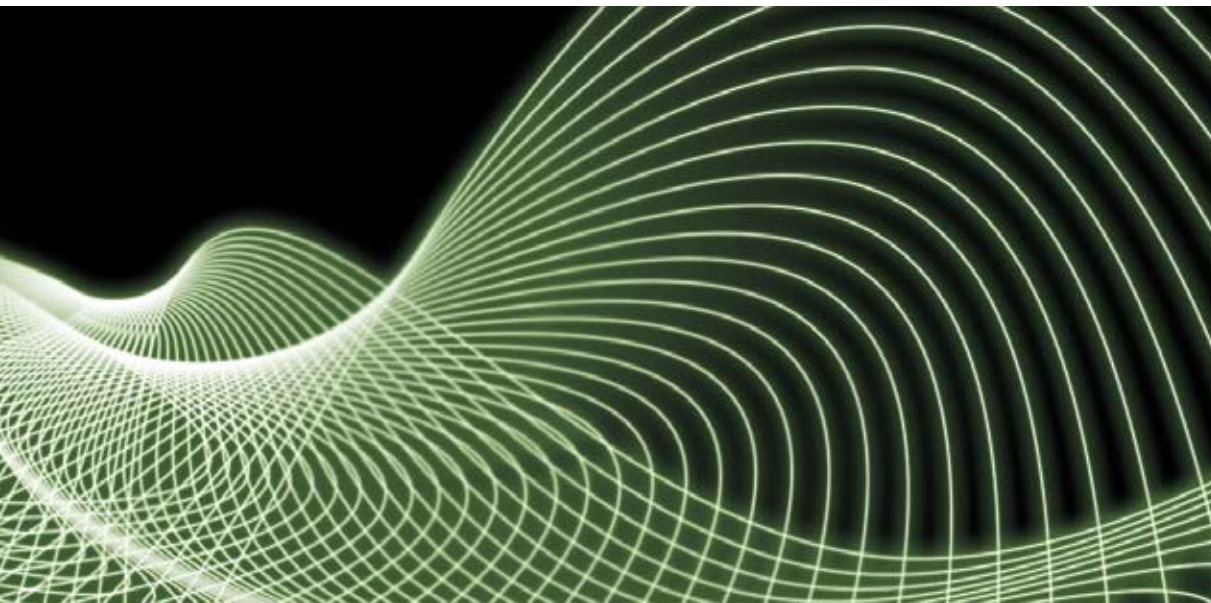




iTaQ™ Supermixes With ROX for qPCR

From the experts in real-time PCR, these supermixes with ROX generate precise, accurate results every time.

- 2 distinct formulas, both validated for use on all ROX-dependent optical thermal cyclers and with any real-time detection chemistry
- Sensitive and accurate detection of low-abundance targets
- Convenient 1-tube formulation preblended with ROX to correct for interwell signal variation

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

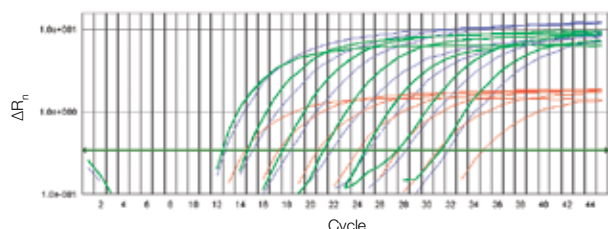


BIO-RAD

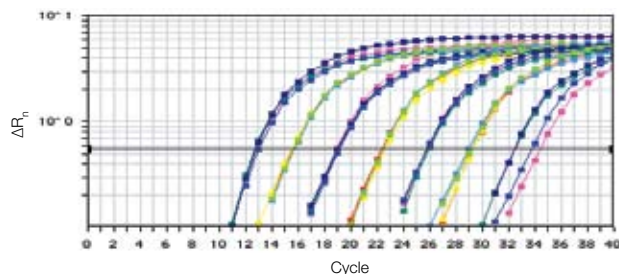
iTaq Supermixes With ROX

The two distinct iTaq supermixes with ROX are formulated to easily achieve optimal results in real-time quantitative PCR assays. These mixes yield high performance with cDNA as well as genomic and plasmid DNA templates over a broad dynamic range, achieving sensitive and specific amplification over at least 7 orders of magnitude. Both supermixes are preblended with hot-start iTaq DNA polymerase, optimized buffer, nucleotides, and ROX passive reference dye. These convenient one-tube formulations will enable you to obtain specific and sensitive amplification every time.

