

SsoFast™ EvaGreen® Supermix with Low ROX

Bio-Rad expands its line of SsoFast supermixes, delivering a reagent capable of superior performance on ROX-dependent real-time PCR systems.

- Convenient 1-tube formulation, preblended with ROX for inter-well signal correction on ABI 7500 and Stratagene Mx series real-time PCR systems
- Unique fusion polymerase and optimized buffer deliver unrivaled speed and performance for a variety of qPCR applications
- Instant polymerase activation and rapid polymerization kinetics for faster qPCR results in less than 40 min

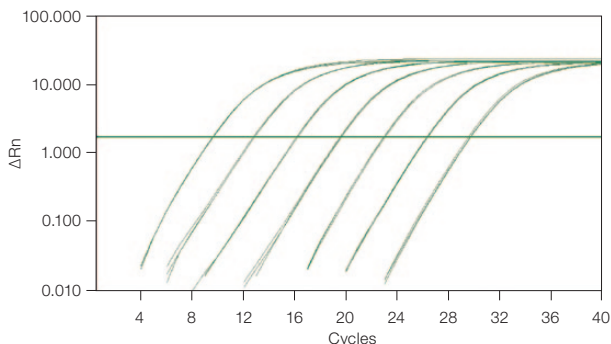
For more information, visit us on the Web at www.bio-rad.com/supermixes.



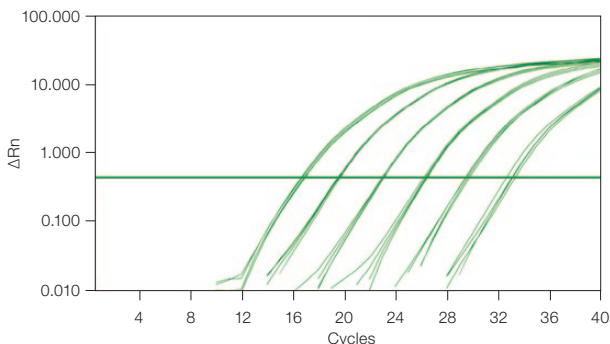
BIO-RAD

SsoFast EvaGreen Supermix with Low ROX

SsoFast EvaGreen supermix with low ROX is part 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 EvaGreen dye and an optimized buffer, robust qPCR results can be generated in less time and with increased speed, reliability, and sensitivity.



SsoFast EvaGreen supermix with low ROX generates linear results over 6 orders of dynamic range on the ABI 7500 fast real-time PCR system. Tenfold serial dilutions of 100 ng to 100 fg of cDNA from HeLa total RNA were used in each 20 μ l reaction to detect 18S rRNA. 18S rRNA efficiency = 98.6%, $r^2 = 0.999$.



The unique fusion polymerase in SsoFast EvaGreen supermix with low ROX generates exceptional qPCR results on the ABI 7500 fast real-time PCR system. Tenfold serial dilutions from 10 ng to 100 fg of cDNA from human spleen were used in each 20 μ l reaction to detect α -tubulin. α -tubulin efficiency = 105.8%, $r = 0.996$. **Total qPCR run time = 39 min** (not including melt curve).

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

Ordering Information

Catalog #	Description	Catalog #	Description
172-5210	SsoFast EvaGreen Supermix with Low ROX , 200 x 20 μ l reactions, 2x mix contains dNTPs, Sso7d fusion polymerase, MgCl ₂ , ROX passive reference dye, stabilizers	172-5212	SsoFast EvaGreen Supermix with Low ROX , 1,000 x 20 μ l reactions
172-5211	SsoFast EvaGreen Supermix with Low ROX , 500 x 20 μ l reactions	172-5213	SsoFast EvaGreen Supermix with Low ROX , 2,000 x 20 μ l reactions, 20 ml bottle

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. Mx is a trademark of Stratagene Corporation. ROX is a trademark of Applied Biosystems Corporation.

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
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 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** 30 210 777 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 **Malaysia** 60 3 2117 5260 **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 3170 **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** 66 2 6518311 **United Kingdom** 020 8328 2000