

SsoFast™ EvaGreen® Control Assay Instructions

About the Assay

The SsoFast EvaGreen control assay is a pre-designed kit for training new users on real-time PCR or verifying the performance of your real-time PCR system.

- Demonstrates a broad dynamic range from 10^2 to 10^8 copies
- Validated for use on the CFX96™, CFX384™, MiniOpticon™, and MyiQ™2 real-time PCR detection systems
- Includes instructions for setting up a dilution series to create a standard curve

Storage and Stability

Store at -20°C . After thawing, remaining stock reagents may be stored at 4°C for up to 6 months. Avoid repeated freeze-thaw cycles. Unused master mix and template DNA dilutions should be discarded.

Kit Contents

The SsoFast EvaGreen control assay includes enough reagents to run the dilution series three times (Table 1).

Table 1. SsoFast EvaGreen control assay kit contents.

| Component | Concentration | Volume |
|---------------------------|-------------------------------------|------------------|
| SsoFast EvaGreen supermix | 2x | Two 1 ml tubes |
| Control assay template | 10^8 copies/ μl | 70 μl |
| Control assay primer mix | 100x (50 μM each primer) | 60 μl |
| Nuclease-free water | NA | Two 1.5 ml tubes |

Additional Materials Required

- Sterile, nuclease-free 1.5 ml tubes for dilutions
- Sterile, nuclease-free Tris-EDTA (TE) buffer for dilutions (10 mM Tris, 0.1 mM EDTA, pH 8.0 is recommended)*
- Compatible 48-well, 96-well, or 384-well PCR plate and sealer, as recommended for your real-time PCR detection system
- Pipets and filtered pipet tips to prevent cross-contamination
- UV hood to prevent cross-contamination (optional)

* If TE buffer is not available, use sterile, nuclease-free water.



IF THIS ASSAY WAS SENT TO YOU BY TECHNICAL SUPPORT, PLEASE FOLLOW THE INSTRUCTIONS EMAILED TO YOU.

Preparing the Template DNA Dilutions

1. Pipet 90 μl of TE buffer into seven 1.5 ml tubes labeled 2–8.
2. Prepare the dilution series according to the directions in Table 2. After adding the template DNA to each tube, mix by repeatedly pipeting up and down ten times. Do not vortex the DNA dilutions.
3. Continue with the 96-Well or 384-Well Plate Standard Curve Setup section appropriate for your instrument.

Table 2. Preparation of dilution series.

| Tube | TE Buffer, μl | Template DNA | Final [DNA], copies/ μl |
|------|--------------------------|------------------------------|------------------------------------|
| 1 | — | Control assay template stock | 10^8 |
| 2 | 90 | 10 μl from Tube 1 | 10^7 |
| 3 | 90 | 10 μl from Tube 2 | 10^6 |
| 4 | 90 | 10 μl from Tube 3 | 10^5 |
| 5 | 90 | 10 μl from Tube 4 | 10^4 |
| 6 | 90 | 10 μl from Tube 5 | 10^3 |
| 7 | 90 | 10 μl from Tube 6 | 10^2 |
| 8 | 90 | — | No template control (NTC) |

Note: Tube 1 does not require dilution. Tube 1 is the control assay template stock that comes with the kit.

96-Well Plate Standard Curve Setup

Preparing the Master Mix

1. Thaw all the components. Gentle rocking will help dissolve the SsoFast EvaGreen supermix. Take care not to generate bubbles or foam.
2. Add the components listed in Table 3 to a 1.5 ml tube and mix thoroughly, but gently.

Table 3. Volumes of components for standard curve setup.

| Component | Volume, μl |
|---------------------------------|-----------------------|
| SsoFast EvaGreen supermix (2x) | 650 |
| Control assay primer mix (100x) | 13 |
| Nuclease-free water | 572 |
| Total volume of master mix | 1,235 |

Preparing the Reactions

1. Label eight 1.5 ml tubes 1–8.
2. Pipet 133 μl of the master mix into each tube.
3. Pipet 7 μl of the corresponding template DNA dilution into each tube.
4. Mix contents thoroughly by vortexing gently.
5. Briefly centrifuge the tubes.

Pipeting the Reactions into a Plate

1. Pipet four replicates of each reaction into a 96-well plate using 20 μl per well (Figure 1). Start pipeting the lowest concentration first, ending with the highest concentration. A 100 μl electronic repeat pipeter is helpful, but is not required.
2. Completely seal the plate with a compatible sealer.
3. Centrifuge the plate to settle its contents.
4. Run the plate using the protocol specified in the Thermal Cycling Conditions section.

