



PrimePCR™ ASSAYS AND PANELS

PrimePCR Assay

Quick Guide

Real-Time PCR Workflow

For a complete guide to PrimePCR Assays, Panels, and Controls, visit bio-rad.com/PrimePCR to download the instruction manual.

Step 1: Isolate RNA

- Aurum™ Total RNA Mini kit
- Aurum Total RNA Fatty and Fibrous Tissue Kit

Step 2: Synthesize cDNA with reverse transcription (RT) control

- Prepare cDNA with an iScript™ reverse transcription kit
- Add 1 μ l RT control template to each 20 μ l cDNA synthesis reaction

Note: Before first use, suspend the RT control template in 200 μ l nuclease-free TE buffer (pH 7.5) or the same reagent used to dilute the RNA samples. Keep the template on ice.

Step 3: Prepare real-time PCR reaction

1. Thaw and mix reagents.
2. Prepare all quantitative PCR (qPCR) reaction mixes according to Table 1 or 2.
3. Transfer the appropriate volume of the PCR reaction mix into each well.
4. Seal the plate with an appropriate seal and briefly centrifuge.
5. Load the PCR plate into a real-time PCR instrument. Follow the PrimePCR cycling protocol in Table 3.

Table 1. Individual assay (SYBR® Green or probe) reaction setup.

Component	Volume per Reaction		Final Concentration
	96-Well	384-Well	
20x PrimePCR Assay or PrimePCR Control Assay	1 μ l	0.5 μ l	1x
2x SsoAdvanced™ Universal Supermix	10 μ l	5 μ l	1x
cDNA sample	1–4 μ l	0.5–2 μ l	100 ng–100 fg
Nuclease-free water	Variable	Variable	—
Total volume	20 μl	10 μl	—

Table 2. Predesigned or custom plate (SYBR® Green) reaction setup.

Component	Volume per Reaction		Final Concentration
	96-Well	384-Well	
20x PrimePCR Assay or PrimePCR Control Assay*	Dried in well	Dried in well	1x
2x SsoAdvanced™ Universal SYBR® Green Supermix	10 μ l	5 μ l	1x
cDNA sample	1–4 μ l	0.5–2 μ l	100 ng–100 fg
Nuclease-free water	Variable	Variable	—
Total volume	20 μl	10 μl	—

* For the PrimePCR PCR Control Assay, add 1 μ l for 96-well plates or 0.5 μ l for 384-well plates to the well(s) designated for PCR control. The total volume will be 21 μ l in a 96-well plate or 10.5 μ l in a 384-well plate for the PrimePCR PCR Control Assay reaction. Additional volume will not affect the qPCR reaction.

BIO-RAD

Step 4: Perform real-time PCR

Table 3. PrimePCR cycling protocol.

Step	Temperature	Time	Number of Cycles
Activation	95°C	2 min	1
Denaturation	95°C	5 sec	40
Annealing/extension	60°C	30 sec	
Melt curve*	65–95°C (0.5°C increments)	5 sec/step	1

* Melt curve step is for SYBR® Green analysis only.

Step 5: Analyze gene expression data

- CFX Manager™ Software
- PrimePCR Analysis Software
- Analyze PrimePCR Controls according to Table 4

Table 4. Analysis of PrimePCR Controls.

PrimePCR Control	Purpose	Pass, Cq	Fail, Cq
PCR control assay	Tests performance of qPCR reaction with sample	<30	≥30
Reverse transcription control assay	Tests performance of reverse transcription reaction	<30	≥30
DNA contamination control assay	Tests for genomic DNA in sample	≥35	<35
RNA quality assay	Tests RNA integrity in sample	$\Delta Cq \leq 3^*$	$\Delta Cq > 3^*$

* $\Delta Cq = |(RQ2 Cq) - (RQ1 Cq)|$

Cq, quantification cycle; RQ, RNA quality assay.

SYBR is a trademark of Life Technologies Corporation. Bio-Rad Laboratories, Inc. is licensed by Life Technologies Corporation to sell reagents containing SYBR Green I for use in real-time PCR, for research purposes only.

The use of SsoAdvanced 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 these products includes a limited, non-transferable 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. These products are 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.