

AMPLIFICATION

SsoAdvanced™ SYBR® Green Supermix

- Robust real-time results under any conditions
- Advanced performance with short run times

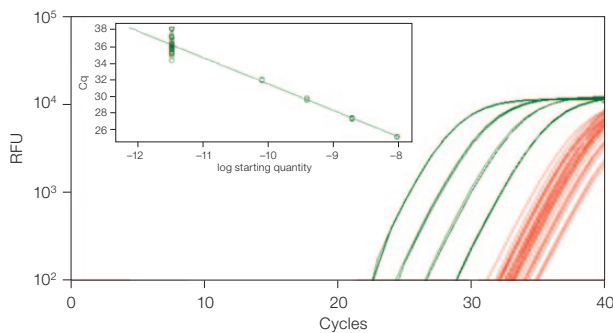
Increased Performance from Your SYBR® Green qPCR

SsoAdvanced™ SYBR® Green supermix is the newest high-performance real-time PCR supermix based on Bio-Rad's patented* Sso7d fusion protein technology. It is formulated for a wide range of qPCR applications. The dsDNA binding protein, Sso7d, stabilizes the polymerase-template complex, providing superior inhibitor tolerance, increased processivity, and greater speed without affecting PCR sensitivity, efficiency, or reproducibility.

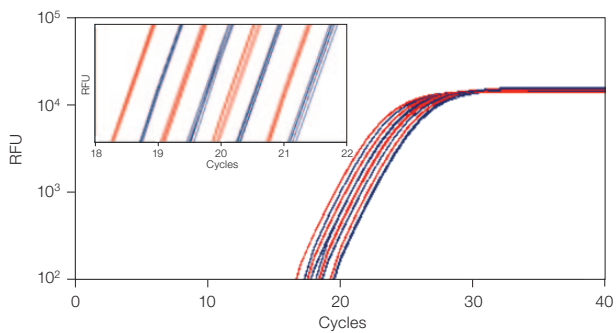
- **Achieve superior gene expression results under various cycling conditions** — robust formulation delivers consistent performance in standard or fast cycling
- **Increase sensitivity and efficiency of detection from compromised samples** — Sso7d fusion polymerase has increased resistance to a wide variety of PCR inhibitors

- **Decrease run times and time to results without compromising qPCR data quality** — Sso7d fusion polymerase and optimized buffer provide rapid polymerization kinetics and instant polymerase activation
- **Obtain better results with predeveloped qPCR assays** — advanced formulation tolerates a broad range of reaction conditions, primer concentrations, and temperature ranges

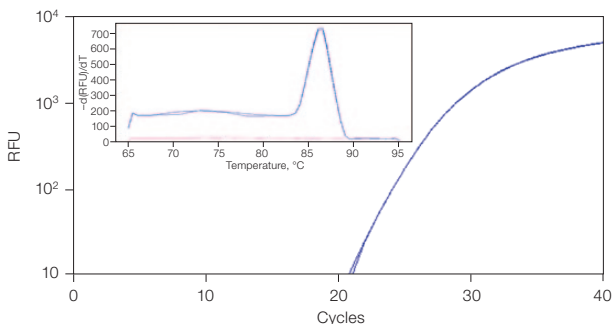




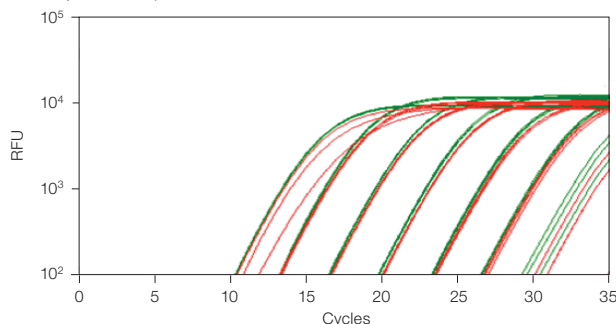
SsoAdvanced™ SYBR® Green supermix provides extreme sensitivity in the detection of a single copy target gene. The cyclin gene was amplified and detected from fivefold serial dilutions of 10 ng–80 pg (■) and 3.2 pg (■) human genomic DNA. Cyclin efficiency = 103%, $R^2 = 1$ (for 10 ng–80 pg). Inset shows the standard curve for the various dilutions. Cq, quantification cycle; RFU, relative fluorescence units.