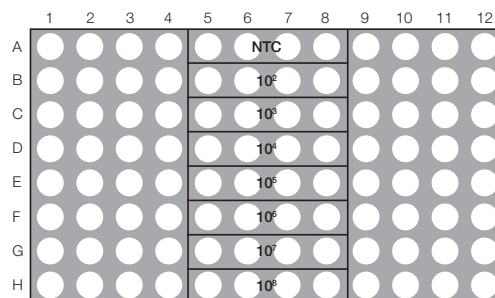


Fig. 1. 96-well plate layout — 20 μl /well.

384-Well Plate Standard Curve Setup

Preparing the Master Mix

1. Thaw all the components. Gentle rocking will help dissolve the SsoFast EvaGreen supermix. Take care not to generate bubbles or foam.
2. Add the components listed in Table 4 to a 1.5 ml tube and mix thoroughly, but gently.

Table 4. Volumes of components for standard curve setup.

| Component | Volume, μ l |
|---------------------------------|-----------------|
| SsoFast EvaGreen supermix (2x) | 450 |
| Control assay primer mix (100x) | 9 |
| Nuclease-free water | 351 |
| Total volume of master mix | 810 |

Preparing the Reactions

1. Label eight 1.5 ml tubes 1–8.
2. Pipet 90 μ l of the master mix into each tube.
3. Pipet 10 μ l of the corresponding template DNA dilution into each tube.
4. Mix contents thoroughly by vortexing gently.
5. Briefly centrifuge the tubes.

Pipeting the Reactions into a Plate

1. Pipet six replicates of each reaction into a 384-well plate using 10 μ l per well (Figure 2). Start pipeting the lowest concentration first, ending with the highest concentration. A 100 μ l electronic repeat pipeter is helpful, but is not required.
2. Completely seal the plate with a compatible sealer.
3. Centrifuge the plate to settle its contents.
4. Run the plate using the protocol specified in the Thermal Cycling Conditions section.

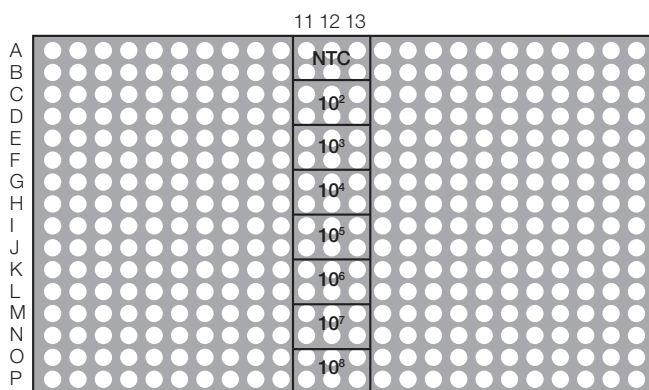


Fig. 2. 384-well plate layout — 10 μ l/well.

Thermal Cycling Conditions

1. Program the real-time PCR instrument using the appropriate plate layout (Figure 1 or 2) and the appropriate thermal cycling protocol for your instrument (Table 5).
2. Place the plate in the real-time PCR instrument and run the program.

Table 5. Thermal cycling protocols for real-time PCR detection systems.

| CFX96 and CFX384 Protocol | MyiQ2 Protocol | MiniOpticon Protocol |
|---------------------------------------------------|----------------------------------------------------|----------------------------------------------------|
| 95°C for 30 sec | 95°C for 30 sec | 95°C for 30 sec |
| 95°C for 1 sec, then 60°C for 5 sec (x 40 cycles) | 95°C for 1 sec, then 60°C for 10 sec (x 40 cycles) | 95°C for 5 sec, then 60°C for 10 sec (x 40 cycles) |
| Default melt curve from 60–95°C | Melt curve from 60–95°C, 30 sec hold | Melt curve from 60–95°C, 5 sec hold |

Analyzing the Data

1. Analyze the data using the default automatic baseline option.
2. With the fluorescence data in log scale view, adjust the threshold such that it is in the linear range of the amplification traces and the reaction efficiency is relatively insensitive to the exact threshold setting (Figure 3).
3. Note the efficiency and R^2 value of the standard curve.

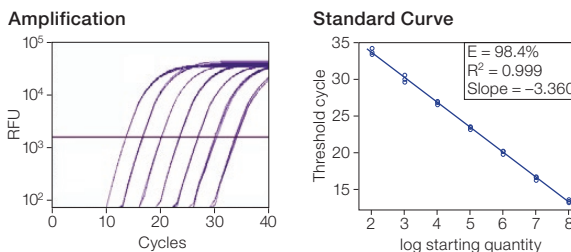


Fig. 3. Sample data for a standard curve on a CFX96 real-time PCR detection system. RFU, relative fluorescence units.

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.

Purchase of this product 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.

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
Laboratories, Inc.**

Life Science
Group

Web site 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