

## SsoFast™ EvaGreen® Supermix

Bio-Rad introduces its next generation of real-time PCR supermixes using our patented\* Sso7d fusion protein technology, delivering a reagent that provides superior qPCR performance in a variety of applications.

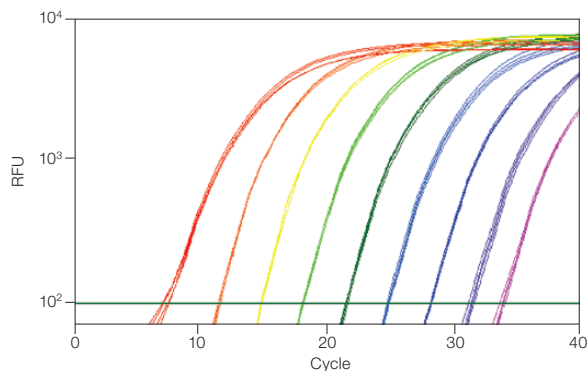
- Unique Sso7d fusion polymerase and optimized buffer deliver unrivaled speed, performance, and tolerance to many common PCR inhibitors
- Minimal PCR inhibition by EvaGreen ensures maximum efficiency, sensitivity, and reproducibility, while generating higher fluorescence compared to SYBR® Green
- Instant polymerase activation and rapid polymerization kinetics for fast qPCR results in less than 30 minutes

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

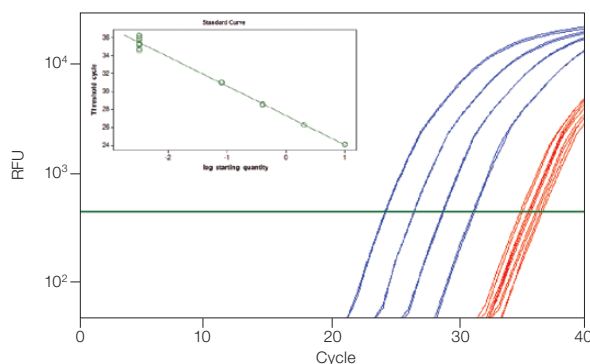
**BIO-RAD**

## SsoFast EvaGreen Supermix

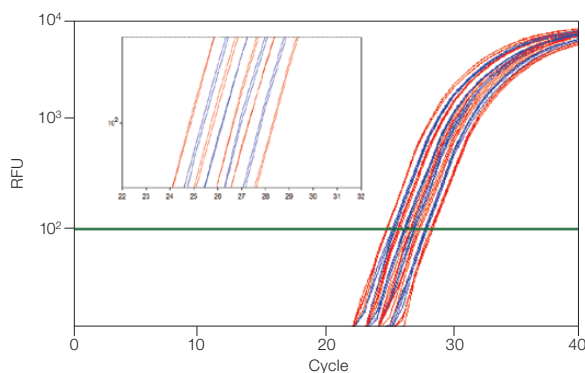
SsoFast EvaGreen supermix is the first member of Bio-Rad's next-generation family of high-performance, real-time PCR reagents. This supermix uses patented\* Sso7d fusion protein technology to deliver excellent performance in a wide range of qPCR applications. By combining a novel engineered hot-start fusion polymerase with optimized buffer and EvaGreen dye, robust qPCR results can be generated in less time and with increased reliability and sensitivity.



The unique fusion polymerase in SsoFast EvaGreen supermix delivers extreme speed and generates exceptional qPCR results in less than 30 min. Tenfold serial dilutions of 10 ng to 100 ag of cDNA from human spleen were used in each 20  $\mu$ l reaction to detect 18S rRNA. 18S rRNA efficiency = 101.8%,  $r = 0.997$ . Total qPCR run time = 29 min. RFU, relative fluorescence units.



SsoFast EvaGreen supermix provides extreme sensitivity in detection of a single copy of target gene. The *ZAP70* gene was amplified and detected from 5-fold serial dilutions of 10 ng to 80 pg (■) and 3.2 pg (■) of human genomic DNA. *ZAP70* efficiency = 102.7%,  $r = 0.991$ . Insert shows the standard curve for the various dilutions. RFU, relative fluorescence units.



Exceptional reproducibility can be achieved with SsoFast EvaGreen supermix. Efficient discrimination and reliable quantification can be obtained from 1.33-fold serial dilutions of input template. The *CBP* gene was amplified from varying amounts of human genomic DNA (5 ng to 500 pg). From left to right: (■) 5 ng, 2.83 ng, 1.60 ng, 903 pg, and 511 pg; (■) 3.76 ng, 2.13 ng, 1.20 ng, and 679 pg. *CBP* efficiency = 96.5%,  $r = 0.996$ . Insert is a magnified view showing robust discrimination and reproducible amplification. RFU, relative fluorescence units.

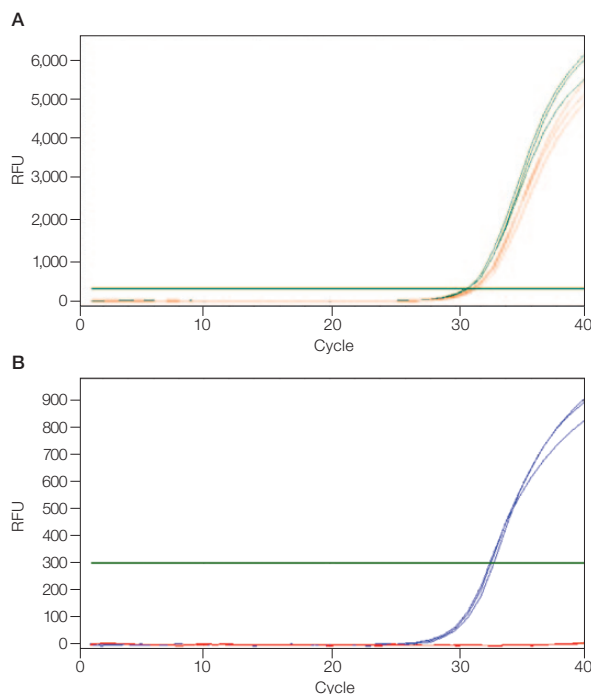
\* U.S. patents 6,627,424; 7,541,170; and 7,560,260.

## Ordering Information

Catalog #	Description
172-5200	SsoFast EvaGreen Supermix, 200 x 20 $\mu$ l reactions, 2x mix contains dNTPs, Sso7d fusion polymerase, MgCl <sub>2</sub> , EvaGreen dye, stabilizers
172-5201	SsoFast EvaGreen Supermix, 500 x 20 $\mu$ l
172-5202	SsoFast EvaGreen Supermix, 1,000 x 20 $\mu$ l
172-5203	SsoFast EvaGreen Supermix, 20 ml bottle, 2,000 x 20 $\mu$ l

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. SYBR is a trademark of Molecular Probes, Inc.

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.



SsoFast EvaGreen supermix demonstrates superior inhibitor tolerance and direct qPCR capability from cell culture. **A**, efficient amplification and detection of a spike-in control template in the absence (■) or presence (■) of conditioned tissue culture medium using SsoFast EvaGreen supermix; **B**, competitor's standard qPCR reagent is able to amplify the spike-in control template only in the absence (■) vs. presence (■) of culture medium. 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 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