

## Advancing qPCR Together

Designed for easy performance verification, the CFX qualification plate gives you confidence in your CFX system, allowing you to focus on your research. Simply thaw, spin, and run the predispensed plate for a quick pass/fail result. The CFX qualification plate combines Bio-Rad's high-performance SsoAdvanced™ SYBR® Green supermix with an optimized assay to create a reliable and efficient PCR instrument validation tool.

The CFX qualification plate offers:

- 3-step run start workflow with predispensed plate, predefined thermal cycling protocol, and plate templates
- 1-click pass/fail performance result with CFX Manager™ software
- 2-fold discrimination with 99.7% confidence level and single product melt peak criteria for determining pass/fail results
- Easy incorporation into an instrument qualification procedure with a printable report



For more information, visit us on the Web at [www.bio-rad.com/cfxqualification](http://www.bio-rad.com/cfxqualification).



## Have Confidence in Your CFX System

The CFX qualification plate allows you to easily verify the performance of your CFX system, giving you confidence in your results. The plate contains a five-point standard curve, two unknown groups, and no template controls (NTCs) for validating the optics and thermal block of your CFX system.

When combined with the user-friendly CFX Manager software, you can run the plate in three easy steps and quickly receive pass/fail performance results with the one-click Qualification Plate Check tool in the software (Figure 1). The tool uses an algorithm to determine if the two unknown groups show a 2-fold discrimination with 99.7% confidence level and a single product melt peak (Figure 2).

Use the CFX qualification plate in your performance qualification or preventive maintenance procedures to ensure the operation of your system. With a printable report listing your CFX system's information along with the bar code of the CFX qualification plate, traceability will no longer be a problem.

The CFX qualification plate can be used on any of Bio-Rad's CFX real-time PCR detection systems, including CFX Connect™, CFX96™, CFX96 Touch™, CFX96 Touch Deep Well™, CFX384™, and CFX384 Touch™ systems.

## Ordering Information

Catalog #	Description
184-5098	<b>CFX Qualification Plate</b> , 96-well format, includes 96-well plate containing template DNA, primers, SsoAdvanced™ SYBR® Green supermix, and nuclease-free water, instruction manual
184-5099	<b>CFX Qualification Plate</b> , 384-well format, includes 384-well plate containing template DNA, primers, SsoAdvanced™ SYBR® Green supermix, and nuclease-free water, instruction manual

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.

Bio-Rad's real-time thermal cyclers are licensed real-time thermal cyclers under Applera's U.S. Patent Number 6,814,934 B1 for use in research, human in vitro diagnostics, and all other fields except veterinary diagnostics.

Bio-Rad's 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.

Purchase of SsoAdvanced™ SYBR® Green supermix includes an immunity from suit under patents specified in the product insert to use only the amount purchased for the purchaser's own internal research. No other patent rights are conveyed expressly, by implication, or by estoppel. 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.

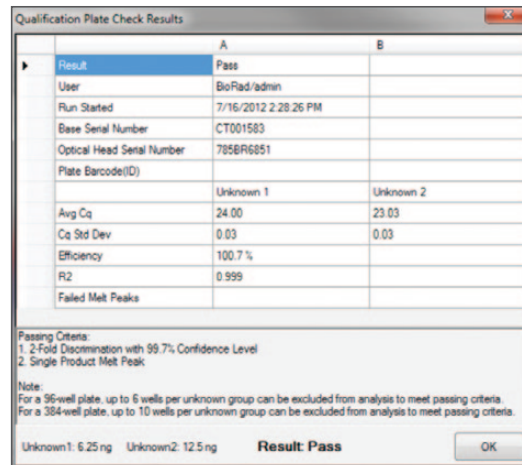


Fig. 1. Qualification Plate Check Results window shows accurate results for easy incorporation into any instrument qualification procedure.

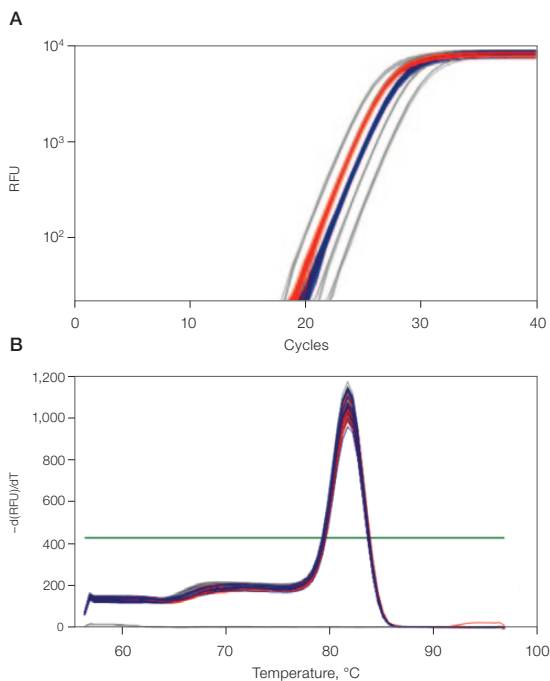


Fig. 2. Amplification (A) and melt peak (B) charts demonstrating a functioning CFX96 Touch real-time PCR detection system. RFU, relative fluorescence units.



**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** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 11 5044 5699  
**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** 089 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** 03 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** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55  
**Taiwan** 886 2 2578 7189 **Thailand** 800 88 22 88 **United Kingdom** 020 8328 2000