows an example of RNA purity and integrity analysis using spectrophotometry and the Experion Automated Electrophoresis System. Mouse liver total RNA samples were subjected to degradation by heating at 90°C for different periods of time and analyzed for both purity and integrity. Green boxes highlight the fact that the total RNA sample appears by absorbance readings alone ( $\geq 1.8$ ) to be of sufficient quality, whereas RQI measurement on the Experion System clearly shows degraded sample with low RQI values. **A**, virtual gel image generated by the Experion Software showing various degradation levels of the RNA samples and a progressively decreasing intensity of the 18S rRNA band. **B**, the  $A_{260}/A_{280}$  ratio measured for all samples on the NanoDrop Spectrophotometer is between 1.8 and 2.0 indicating that the samples are devoid of protein contamination. **C**, summary of RNA integrity analysis using the Experion Software. The 28S/18S rRNA ratio and the RQI indicate a decreasing integrity of the heat-treated RNA samples. A color classification allows for easy identification of the samples that are not adequate for qPCR. **D**, example of qPCR analysis of *GAPDH* expression performed on degraded samples (Gingrich et al. 2008). The profiles show increasing quantification cycle (Cq) with the degraded samples. No degradation (—); 1 hr degradation (—); 3 hr degradation (—); 5 hr degradation (—); 8 hr degradation (—). RFU, relative fluorescence units.

#### 4. Reverse Transcription

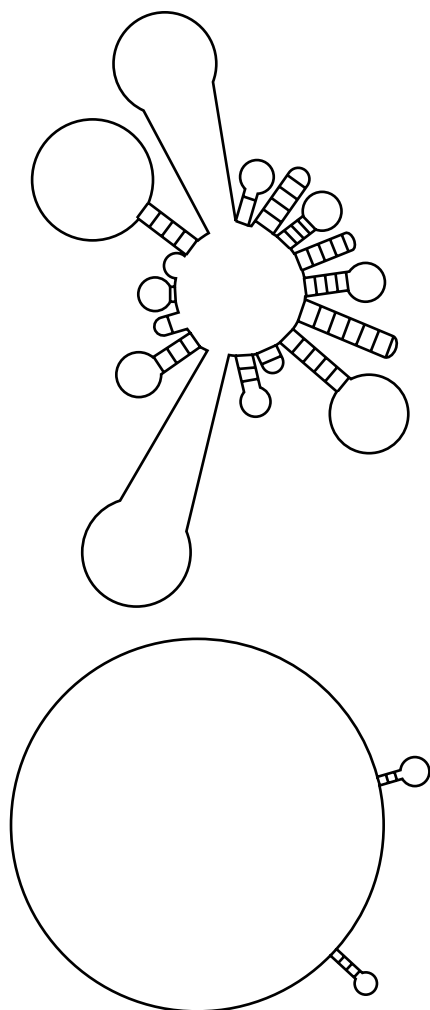
Given the prevalence of RNase in the environment, we recommend performing the reverse transcription of total RNA samples to cDNA immediately following the quality control assessment. This will avoid the risk of RNA sample degradation from multiple freeze/thaws before conversion to cDNA. For the RT step, the key is to ensure consistent and complete coverage of the transcribed genome in the extracted RNA sample. Some genes are very long, but the associated RT products for these sequences cannot be as large, especially if the RT primers anneal only at the ends of each mRNA. By annealing primers at the end of the mRNA plus at random points within each mRNA sequence, a good sampling of the population of each gene is obtained. This method, which is more representative than just annealing at the ends or at random sites within each transcript, provides better coverage of the transcribed genome.

A reverse transcription buffer should contain a mix of primers that are random in sequence, allowing for a better sampling of the mRNA; RNase H, the enzyme that specifically degrades RNA in DNA/RNA duplexes; a specific and sensitive reverse transcriptase with broad dynamic range for RNA amounts from 1  $\mu$ g to 1 pg and; a simple and fast protocol. The iScript™ cDNA Synthesis Kits (Bio-Rad) meet all of these criteria and are well suited for reverse transcription of any total RNA sample. We recommend that the same amount of total RNA be used and that reaction time be maintained for reverse transcription for all experimental samples to minimize variability between biological replicates. Reverse transcribed RNA can be stored frozen at  $-20^{\circ}\text{C}$  or  $-80^{\circ}\text{C}$  until use and diluted before use in the downstream qPCR reaction.

