

## Convenient Multiplex PCR Assays With iQ™ Multiplex Powermix

Liz Jordan and Naheed Aslam, Bio-Rad Laboratories, Inc.,  
2000 Alfred Nobel Drive, Hercules, CA 94547 USA

### Multiplex Quantitative Studies of Gene Expression

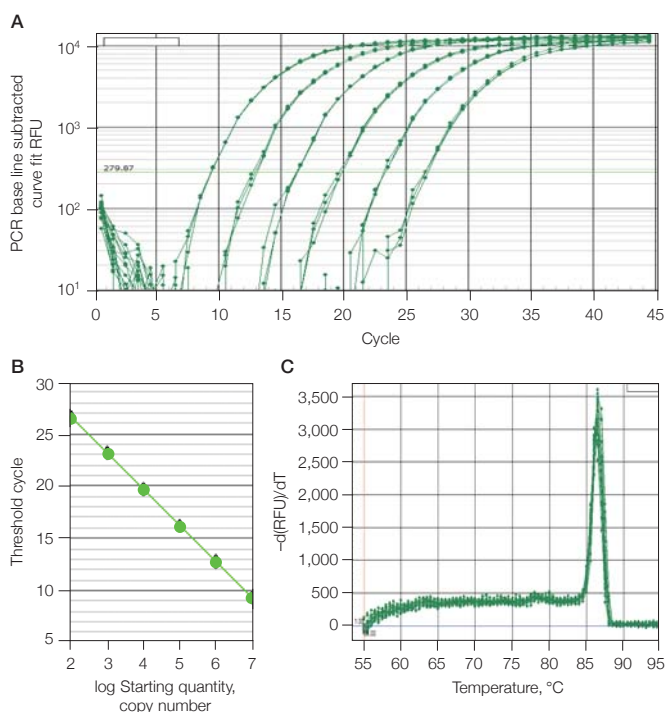
Gene expression analysis is a fundamental element of research on disease and drug development. To study gene expression, the corresponding mRNA levels must be assessed either quantitatively or qualitatively. One reliable method of assessment is reverse transcription of the mRNA into copies of DNA (cDNA) using reverse transcriptase, followed by quantitation of the cDNA using quantitative PCR. Reverse transcription-quantitative PCR (RT-qPCR) has been used for validating expression results obtained from microarrays as well as for more detailed quantitation.

Absolute quantitation of gene expression results in limited information about the transcripts of a specific gene. More relevant information can be obtained from relative gene expression studies in which one looks at internal reference gene(s) within the sample to normalize the relative change in the gene of interest (Bustin 2000, Vandesompele et al. 2002). Determination of relative gene expression can be accomplished by setting up separate reactions to analyze the gene of interest and the internal reference gene (this is required when using SYBR Green I); alternatively, the target genes can be amplified concurrently in a single reaction tube (referred to as multiplexing). In addition to gene expression analysis, multiplex real-time PCR has been used to detect different bacteria in the same sample (Templeton et al. 2003, 2004, Khanna et al. 2005), to detect mutations in more than one PCR product (Ugozzoli et al. 2002), and to analyze multiple transcripts in the same sample (Persson et al. 2005). The advantage of performing multiplex qPCR assays is that one can assess up to five different PCR products in the same tube (using the iQ™5 multicolor real-time PCR detection system). Thus, it is possible to assess several internal reference genes and genes of interest in the same tube, using much less sample than when each reaction is set up individually.

To multiplex, each PCR product requires a gene-specific primer set and a fluorophore-labeled gene-specific probe along with standard PCR reagents. Until recently, qPCR multiplexing was accomplished by supplementing a real-time 2x supermix with additional dNTPs, MgCl<sub>2</sub>, and Taq polymerase, so that limiting

reagents would not influence any of the reactions. Another approach sometimes used in optimizing multiplex reactions is to determine the limiting primer concentration for the more highly expressed targets by testing a series of concentrations. This approach conserves reagents so that amplification of lower-abundance targets can occur unhindered. This practice not only is time consuming, but also uses more of the required reagents. A more practical approach is to use iQ multiplex powermix, which allows multiplexing of up to five unique targets with minimal need for optimizing reaction conditions, allowing considerable savings in cost and time.

This tech note describes several points that must be addressed for optimal multiplexing of reactions.



**Fig. 1. SYBR Green I assay validation and melt curve for the  $\beta$ -actin primer set.** **A**, amplification curve for 10-fold serial dilutions of  $\beta$ -actin plasmid ( $10^7$ – $10^2$  copies per reaction); **B**, resulting standard curve with efficiency = 94.6%, slope =  $-3.459$ ,  $R^2 = 1.00$ ; **C**, melt curve run immediately after amplification.

