

## SsoFast™ EvaGreen® Control Assay

The SsoFast EvaGreen control assay is a pre-designed control kit for real-time PCR. The assay includes control DNA template, SsoFast EvaGreen supermix, primer mix, and nuclease-free water.

- Quickly verifies the performance of your real-time PCR instrument
- Trains new users on real-time PCR setup, instrument operation, and data analysis
- Provides a convenient format with template, pre-designed primers, and a recommended thermal cycling protocol
- Is validated for use on the CFX96™, CFX384™, MiniOpticon™, and MyiQ™2 real-time PCR detection systems



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

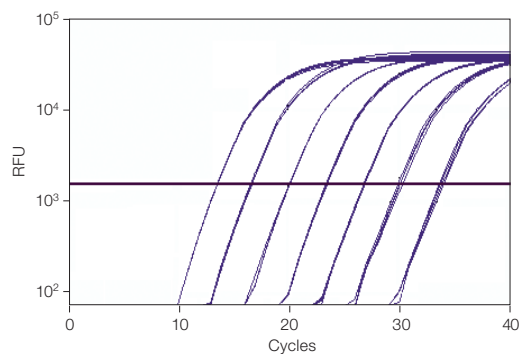
**BIO-RAD**

## Take Control with the SsoFast EvaGreen Control Assay

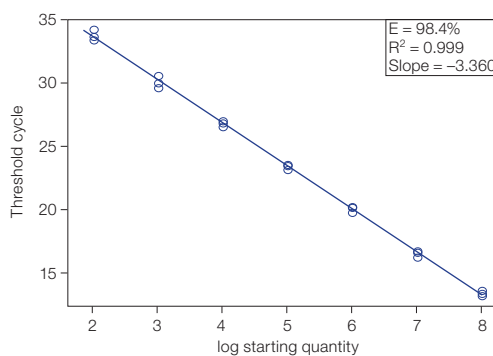
As your research changes, you need to maintain control of your results. The SsoFast EvaGreen control assay allows you to control your experimental conditions and instrument verification and training.

With a reliable assay that is specifically validated for your real-time PCR system, you can easily train new laboratory personnel and confirm that everyone can obtain consistent results. You can verify the consistency of your results from run to run or day to day. You can verify the thermal and optical performance of your real-time PCR system to meet your regulatory needs.

Results obtained using the SsoFast EvaGreen control assay demonstrate high efficiency over a broad dynamic range of  $10^2$ – $10^8$  copies. The assay is validated for use on the CFX96, CFX384, MiniOpticon, and MyiQ2 real-time PCR detection systems. Enough reagents are provided to run a standard curve three times, or up to two hundred 20  $\mu$ l reactions.



Sample amplification plot from a CFX96 real-time PCR detection system. RFU, relative fluorescence units.



Standard curve of  $10^2$ – $10^8$  copies of plasmid DNA template detected on a CFX96 real-time PCR system.

## Ordering Information

Catalog #	Description
172-5996	<b>SsoFast EvaGreen Control Assay</b> , pre-designed kit for verifying performance of real-time PCR systems or training new users, includes template DNA, primers, supermix, and water

EvaGreen is a trademark of Biotium, Inc. Bio-Rad Laboratories, Inc. is licensed by Biotium, Inc. to sell reagents containing EvaGreen dye 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 United States 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.

**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 01 877 89 01 Belgium 09 385 55 11 Brazil 55 31 3689 6600  
Canada 905 364 3435 China 86 20 8732 2339 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 30 210 777 4396 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 0508 805 500  
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 United Kingdom 020 8328 2000