

iQ™ Supermixes for qPCR

Always crowd-pleasers, Bio-Rad's iQ supermixes deliver memorable performances every time in real-time PCR assays.

- iTaq™ hot-start DNA polymerase allows sensitive and accurate detection of as few as 10 copies of template
- 2 unique formulas afford compatibility with any fluorescent detection chemistry
- These robust supermixes provide precise linear detection over 6 orders of magnitude



iQ™ SYBR® Green Supermix

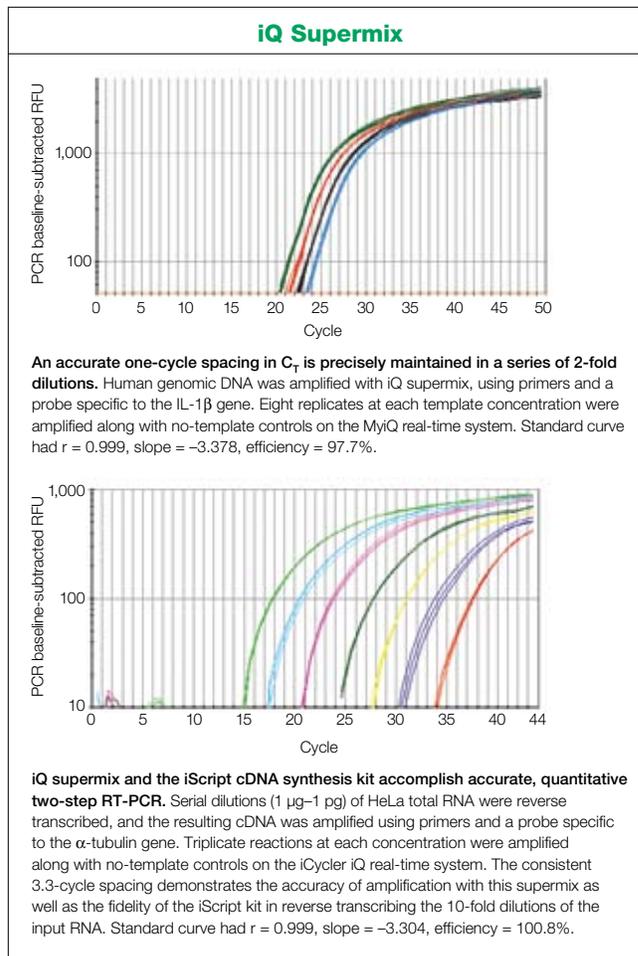
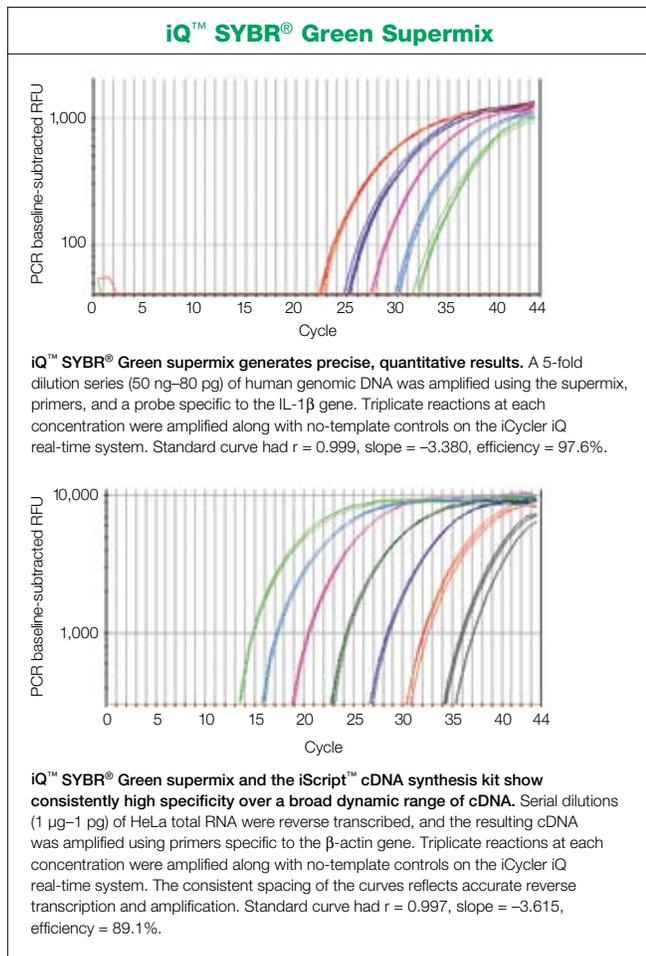
iQ Supermix

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

BIO-RAD

iQ Supermixes for Real-Time PCR

Our versatile iQ supermixes provide the ultimate convenience in preblended solutions for a wide array of real-time PCR applications. The two unique formulas have been specifically optimized for maximum flexibility and performance with any fluorescent detection chemistry, whether through the use of sequence-specific probes or the SYBR® Green I DNA binding dye. Both supermixes contain iTaq DNA polymerase, a highly sensitive hot-start polymerase that minimizes the amplification of nonspecific PCR products. In addition, both supermixes contain buffering agents and qPCR-qualified dNTPs, to afford the most sensitive amplification available. The iQ™ SYBR® Green supermix also contains fluorescein for the collection of well factors on the iCycler iQ®, MyiQ™, and iQ™5 real-time PCR detection systems. With these mixes, you can be assured of optimal qPCR results each and every time.



Ordering Information

Catalog #	Description
170-8862	iQ Supermix , 500 x 50 μl reactions, 2x mix contains 100 mM KCl, 40 mM Tris-HCl, pH 8.4, 0.4 mM each dNTP (dATP, dCTP, dGTP, dTTP), 50 U/ml iTaq DNA polymerase, 6 mM MgCl_2 , stabilizers
170-8864	iQ Supermix , 1,000 x 50 μl reactions
170-8882	iQ™ SYBR® Green Supermix , 500 x 50 μl reactions, 2x mix contains 100 mM KCl, 40 mM Tris-HCl, pH 8.4, 0.4 mM each dNTP (dATP, dCTP, dGTP, dTTP), 50 U/ml iTaq DNA polymerase, 6 mM MgCl_2 , SYBR® Green I, 20 nM fluorescein, stabilizers
170-8884	iQ™ SYBR® Green Supermix , 1,000 x 50 μl reactions
170-8885	iQ™ SYBR® Green Supermix , 2,000 x 50 μl reactions (50 ml bottle)
170-8890	iScript cDNA Synthesis Kit , 25 x 20 μl reactions
170-8891	iScript cDNA Synthesis Kit , 100 x 20 μl reactions

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

Practice of the patented 5' Nuclease Process requires a license from Applied Biosystems. The purchase of these products 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 when used with the separate purchase of Licensed Probe. No other patent rights are conveyed expressly, by implication, or by estoppel. Further information on purchasing licenses may be obtained from 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 Applied's United States Patent No. 6,814,934 B1 for use in research and for all other fields except the fields of human diagnostics and 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 Nos. 6,767,512 and 7,074,367.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA 800 4BIORAD Australia 61 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400 Canada 905 364 3435 China 86 21 6426 0808 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65 Germany 089 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 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