#### 5. Primer and Amplicon Design

Both primer design and careful choice of target sequence are essential to ensure specific and efficient amplification of the products. Target sequences should be unique, 75–150 bp long with a GC content between 50 and 60%, and should not contain secondary structures. It is recommended that primers should have a GC content of 50–60% and a melting temperature ( $T_m$ ) of 55–65°C. Long G or C stretches in the primer should be avoided, but it is recommended to have G or C at the end of the primers.

A number of programs are available to help design primer pairs and pick target sequences. We recommend designing oligonucleotides using Primer-Blast ([www.ncbi.nlm.nih.gov/tools/primer-blast/index.cgi?LINK\\_LOC=BlastHomeAd](http://www.ncbi.nlm.nih.gov/tools/primer-blast/index.cgi?LINK_LOC=BlastHomeAd)), a program developed by NCBI that uses the algorithm Primer3 (Rozen and Skaletsky 2000). Primer sequences are compared (blasted) to the user-selected databases to ensure they are unique and specific for the gene of interest. The program MFOLD (<http://unafold.rna.albany.edu/?q=mfold/dna-folding-form>) can then be used to analyze the amplicon for potential secondary structures that may prevent efficient amplification (Zuker 2003) (Figure 2). Ideally, two sets of oligonucleotides should then be ordered and tested for their performance in a qPCR reaction.



**Fig. 2. Effect of annealing temperature on amplicon secondary structure.** Analysis of the amplicon sequence using MFOLD predicts a high secondary structure at 60°C (top) and low secondary structure at 65°C (bottom).

## 6. qPCR Validation

A validated qPCR assay is one that has been assessed for the optimal range of primer annealing temperatures, reaction efficiency, and specificity using a standard set of samples (Bustin et al. 2009). This will ensure that the reaction conditions, buffers, and primers have been optimized and that the cDNA samples are not contaminated with inhibitors of Taq polymerase. Bio-Rad has created a practical web resource ([bio-rad.com/genomics/pcrsupport](http://bio-rad.com/genomics/pcrsupport)) for qPCR design and validation. The major points in assay validation are summarized below.

### Determination of Annealing Temperature, Melt Curve Analysis, Gel Analysis of Amplicon, and No Template Control

A critical step in a PCR reaction is the annealing of the primers to their target sequences. It has to be performed at the right temperature for the primers to anneal efficiently to their targets, while preventing nonspecific annealing and primer-dimer formation. The fastest way to determine optimal annealing

temperature is to use a thermal cycler equipped with a temperature gradient feature. All Bio-Rad thermal cyclers and real-time PCR instruments are available with a thermal gradient block option. A range of temperatures around the calculated  $T_m$  of the primers should be tested (Figure 3A).

