

SsoFast™ Probes Supermix

Bio-Rad expands its next-generation line of SsoFast supermixes, delivering a reagent that provides superior qPCR performance for gene expression analysis.

- Unique fusion polymerase and optimized buffer deliver unrivaled speed and performance
- Robust, simultaneous detection of up to 2 different gene targets
- Instant polymerase activation and rapid polymerization kinetics for results in less than 30 min
- Compatible with fast and standard cycling conditions under a broad range of primer/probe concentrations

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

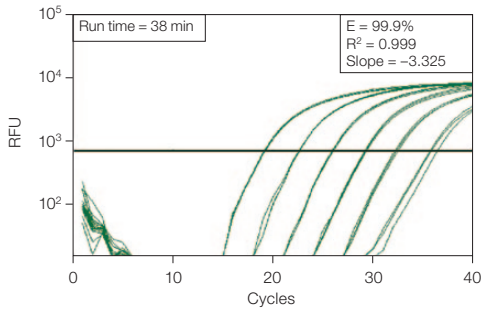


BIO-RAD

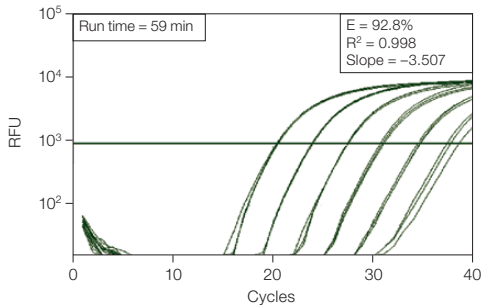
SsoFast Probes Supermix

SsoFast probes supermix 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 an optimized buffer, robust qPCR results can be generated in less time and with increased speed, reliability, and sensitivity.

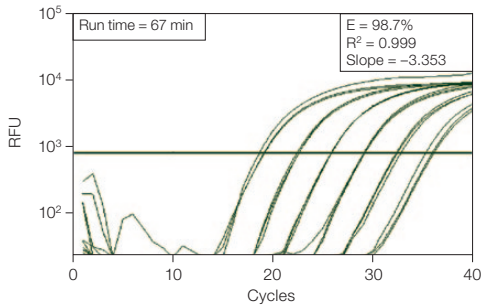
A. Fast 2-Step Cycling



B. Standard 2-Step Cycling



C. Standard 3-Step Cycling



SsoFast probes supermix delivers robust results in a broad range of cycling conditions. Tenfold serial dilutions of 100 ng to 1 pg of cDNA from HeLa cells were used in each 20 µl reaction to detect α -tubulin. RFU, relative fluorescence units.

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

Ordering Information

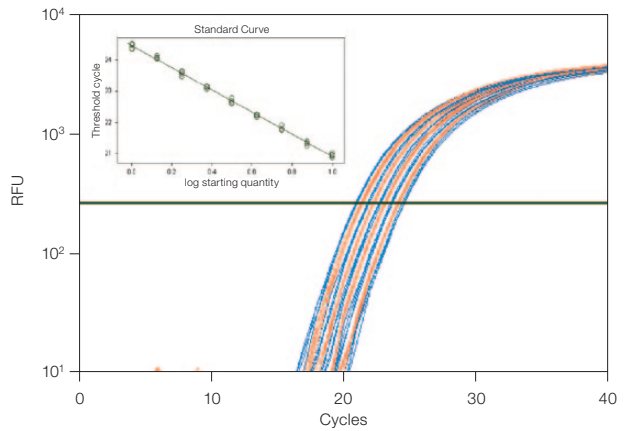
Catalog #	Description
172-5230	SsoFast Probes Supermix , 200 x 20 µl reactions, 2x mix contains dNTPs, Sso7d fusion polymerase, MgCl ₂ , stabilizers
172-5231	SsoFast Probes Supermix , 500 x 20 µl reactions
172-5232	SsoFast Probes Supermix , 1,000 x 20 µl reactions
172-5233	SsoFast Probes Supermix , 2,000 x 20 µl reactions, 20 ml bottle

Texas Red is a trademark of Invitrogen Corporation.

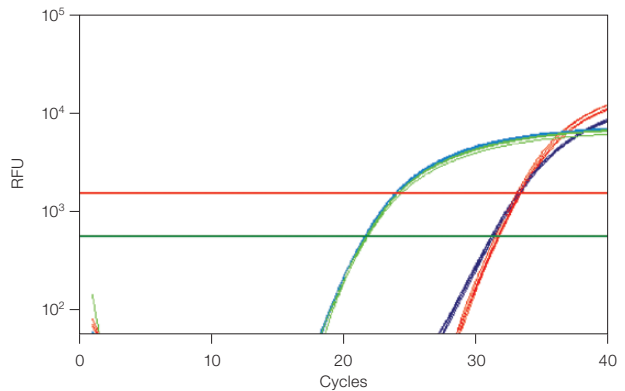
Purchase of this product includes an immunity from suit under patents specified in the product insert to use only the amount purchased for the purchaser's own internal research. No other patent rights are conveyed expressly, by implication, or by estoppel. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

Bio-Rad'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.



Exceptional reproducibility can be achieved on the CFX384™ real-time PCR detection system with SsoFast probes supermix. Efficient discrimination and reliable quantification can be obtained from 1.33-fold serial dilutions of input template. The *GAPDH* gene was amplified from varying amounts of HeLa cDNA (1 ng to 100 pg). From left to right: (■) 1 ng, 565 pg, 320 pg, 181 pg, and 102 pg; (■) 752 pg, 425 pg, 240 pg, and 136 pg. *GAPDH* efficiency = 91.5%, R² = 0.997. Inset shows the standard curve for the various dilutions. **Total qPCR run time = 50 min.** RFU, relative fluorescence units.



SsoFast probes supermix delivers superior results for gene expression analysis of two targets on the CFX96™ real-time PCR detection system, with no difference in detection of a low-expressing gene in duplex or simplex. cDNA from human liver (100 ng) was used in each 20 µl reaction. (■) HEX-labeled *GAPDH* probe duplex reaction; (■) Texas Red-labeled IL-2 probe duplex reaction; (■) HEX-labeled *GAPDH* probe simplex reaction; (■) Texas Red-labeled IL-2 probe simplex reaction. **Total qPCR run time = 38 min.** RFU, relative fluorescence units.



**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