## Guidelines for Experimental Design

### Primers, Probes, and Internal Reference Genes

Primers and their associated probes should be designed at the same time, but only the primers should be ordered for the initial validation experiments (see next section). It is often advisable to design several primer and probe sets per PCR product, and order and validate the primer sets first.

It is critically important in qPCR multiplex assays that there be no complementarity between any of the primers and probes that will be included in the reaction. During the experimental design stage, it is important to assess all the primer and probe sequences for complementarity. A convenient rule of thumb is to have no more than four complementary bases in a row, to avoid any possible cross-hybridization. Many tools are available to test these interactions and assist in the design of multiplex reactions, including oligonucleotide design web sites and more extensive software packages such as Beacon Designer. This software will perform all pairwise comparisons and will suggest only primers and probes that will not cross-hybridize with each other.

Another important experimental design consideration is the choice of appropriate internal reference genes for normalizing sample input across treatment conditions or samples (Bustin 2000). Several researchers have investigated the use of appropriate housekeeping genes in detail, suggesting that up to three internal reference genes be used to reduce the likelihood of incorporating large errors into the results (Vandesompele et al. 2002). These internal reference genes must be shown to be stably expressed across the proposed experimental conditions (Livak and Schmittgen 2001, Guo et al. 2002).

### Validation of Primer and Probe Specificity and Efficiency

Each real-time reaction should be characterized individually using SYBR Green I before ordering the probe. By doing so, if primers need to be redesigned, one will not incur the added potential cost of probe redesign. Figure 1 shows an example of a reaction run with SYBR Green I to determine reaction efficiency, with melt-curve analysis to show specificity. The melt curve shows the absence of primer-dimers even at low template concentrations. The reaction efficiency is calculated from the slope of the standard curve (expected efficiency >90%). In the case shown, the efficiency was 94.6%.

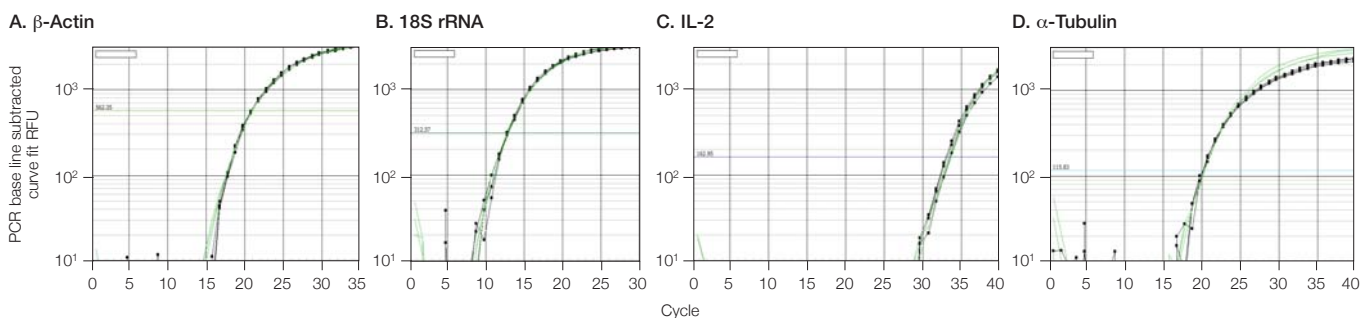
Specificity (amplification of only one PCR product) is verified using melt-curve analysis as well as by running representative samples on a high-resolution gel. The optimal annealing temperature of the primers can be determined using the gradient feature included in iCycler iQ®, MyiQ™, iQ5, and Opticon Monitor™ software. Primer and MgCl<sub>2</sub> concentrations can also be titrated to improve reaction efficiencies. If none of these methods results in adequate reaction optimization, we recommend redesigning the primers.

When both the reaction efficiency and the specificity are acceptable, probes should be ordered and validated over the relevant template concentration range. These assays should display the same performance with the probe as with SYBR Green I (>90% efficiency for each reaction). Because the goal of a multiplex gene expression experiment is to compare expression levels of a gene of interest to a reference gene, it is important that the efficiencies of all reactions be similar to one another (generally within 5%).