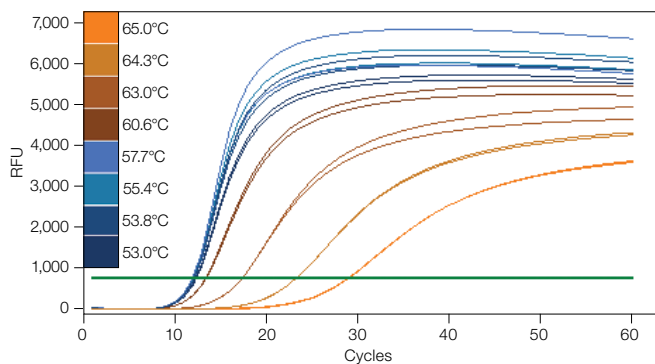
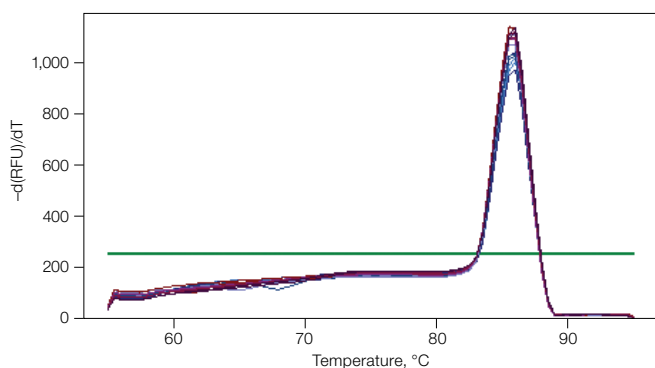
It is important to check the specificity of the reaction by analyzing the PCR product. A melt curve analysis, performed at the end of the PCR cycles, will confirm specificity of primer annealing. The melt curve should display a single sharp peak (Figure 3B). In addition, at least one sample per primer pair should be run on an appropriate gel (agarose gel or polyacrylamide gel to resolve smaller DNA fragments) to confirm that the amplicon is of the expected size (Figure 4). Alternatively, an automated electrophoresis system such as the Experion Automated Electrophoresis System (using the Experion DNA 1K Analysis Kit, Bio-Rad) can be used to rapidly and precisely confirm the size of the amplicon.

Since most qPCR experiments involve the study of multiple genes in a single protocol it is important to design primers with similar annealing temperatures and to run the reactions at a temperature that is adequate for all the primer pairs to be studied during the same run. A duplicate, no template control (NTC) reaction should be included in every run for each primer pair to test buffers and solutions for DNA contamination and to assess for primer-dimers.

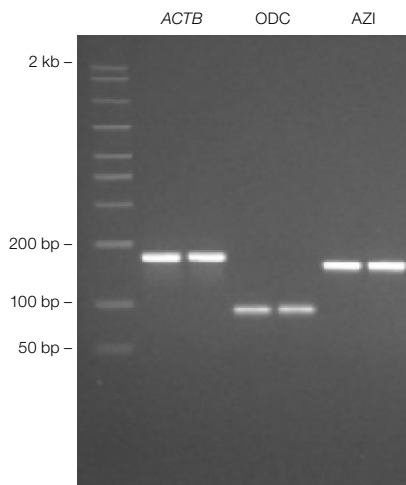
### Establishment of a Standard Curve to Evaluate PCR Efficiency

The efficiency of a PCR reaction is a measure of the rate at which the polymerase converts the reagents (dNTPs, oligonucleotides, and template cDNA) to amplicon. The maximum increase of amplicon per cycle is twofold, representing a reaction that is 100% efficient. It is important to measure reaction efficiency because it is indicative of problems with the qPCR reaction that can cause artifactual results. Low reaction efficiencies (<90%) may be caused by contaminating Taq inhibitors, high or suboptimal annealing temperature, old or inactive Taq, poorly designed primers, or amplicons with secondary structures. High reaction efficiency (>110%) is generally the result of primer-dimers or nonspecific amplicons. The most common causes of both high and low reaction efficiencies are poorly calibrated pipets or poor pipetting technique.

A standard curve is generally used to determine the reaction efficiency for any qPCR reaction. The template for this typically is a sample of cDNA or spiked plasmid cDNA in a sample extract. We recommend initially producing a tenfold dilution series over eight points starting from the most concentrated cDNA sample, to ensure the standard curve covers all potential template concentrations that may be encountered during the study (i.e., broad dynamic range). For each dilution, a standard qPCR protocol should be performed in triplicate for all the primer pairs to be used in the experiment and  $C_T$  values determined. CFX Manager™ Software versions 1.6 and

**A. Amplification****B. Melt Peak**

**Fig. 3. Validation of qPCR primers.** **A**, qPCR is performed at a range of annealing temperatures using a thermal gradient block. Amplification profiles indicate that the most efficient amplification occurs at the four lowest annealing temperatures between 53 and 57.7°C where the curves are at the lowest Cq. **B**, a single peak on the melt curve analysis indicates a single PCR product. RFU, relative fluorescence units.