iTaq SYBR Green Supermix With ROX

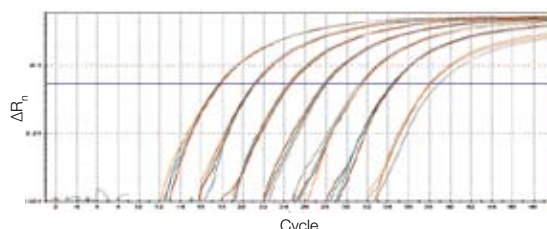


iTaq SYBR Green supermix with ROX gives superior results on the ABI PRISM 7000 sequence detection system. The Bio-Rad supermix (green) or two other ROX-containing supermixes (red and blue) were used to amplify 10-fold serial dilutions (10^7 to 10 copies) of the GAPDH gene from a plasmid template. Reactions were performed in quadruplicate (replicates not shown). The Bio-Rad supermix consistently generated earlier C_T values than the other mixes, and produced high ΔR_n values.

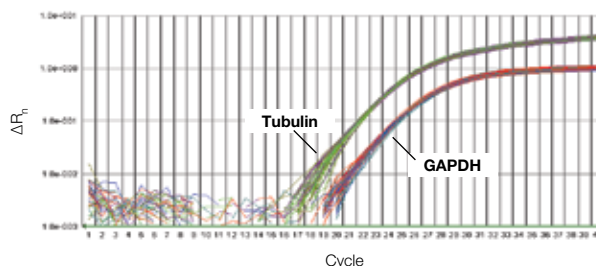


iTaq SYBR Green supermix with ROX performs well over 6 orders of magnitude on the ABI PRISM 7700 sequence detection system. The Bio-Rad supermix was used to amplify 10-fold serial dilutions (10^7 to 10 copies) of the β -actin gene from a plasmid template. Reactions were performed in quadruplicate. Standard curve had $r = 0.996$, slope = -3.40 , efficiency = 96.8%.

iTaq Supermix With ROX



iTaq supermix with ROX generates linear results over 6 orders of magnitude on the Stratagene Mx3000P real-time PCR system. The Bio-Rad supermix was used to amplify 10-fold serial dilutions of 100 ng to 100 fg of cyclophilin cDNA from HeLa RNA. An Applied Biosystems PDAR probe (FAM/MGB) was used for detection. Standard curve had $r = 0.999$, slope = -3.402 , efficiency = 96.8%.



iTaq supermix with ROX produces highly uniform results in a duplex reaction on the ABI PRISM 7000 sequence detection system. The Bio-Rad supermix was used to amplify tubulin and GAPDH targets from 1 ng of cDNA. A FAM-labeled probe was used for tubulin detection, while an Applied Biosystems PDAR probe (VIC/MGB) was used to detect GAPDH. In both dye layers, the Bio-Rad mix generated highly uniform results. The tubulin gene target had an average C_T of 22.83 and a standard deviation of 0.130 over the 47 replicates tested; the GAPDH gene target had an average C_T of 26.01 and standard deviation of 0.145.

Ordering Information

Catalog #	Description
172-5850	iTaq SYBR Green Supermix With ROX, 200 x 50 μ l reactions, 2x mix contains 0.4 mM each of dATP, dCTP, dGTP, and dTTP, 50 U/ml iTaq DNA polymerase, 6 mM Mg^{2+} , SYBR Green I, ROX reference dye, stabilizers
172-5851	iTaq SYBR Green Supermix With ROX, 1,000 x 50 μ l reactions
172-5852	iTaq SYBR Green Supermix With ROX, 2,000 x 50 μ l reactions (50 ml bottle)
172-5854	iTaq Supermix With ROX, 200 x 50 μ l reactions, 2x mix contains 0.4 mM each of dATP, dCTP, dGTP, and dTTP, 50 U/ml iTaq DNA polymerase, 10 mM Mg^{2+} , ROX reference dye, stabilizers
172-5855	iTaq Supermix With ROX, 500 x 50 μ l reactions
172-5856	iTaq Supermix With ROX, 1,000 x 50 μ l reactions
172-5857	iTaq Supermix With ROX, 2,000 x 50 μ l reactions, 50 ml bottle

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. 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.

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 by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

ABI PRISM and VIC are trademarks of Applied Biosystems. Mx3000P is a trademark of Stratagene Corporation. 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.

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