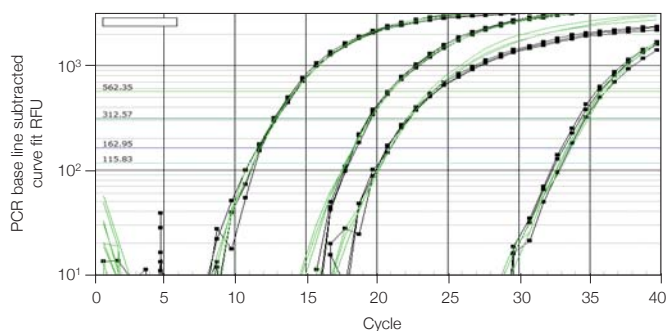
### Multiplex Reactions

Once all individual reactions have been tested, it is important to compare the individual reactions with the multiplex reaction. This comparison will expose any effects of one reaction on the others in the multiplex reaction. There should be no significant difference between the threshold cycle (C<sub>T</sub>) value obtained for a reaction in the singleplex format and the value obtained in the multiplex format. If this is not the case, further investigation of primer interactions will be necessary.

Figures 2 and 3 display a validation experiment showing that the C<sub>T</sub> of each singleplex reaction does not shift when the same assay is performed in a four-target multiplex reaction. Each of the singleplex reactions was previously validated following the strategies described above. Amplification was performed using iQ multiplex powermix and the same primer (300 nM) and probe (200 nM) concentrations for singleplex and multiplex reactions. We recommend running a two-step PCR protocol (combining the annealing and extension steps) and allowing adequate time (about 1 min) for the instrument to collect data through all four or five filter sets.



**Fig. 2. Comparison of singleplex and multiplex results.** Template was human spleen cDNA prepared from purchased RNA, using iScript™ reverse transcriptase. One-tenth of a 1 µg cDNA synthesis reaction was added to each 50 µl reaction. (■), multiplex reactions; (—), singleplex reactions. **A**, β-actin primers and FAM-labeled probe; **B**, 18S rRNA primers with HEX-labeled probe; **C**, interleukin-2 (IL-2) primers and Texas Red-labeled probe; **D**, α-tubulin primers with Cy5-labeled probe.



**Fig. 3. Multiplex reactions showing all four fluorophore traces on the same graph.** From left to right: 18S rRNA (HEX),  $\beta$ -actin (FAM),  $\alpha$ -tubulin (Cy5), and IL-2 (Texas Red). Data set is the same as in Figure 2.

These data demonstrate that each of the singleplex reactions performs similarly whether run independently or in the four-target multiplex assay. Additionally, the difference in  $C_T$  values of 20.6 between the 18S rRNA (HEX) and IL-2 (Texas Red) traces (Figure 3) represents over 6 orders of magnitude difference in template concentration, yet the presence of more highly expressed targets had no effect on the amplification of IL-2. From these data, we have determined that multiplexing these reactions will not skew our data on the relative expression of one gene compared with the others. Furthermore, no additional optimization was required when using iQ multiplex powermix.

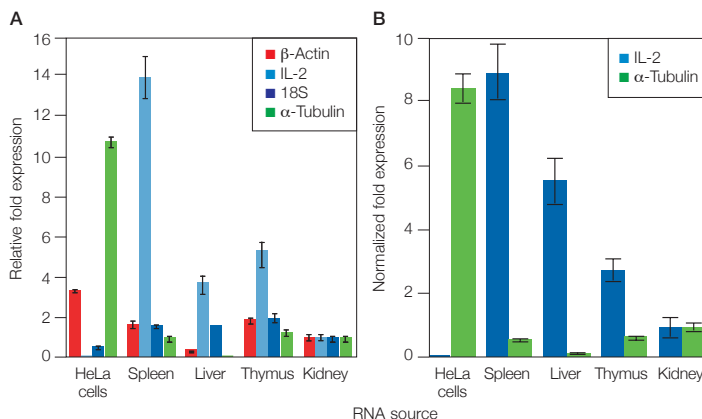
Table 1 shows the average  $C_T$  values obtained for four replicate amplifications using each primer/probe set in singleplex and corresponding multiplex reactions. The  $C_T$  values are essentially the same when including the SD; there is a slight shift of  $\sim 0.12 C_T$  for  $\alpha$ -tubulin, which is acceptable.