**Fig. 4. Gel analysis of PCR products.** Gel electrophoresis analysis of the amplicon confirms the presence of a single fragment of the expected size. ODC, ornithine decarboxylase; AZI, antizyme inhibitor.

earlier provide  $C_T$  values for threshold cycle; however, MIQE guidelines recommend the more generic quantification cycle ( $C_q$ ) term. Since these terms are used interchangeably and refer to the same value,  $C_q$  will be used here. The standard curve is constructed by plotting the log of the starting quantity of the template against the  $C_q$  values obtained. The equation of the linear regression line, along with Pearson correlation coefficient ( $r$ ) or the coefficient of determination ( $r^2$ ), can then be used to evaluate whether the qPCR assay is optimized.

Ideally, the dilution series will produce amplification curves with tight technical replicates that are evenly spaced (Figure 5A). If perfect doubling occurs with each amplification cycle, the spacing of the fluorescence curves will be determined by the equation  $2^n = \text{dilution factor}$ , where  $n$  is the number of cycles between curves at the fluorescence threshold (in other words, the difference between the  $C_q$  values of the curves). For example, with a tenfold serial dilution of DNA,  $2^n = 10$ . Therefore,  $n = 3.32$ , and the  $C_q$  values should be separated by 3.32 cycles. Evenly spaced amplification curves (Figure 5A) will produce a linear standard curve (Figure 5B) with the goal of 90–110% reaction efficiency.

The  $r$  or  $r^2$  value of a standard curve represents how well the experimental data fit the regression line, that is, how linear the data are. Linearity, in turn, gives a measure of the variability across assay replicates and whether the amplification efficiency is the same for different starting template copy numbers. A significant difference in observed  $C_q$  values between replicates will lower the  $r$  or  $r^2$  value. An  $r$  with an absolute value  $>0.990$  or an  $r^2$  value  $>0.980$  is desirable for RT-qPCR reactions. Deletion of points at both ends of the standard curve may be required to obtain an acceptable slope (efficiency) and  $r^2$  value. This will ultimately define the dynamic range of cDNA concentration for each primer pair with respect to sample dilution.

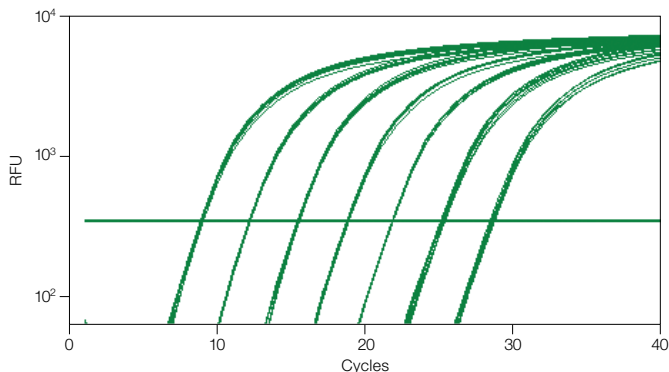
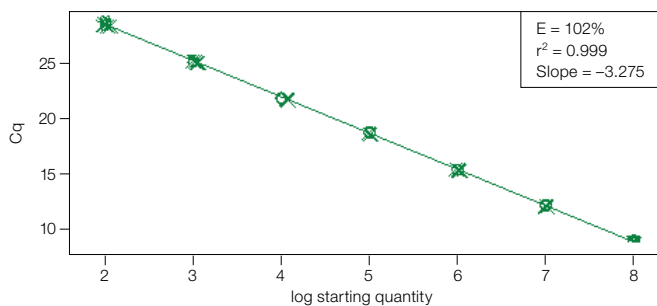
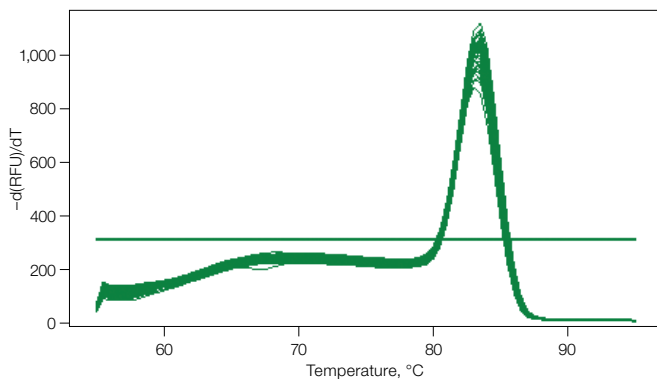
**qPCR Reagents, Instrument, and Analysis Program**

