

iTaq™ DNA Polymerase

Bring your results to the forefront with iTaq DNA polymerase.

- Antibody-mediated hot-start DNA polymerase for sensitive and specific amplification
- Performance tested over a broad range of genomic and plasmid DNA targets
- Qualified for use in both standard and real-time PCR

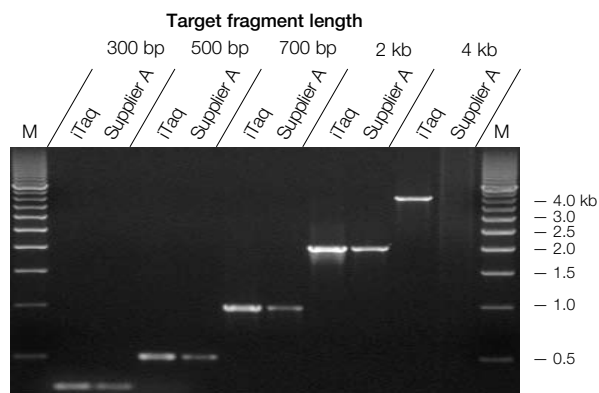
For more information, visit us on the Web at www.bio-rad.com/pcrreagents.



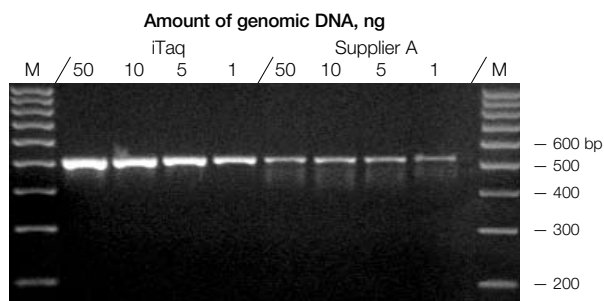
BIO-RAD

High Quality and High Specificity

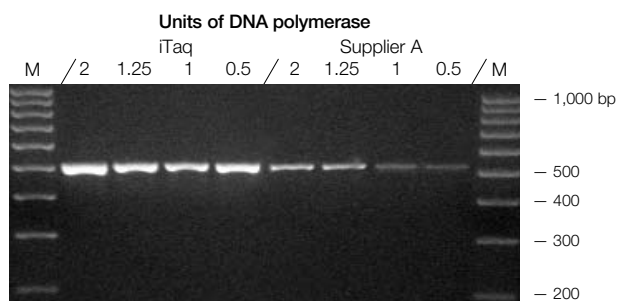
iTaq DNA polymerase is a hot-start DNA polymerase suitable for many PCR applications. The hot-start attribute is mediated through a highly specific antibody. The enzyme is activated after an initial 3 minute denaturation step at 95°C to ensure ease of use and high specificity.



iTaq DNA polymerase outperforms the gold standard chemically modified hot-start DNA polymerase over a broad range of targets. Human genomic DNA fragments were amplified with both polymerases according to manufacturers' recommendations. Lane M, EZ Load™ 500 bp molecular ruler.



iTaq DNA polymerase performs better than the competition, even at low target concentrations. A 510 bp target in human genomic DNA was amplified with iTaq DNA polymerase or supplier A's hot-start DNA polymerase according to manufacturers' recommendations. Lane M, EZ Load 100 bp molecular ruler.



The high sensitivity of iTaq DNA polymerase translates to higher yields per unit of enzyme. A 510 bp target in human genomic DNA was amplified with iTaq DNA polymerase or supplier A's hot-start DNA polymerase according to manufacturers' recommendations. Lane M, EZ Load 100 bp molecular ruler.

Ordering Information

Catalog #	Description
Reagents	
170-8870	iTaq DNA Polymerase , 5 U/μl, includes 250 U polymerase, 1.25 ml 10x PCR buffer, 1.25 ml 50 mM MgCl ₂ solution
170-8872	MgCl₂ Solution , 50 mM, 1.25 ml
170-8874	dNTP Mix , 200 μl, 10 mM each dNTP
170-8860	iQ™ Supermix , 100 x 50 μl reactions, 2x mix
170-8862	iQ Supermix , 500 x 50 μl reactions
170-8864	iQ Supermix , 1,000 x 50 μl reactions
170-8880	iQ™ SYBR® Green Supermix , 100 x 50 μl reactions, 2x mix
170-8882	iQ™ SYBR® Green Supermix , 500 x 50 μl reactions
170-8884	iQ™ SYBR® Green Supermix , 1,000 x 50 μl reactions
170-8890	iScript™ cDNA Synthesis Kit , 25 x 20 μl reactions, includes 5x iScript reaction mix, iScript reverse transcriptase, nuclease-free water
170-8891	iScript cDNA Synthesis Kit , 100 x 20 μl reactions

Thermal Cyclers

185-1096R	C1000™ Thermal Cycler With 96-Well Fast Reaction Module , includes C1000 thermal cycler chassis, 96-well fast reaction module, USB key, power cord, reagents, consumables, instructions
185-1048R	C1000 Thermal Cycler With Dual 48/48 Fast Reaction Module , includes C1000 thermal cycler chassis, dual 48/48 fast reaction module, USB key, power cord, reagents, consumables, instructions
185-1384R	C1000 Thermal Cycler With 384-Well Reaction Module , includes C1000 thermal cycler chassis, 384-well reaction module, USB key, power cord, reagents, consumables, instructions
185-2096R	S1000™ Thermal Cycler With 96-Well Fast Reaction Module , includes S1000 thermal cycler chassis, 96-well fast reaction module, power cord, reagents, consumables, instructions
185-2048R	S1000 Thermal Cycler With Dual 48/48 Fast Reaction Module , includes S1000 thermal cycler chassis, dual 48/48 fast reaction module, power cord, reagents, consumables, instructions
185-2384R	S1000 Thermal Cycler With 384-Well Reaction Module , includes S1000 thermal cycler chassis, 384-well reaction module, power cord, reagents, consumables, instructions

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.

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.

Notice regarding Bio-Rad thermal cyclers and real-time systems: Purchase of this instrument conveys a limited non-transferable immunity from suit for the purchaser's own internal research and development and for use in human in vitro diagnostics and all other applied fields under one or more of U.S. Patent Numbers 5,656,493; 5,333,675; 5,475,610 (Claims 1, 44, 158, 160-163, and 167 only); and 6,703,236 (Claims 1-7 only), or corresponding claims in their non-U.S. counterparts, owned by Applied Biosystems. No right is conveyed expressly, by implication or by estoppel under any other patent claim, such as claims to apparatus, reagents, kits, or methods such as 5' nuclease methods. 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 thermal cyclers and 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.

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 20 8732 2339 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 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