**Table 1. Comparison of  $C_T$  of each reaction in singleplex and multiplex.** Values shown are averages  $\pm$  SD.

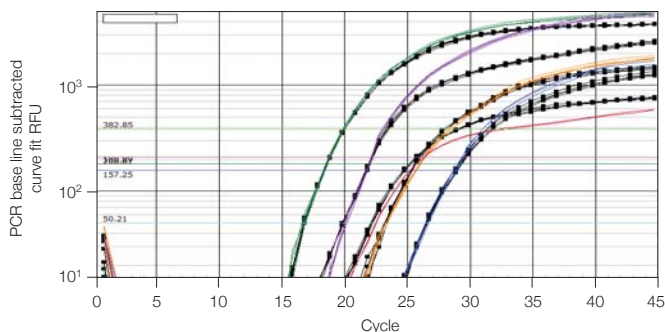
Target	Singleplex	Multiplex
$\beta$ -Actin	20.89 $\pm$ 0.07	20.82 $\pm$ 0.08
18S rRNA	12.85 $\pm$ 0.07	12.82 $\pm$ 0.08
IL-2	33.46 $\pm$ 0.18	33.21 $\pm$ 0.30
$\alpha$ -Tubulin	20.27 $\pm$ 0.04	20.09 $\pm$ 0.02

At this point, the validation of our model system has led to an optimized experimental design for gene expression studies. Modifications of the  $2^{-\Delta\Delta C_T}$  method (Livak and Schmittgen 2001, Vandesompele et al. 2002) are most often used to present relative gene expression data. Figure 4 illustrates the type of relative and normalized gene expression graphs that can be displayed with iQ5 software. The data suggest that HeLa cells express about 8 times more  $\alpha$ -tubulin than kidney, and spleen expresses about 8 times more IL-2 than kidney.

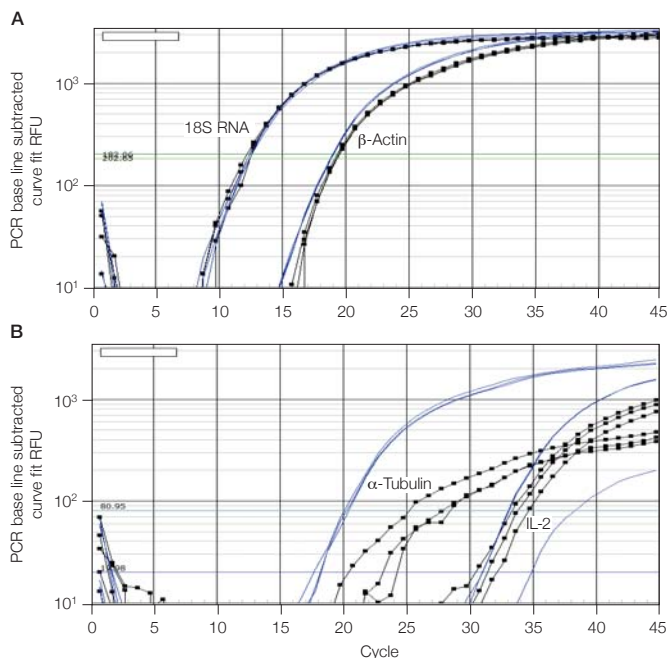
iQ multiplex powermix permits multiplexing of five PCR products in a single tube (Figure 5). Singleplex assays run alongside the five-color multiplex assay verified that there was no significant difference in  $C_T$  values between the two formats. In contrast, when another commercially available supermix designed for multiplex qPCR was used, there was a shift in  $C_T$  values between singleplex and multiplex reactions for targets expressed at low levels (Figure 6B). Therefore, the reaction would require further optimization. Analysis of this suboptimal multiplex data might cause one to conclude that  $\alpha$ -tubulin is expressed over 100-fold lower relative to singleplex data, and normalized to  $\beta$ -actin and 18S rRNA expression levels.



**Fig. 4. Four-color multiplex reactions carried out using cDNAs prepared from different RNA sources.** Reactions were carried out in triplicate with one-tenth of a 1  $\mu$ g cDNA synthesis reaction using total human RNA in each 50  $\mu$ l reaction. Genes analyzed were  $\beta$ -actin, 18S rRNA, IL-2, and  $\alpha$ -tubulin. Kidney was set as the control sample. **A**, relative quantity ( $\Delta C_T$ ). **B**, normalized expression data ( $\Delta\Delta C_T$ ) using  $\beta$ -actin and 18S as reference genes.



**Fig. 5. Five-target multiplex reaction.** One-tenth of a 1  $\mu$ g cDNA synthesis reaction of human spleen total RNA was used in each 50  $\mu$ l reaction. ( $\blacksquare$ ), multiplex; singleplex reactions: ( $\bullet$ ), FAM-labeled  $\beta$ -actin probe; ( $\blacksquare$ ), Cy5-labeled  $\alpha$ -tubulin probe; ( $\blacktriangle$ ), TET-labeled GAPDH probe; ( $\blacklozenge$ ), TAMRA-labeled cyclophilin probe; ( $\blacktriangleleft$ ), Texas Red-labeled IL-2 probe.