There are a wide variety of commercial qPCR reagent kits available. We recommend SsoAdvanced™ Universal Supermixes from Bio-Rad, which are ready-to-use reagents employing Bio-Rad's patented\* Sso7d fusion polymerase and advanced buffer formulation. These uniquely formulated supermixes can be used for a wide range of real-time PCR applications and with all real-time PCR systems.

A large number of qPCR detection instruments are commercially available and the following guidelines can be employed for selecting an appropriate instrument:

1. Sample volumes with a range of 10 to 50  $\mu$ l in standard 96-well format, 0.2 ml, low-profile thermal cycler plates, strips, or tubes from any manufacturer.
2. Gradient-enabled thermal block with fast ramping.
3. No requirement for proprietary reagents, consumables, or fluorophores.

\* U.S. patents 6,627,424; 7,541,170; and 7,560,260.

**A. Amplification****B. Standard Curve****C. Melt Peak**

**Fig. 5. Standard curve.** **A**, qPCR is performed on serial dilutions of the template and Cq is determined for each dilution. Three replicates are used for each concentration of template. Note that the Cq values for the replicates are very tight and nearly overlay perfectly indicating low technical variability. **B**, measured Cq values are plotted against the log of the copy number (or relative concentration) of the template to establish a standard curve that permits the calculation of the efficiency from the slope and  $r^2$  value. Ideally the three replicates per dilution should give very tightly grouped amplification curves with efficiency close to 100% (90–110%),  $r^2$  values above 0.98, and single, sharp melt curve peak per gene. **C**, melt curve analysis of the amplicons shows a single peak. Cq, quantification cycle; RFU, relative fluorescence units.

The CFX96 Touch™ and CFX384 Touch™ Real-Time PCR Detection Systems from Bio-Rad meet all of these criteria.

All qPCR instruments are packaged with data analysis software and we recommend that a good software package include the following features:

1. Flexibility to enter plate setup information such that well identifiers can be loaded and edited before, during, or after a run.
2. Ability to group wells from multiple experiments on one plate.
3. Built-in gene expression analysis that can normalize data to multiple reference genes (Vandesompele et al. 2002) and individual reaction efficiencies (Pfaffl 2001).
4. The ability to combine multiple plates of experimental data into a large gene study.

The CFX Manager Software from Bio-Rad meets all of these criteria and comes packaged with the CFX96 Touch and CFX384 Touch Real-Time PCR Detection Systems.

**7. Choice of Reference Genes**

In RT-qPCR experiments, reference genes are used as controls to normalize the data by correcting for differences in quantities of cDNA used as a template (Gutierrez et al. 2008, Huggett et al. 2005, Vandesompele et al. 2002). A perfect reference gene is therefore one that does not exhibit changes in expression between samples from various experimental conditions or time points. Several genes such as *GADPH*, *ACTB*, or rRNA are often used as reference genes. However, a number of studies have indicated that the expression of these genes may vary considerably between tissues or between treatments, which may make them unsuitable for use as reference genes. Reference genes must therefore be carefully selected based on experimental data and we recommend the following protocol:

1. Extract the total RNA from at least one or two samples from each experimental condition or time point and confirm their purity and quality (see sections 2 and 3).
2. Normalize the sample concentration and perform reverse transcription PCR (see section 4) from the same volume of each sample.
3. Perform the qPCR experiment using the same volume of each cDNA sample as a template.
4. Use the geNorm method (Hellemans et al. 2007) to calculate the target stability between the different conditions (available at <https://genorm.cmgg.be> or automatically calculated with CFX Manager Software).

A good reference gene should have an M value below 0.5 or 1 in homogeneous and heterogeneous sample sets, respectively (Vandesompele et al. 2002). geNorm also helps in the selection of the optimal number of reference genes. Typically, between three and five good reference genes are required to achieve the most accurate normalization (Vandesompele et al. 2002).