Exceptional reproducibility can be achieved with SsoAdvanced™ SYBR® Green supermix. Efficient discrimination and reliable quantification can be obtained from a 1.33-fold serial dilution of input template. The *GAPDH* gene was amplified from varying amounts of human genomic DNA (1 ng–136 pg). From left to right: 1 ng, 753 pg, 565 pg, 425 pg, 320 pg, 240 pg, 181 pg, and 136 pg. *GAPDH* efficiency = 96.2%, $R^2 = 0.999$. Inset is a magnified view showing robust discrimination and reproducible amplification (six replicates for each input amount). RFU, relative fluorescence units.



SsoAdvanced™ SYBR® Green supermix demonstrates superior inhibitor tolerance. The *ADAR* gene was amplified from HeLa cDNA in the presence of water alone, or in the presence of a known PCR inhibitor, Eagle's minimal essential medium (EMEM) with fetal bovine serum (FBS; 0, 2.5, 5, 10, and 20%), added to SsoAdvanced™ SYBR® Green supermix (■) or a traditional Taq DNA polymerase–based qPCR master mix (■). SsoAdvanced™ SYBR® Green supermix showed quality amplification in all reactions (EMEM with 20% FBS data shown), while the Taq DNA polymerase–based qPCR master mix failed to amplify in all EMEM with FBS combinations (shown in the inset melt curve). RFU, relative fluorescence units.



Start Time, hr	<i>GAPDH</i> Efficiency	R^2	Slope
0	100.9	0.998	-3.30
24	101.0	0.999	-3.30
48	98.6	0.998	-3.36
72	99.5	0.999	-3.33
144	101.2	0.993	-3.29

SsoAdvanced™ SYBR® Green supermix maintains exceptional stability on the CFX automation system. Tenfold serial dilutions of 100 ng–1 pg cDNA from human spleen were used in each 20 μ l reaction to detect *GAPDH*. All reactions were assembled and loaded onto the CFX automation system and run after varying times (0–144 hr) on the CFX96™ real-time PCR detection system. The thermal cycling conditions were 95°C for 30 sec, then 35 cycles of 95°C for 10 sec and 60°C for 30 sec. Results (*GAPDH* efficiency, R^2 , and slope) for five time points are shown in the table above. The amplification plot shows real-time traces from 0 hr (■) and 144 hr (■). RFU, relative fluorescence units.

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

SYBR is a trademark of Molecular Probes, Inc. Bio-Rad Laboratories, Inc. is licensed by Molecular Probes, Inc. to sell reagents containing SYBR Green I for use in real-time PCR, for research purposes only.

Bio-Rad's real-time thermal cyclers are licensed real-time thermal cyclers under Applied's U.S. 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.

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.



Ordering Information

Catalog #	Description
172-5260	SsoAdvanced™ SYBR® Green Supermix , 200 x 20 μ l reactions, 2 x 1 ml, 2x real-time PCR mix contains dNTPs, Sso7d fusion polymerase, $MgCl_2$, SYBR® Green I, stabilizers
172-5261	SsoAdvanced™ SYBR® Green Supermix , 500 x 20 μ l reactions, 5 x 1 ml
172-5262	SsoAdvanced™ SYBR® Green Supermix , 1,000 x 20 μ l reactions, 10 x 1 ml
172-5264	SsoAdvanced™ SYBR® Green Supermix , 2,500 x 20 μ l reactions, 5 x 5 ml
172-5265	SsoAdvanced™ SYBR® Green Supermix , 5,000 x 20 μ l reactions, 10 x 5 ml



**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