**Fig. 6. Four-target multiplex reactions with another supplier's reagent are not as robust for low-expression targets.** One-tenth of a 1  $\mu$ g cDNA synthesis reaction was used in each 50  $\mu$ l reaction. Genes analyzed were  $\beta$ -actin (FAM), 18S rRNA (HEX),  $\alpha$ -tubulin (Cy5), and IL-2 (Texas Red), and the experiment was performed according to the supplier's protocol. (■), multiplex; (□), singleplex reactions. The four traces are plotted in two panels for clarity: Higher-expressing targets are shown in **A**, lower-expressing targets in **B**.

## Conclusions

Multiplexing requires some initial optimization of individual reactions. Once the assays have been validated, multiplexing is a powerful tool that can provide a large amount of data from small sample quantities. The versatility of iQ multiplex powermix reduces the need to optimize multiplex reaction conditions.

## References

- Bustin SA, Absolute quantification of mRNA using real-time reverse transcription polymerase chain reaction assays, *J Mol Endocrinol* 25, 169–193 (2000)
- Guo D et al., The iCycler iQ detection system for evaluating reference gene expression in normal human tissue, *Bio-Rad bulletin* 2804 (2002)
- Khanna M et al., The pneumoplex assays, a multiplex PCR-enzyme hybridization assay that allows simultaneous detection of five organisms, *Mycoplasma pneumoniae*, *Chlamydia (Chlamydothila) pneumoniae*, *Legionella pneumophila*, *Legionella micdadei*, and *Bordetella pertussis*, and its real-time counterpart, *J Clin Microbiol* 43, 565–571 (2005)
- Livak KJ and Schmittgen TD, Analysis of relative gene expression data using real-time quantitative PCR and the  $2^{-\Delta\Delta C_t}$  method, *Methods* 25, 402–408 (2001)
- Persson K et al., Four-color multiplex reverse transcription polymerase chain reaction — overcoming its limitations, *Anal Biochem* 344, 33–42 (2005)
- Templeton KE et al., Rapid and sensitive method using multiplex real-time PCR for diagnosis of infections by influenza A and influenza B viruses, respiratory syncytial virus, and parainfluenza viruses 1, 2, 3, and 4, *J Clin Microbiol* 42, 1564–1569 (2004)
- Templeton KE et al., Development and clinical evaluation of an internally controlled, single-tube multiplex real-time PCR assay for detection of *Legionella pneumophila* and other *Legionella* species, *J Clin Microbiol* 41, 4016–4021 (2003)
- Ugozzoli LA et al., Fluorescent multicolor multiplex homogeneous assay for the simultaneous analysis of the two most common hemochromatosis mutations, *Anal Biochem* 307, 47–53 (2002)
- Vandesompele J et al., Accurate normalization of real-time quantitative RT-PCR data by geometric averaging of multiple internal control genes, *Genome Biol* 3, research0034.1–0034.11 (2002)

Notice regarding Bio-Rad thermal cyclers and real-time systems.

Purchase of this instrument conveys a limited non-transferable immunity from suit for the purchaser's own internal research and development and for use in applied fields other than Human In Vitro Diagnostics under one or more of U.S. Patents Nos. 5,656,493, 5,333,675, 5,475,610 (claims 1, 44, 158, 160–163, and 167 only), and 6,703,236 (claims 1–7 only), or corresponding claims in their non-U.S. counterparts, owned by Applied Biosystems. No right is conveyed expressly, by implication or by estoppel under any other patent claim, such as claims to apparatus, reagents, kits, or methods such as 5' nuclease methods. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

Bio-Rad's real-time thermal cyclers are licensed real-time thermal cyclers under Applied's United States Patent No. 6,814,934 B1 for use in research and for all other fields except the fields of human diagnostics and veterinary diagnostics.

Beacon Designer is a trademark of PREMIER Biosoft International. SYBR and Texas Red are trademarks of Molecular Probes, Inc. Cy is a trademark of GE Healthcare.

Information in this tech note was current as of the date of writing (2006) and not necessarily the date this version (rev A, 2006) was published.

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://www.bio-rad.com) USA 800 4BIORAD Australia 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400  
Canada 905 712 2771 China 86 21 6426 0808 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65  
Germany 089 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 India 91 124 4029300/5013478 Israel 03 963 6050  
Italy 39 02 216091 Japan 03 5811 6270 Korea 82 2 3473 4460 Mexico 55 5200 05 20 The Netherlands 0318 540666 New Zealand 64 9415 2280  
Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 095 721 14 04 Singapore 65 6415 3188 South Africa 27 0861 246 723  
Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189/2578 7241 United Kingdom 020 8328 2000