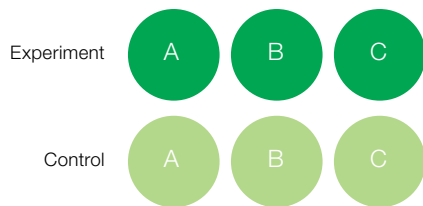
### 8. Experimental Reproducibility

There are two sources of variability in a gene expression experiment that may affect the results:

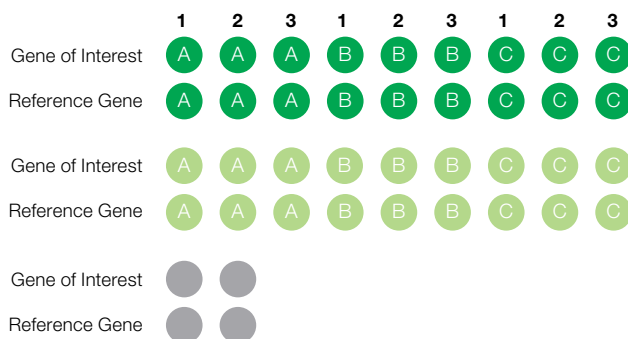
1. Biological variability which is due to inherent differences in gene expression levels between individual organisms, tissues, or cell culture samples.
2. Technical variability in the experimental process itself, which is typically associated with pipetting, poorly calibrated pipets, and sample quality and quantity.

To mitigate the effect of biological and technical variability, it is generally accepted that at least three biological and two technical replicates per biological replicate be performed for each experiment (Figure 6). If the experiment compares gene expression levels between control and treated samples, the three biological replicates should be samples that were treated in separate and independent experiments.

#### Biological Replicates



#### Technical Replicates



**Fig. 6. Experimental replicates.** All experiments should be designed with a combination of biological and technical replicates. This illustrates a simple experiment with triplicate biological samples from control (■) and treatment/experimental (■) conditions. For each biological sample, three technical replicates are recommended for the gene of interest as well as for the reference gene(s). This results in a total of at least 36 samples plus the duplicate NTC (■) for a total of >40 wells.

### Conclusions

RT-qPCR is the method of choice for gene expression analysis because of its high sensitivity from samples of very low RNA concentrations. However, in order to ensure accurate and reproducible quantitative data, strict standard operating procedures should be followed. All experimental details and controls should be accurately reported when publishing gene expression experiments. This will allow proper assessment of the data by the scientific community, and enable informed comparison of expression data between laboratories and between experiments.

In summary, the key steps for most RT-qPCR experiments include:

1. Experimental design with appropriate number of biological replicates and proper control samples.
2. Sample procurement which requires adherence to strict experimental protocols for acquisition, processing, and storage to ensure reproducibility and minimize standard deviations between replicates.
3. Quality control of RNA for purity and integrity.
4. Reverse transcription to convert the total RNA to cDNA.
5. qPCR experiments with proper primer design, choice of target sequence and reference genes, and technical replicates.

The MIQE guidelines were written to provide all the parameters that should be met to publish acceptable results from qPCR experiments. Here we have shown how the most important items in the MIQE checklist can be addressed in practice to ensure high-quality results.

