

SsoFast™ Probes Supermix with ROX

SsoFast probes supermix with ROX is a ROX-blended version of the high-performing SsoFast probes supermix. It employs patented* Sso7d fusion protein technology for use in probe-based real-time quantitative PCR assays on **Applied Biosystems 7900HT** and **7900HT fast real-time PCR systems**.

- **Fast qPCR results** — Sso7d fusion polymerase's stringent hot-start and rapid polymerization kinetics provide fast results in less than 30 min
- **Universal conditions** — fully compatible with your pre-existing cycling protocols and primer/probe conditions without the need for additional optimization
- **Minimal PCR inhibition** — highly tolerant polymerase exhibits low PCR inhibition for better reproducibility
- **Distinct target detection** — robust, simultaneous detection of up to 2 different gene targets using fluorogenic probes under fast qPCR conditions



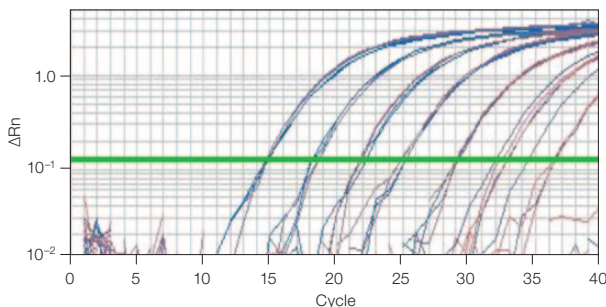
For more information, visit us on the Web at www.bio-rad.com/supermixes.

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

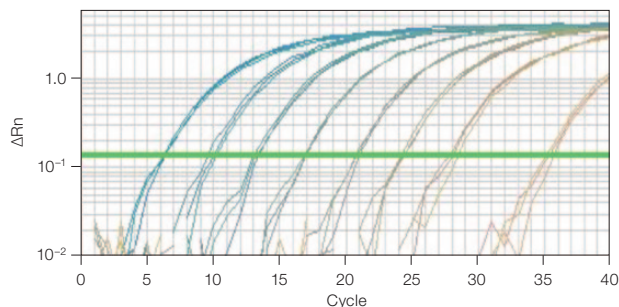


SsoFast Probes Supermix with ROX

SsoFast probes supermix with ROX combines Sso7d fusion protein technology and an optimized formulation to help perform fast duplex qPCR on **Applied Biosystems 7900HT** and **7900HT fast real-time PCR systems** while maintaining a broad dynamic range, reliability, and accurate quantification for robust gene expression analysis.

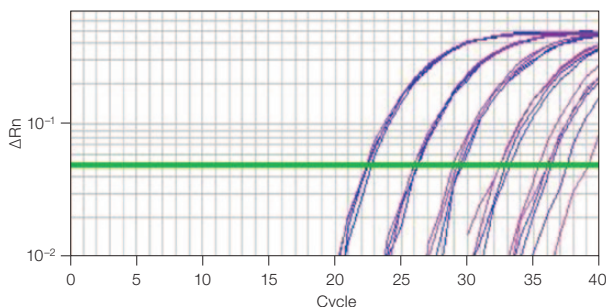


SsoFast probes supermix with ROX demonstrates excellent sensitivity and efficiency for simplex qPCR from cDNA using an ABI 7900HT fast real-time PCR system (384-well standard block). cDNA from HeLa cells (100 ng to 0.1 pg) was used in a 10 μ l reaction. Simplex reaction using a VIC-labeled tubulin probe ($R^2 = 0.9966$, efficiency = 92.2%, slope = -3.53).

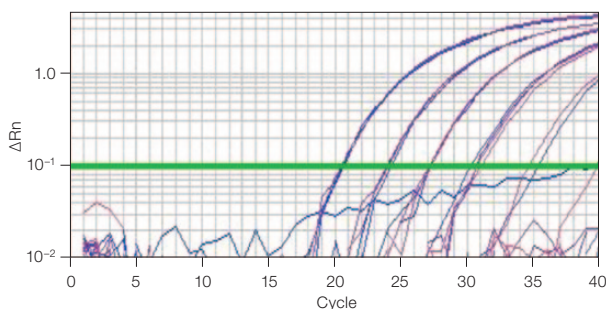


SsoFast probes supermix with ROX demonstrates excellent sensitivity and efficiency for duplex qPCR from targets with approximately a 1,000-fold difference in expression levels using an ABI 7900HT fast real-time PCR system (96-well standard block). cDNA from HeLa cells (100 ng to 0.1 pg) was used in a 20 μ l reaction. Duplex reaction using a FAM-labeled TFRC probe (\blacksquare ; $R^2 = 0.9970$, efficiency = 90.3%, slope = -3.58) and a VIC-labeled 18S probe (\blacksquare ; $R^2 = 0.9990$, efficiency = 88.6%, slope = -3.63).

A



B



SsoFast probes supermix with ROX demonstrates excellent sensitivity and efficiency for simplex qPCR from genomic DNA using an ABI 7900HT fast real-time PCR system (384-well standard block). Human gDNA (250 ng to 2.5 ng) was used in a 10 μ l reaction. **A**, simplex reaction using a TAMRA-labeled IL-1 β probe ($R^2 = 0.9925$, efficiency = 103.9%, slope = -3.23); **B**, simplex reaction using a TET-labeled CBP probe for a 70.9% GC-rich amplicon ($R^2 = 0.9959$, efficiency = 92.7%, slope = -3.51).

Ordering Information

Catalog #	Description
172-5250	SsoFast Probes Supermix With ROX , 200 x 20 μ l reactions, 2x real-time PCR mix contains dNTPs, Sso7d fusion polymerase, MgCl ₂ , ROX passive reference dye, enhancers, stabilizers
172-5251	SsoFast Probes Supermix With ROX , 500 x 20 μ l reactions
172-5252	SsoFast Probes Supermix With ROX , 1,000 x 20 μ l reactions
172-5253	SsoFast Probes Supermix With ROX , 2,000 x 20 μ l reactions, 20 ml bottle

VIC is a trademark of Applied Biosystems.

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

**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 Vietnam 84 8 38131140