## References

- Bustin SA et al. (2009). The MIQE guidelines: Minimum information for publication of quantitative real-time PCR experiments. *Clin Chem* 55, 611–622.
- Denisov V et al. (2008). Development and validation of RQI: An RNA quality indicator for the Experion™ automated electrophoresis system. *Bio-Rad Bulletin* 5761.
- Fleige S and Pfaffl MW (2006). RNA integrity and the effect on the real-time qRT-PCR performance. *Mol Aspects Med* 27, 126–139.
- Garson JA et al. (2009). Unreliable real-time PCR analysis of human endogenous retrovirus-W (HERV-W) RNA expression and DNA copy number in multiple sclerosis. *AIDS Res Hum Retroviruses* 25, 377–378; author reply 379–381.
- Gingrich J et al. (2008). Effect of RNA degradation on data quality in quantitative PCR and microarray experiments. *Bio-Rad Bulletin* 5452.
- Gutierrez L et al. (2008). The lack of a systematic validation of reference genes: A serious pitfall undervalued in reverse transcription-polymerase chain reaction (RT-PCR) analysis in plants. *Plant Biotechnol J* 6, 609–618.
- Hellemans J et al. (2007). qbase relative quantification framework and software for management and automated analysis of real-time quantitative PCR data. *Genome Biol* 8, R19.
- Huggett J et al. (2005). Real-time RT-PCR normalisation; strategies and considerations. *Genes Immun* 6, 279–284.
- Imbeaud S et al. (2005). Towards standardization of RNA quality assessment using user-independent classifiers of microcapillary electrophoresis traces. *Nucleic Acids Res* 33, e56.
- Lefever S et al. (2009). RDML: Structured language and reporting guidelines for real-time quantitative PCR data. *Nucleic Acids Res* 37, 2,065–2,069.
- Pfaffl MW (2001). A new mathematical model for relative quantification in real-time RT-PCR. *Nucleic Acids Res* 29, e45.
- Rozen S and Skaletsky HJ (2000). Primer3 on the WWW for general users and for biologist programmers. In *Bioinformatics Methods and Protocols: Methods in Molecular Biology*, Krawetz S, Misener S, eds. (Totowa, NJ: Humana Press), pp. 365–386.
- Vandesompele J et al. (2002). Accurate normalization of real-time quantitative RT-PCR data by geometric averaging of multiple internal control genes. *Genome Biol* 3, research 0034.1–0034.11.
- Zuker M (2003). Mfold web server for nucleic acid folding and hybridization prediction. *Nucleic Acids Res* 31, 3,406–3,415.
- NanoDrop is a trademark of Thermo Fisher Scientific Inc.
- The use of SsoAdvanced Universal Supermixes is covered by one or more of the following U.S. patents and corresponding patent claims outside the U.S.: 5,804,375; 5,994,056; and 6,171,785. The purchase of this product includes a limited, nontransferable immunity from suit under the foregoing patent claims for using only this amount of product for the purchaser's own internal research. No right under any other patent claim and no right to perform commercial services of any kind, including without limitation reporting the results of purchaser's activities for a fee or other commercial consideration, is conveyed expressly, by implication, or by estoppel. This product is for research use only. Diagnostic uses under Roche patents require a separate license from Roche. Further information on purchasing licenses may be obtained from the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.
- Bio-Rad's thermal cyclers and real-time thermal cyclers are covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Numbers 6,767,512 and 7,074,367.



LabChip and the LabChip logo are trademarks of Caliper Life Sciences, Inc. Bio-Rad Laboratories, Inc. is licensed by Caliper Life Sciences, Inc. to sell products using the LabChip technology for research use only. These products are licensed for internal research and development use only in detecting, quantitating, and sizing macromolecules, in combination with microfluidics, where internal research and development use expressly excludes the use of this product for providing medical, diagnostic, or any other testing, analysis, or screening services, or providing clinical information or clinical analysis, in any event in return for compensation by an unrelated party.

Information in this tech note was current as of the date of writing (2009) and not necessarily the date this version (Ver E, 2015) was published.

## Acknowledgements

We are grateful to Dr. Jo Vandesompele for his insightful comments and helpful contributions to this paper.

Visit [bio-rad.com/web/PCRampl](http://bio-rad.com/web/PCRampl) for more information.



**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [www.bio-rad.com](http://www.bio-rad.com) **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 1 877 89 01 **Belgium** 03 710 53 00 **Brazil** 55 11 3065 7550  
**Canada** 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00  
**France** 01 47 95 69 65 **Germany** 49 89 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300  
**Israel** 03 963 6050 **Italy** 39 02 216091 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666  
**New Zealand** 64 9 415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04  
**Singapore** 65 6415 3188 **South Africa** 27 (0) 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 026674 55 05  
**Taiwan** 886 2 2578 7189 **Thailand** 1800 88 22 88 **United Kingdom** 020 8328 2000