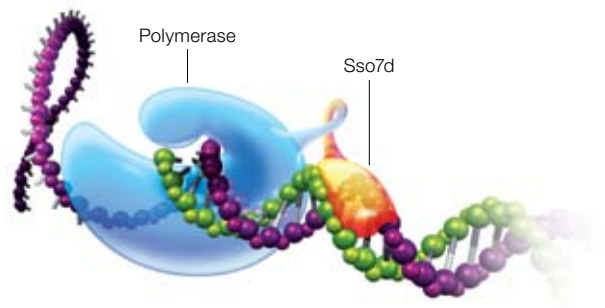


iProof™ High-Fidelity DNA Polymerase

Bio-Rad's unique iProof high-fidelity DNA polymerase delivers unsurpassed speed and fidelity, with higher yields and fewer reaction failures, for all your PCR needs.

- **Fidelity** — novel proofreading enzyme is the most accurate thermostable polymerase (52-fold more accurate than *Taq*)
- **Speed** — increased processivity dramatically reduces extension steps (15–30 sec/kb) and overall reaction times
- **Length** — large fragments (up to 37 kb) amplified in less time and with less enzyme (0.25–1.0 unit/reaction)



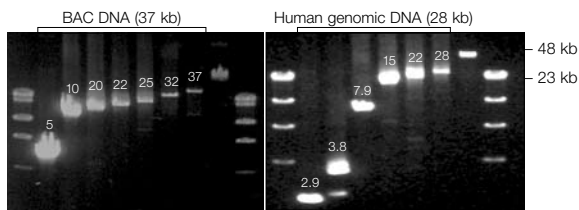
Fusing the double-stranded DNA binding protein Sso7d to iProof polymerase gives it a powerful sliding grip on the replicated DNA.

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

BIO-RAD

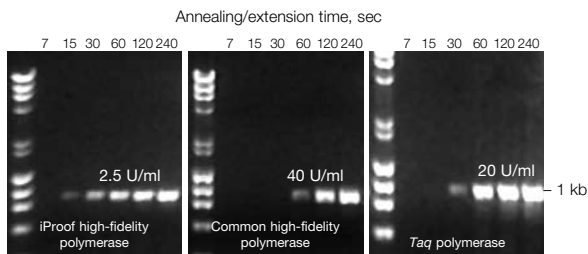
Unsurpassed Speed and Fidelity for All Your PCR Needs

iProof high-fidelity DNA polymerase consists of a unique *Pyrococcus*-like proofreading enzyme fused to a dsDNA binding protein, Sso7d. This novel technology results in a thermostable polymerase capable of amplifying long products from a variety of DNA templates, while providing the highest fidelity of any available polymerase (52-fold more accurate than *Taq*). iProof polymerase is available in three convenient formats: as a stand-alone enzyme, as an easy-to-use master mix, and in a PCR kit complete with controls.

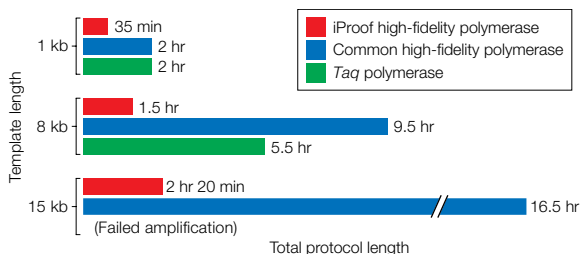


iProof DNA polymerase amplifies long templates with high yields.

Left, various fragments up to 37 kb in length were amplified from BAC DNA using a combined annealing/extension step of 10 min per cycle and 30 U/ml of iProof polymerase. Right, various sequences up to 28.8 kb were amplified directly from human genomic DNA using 30 U/ml of iProof polymerase in GC buffer with a combined annealing/extension time of 10 min per cycle.



iProof DNA polymerase delivers high yields with significantly less enzyme. A 1 kb fragment was amplified from 1.25 pM of λ template DNA in 20 cycles using a combined annealing/extension step of 7–240 sec per cycle. All enzymes were used according to manufacturer recommendations.



iProof DNA polymerase demonstrates unrivaled speed, leading to dramatically shorter overall reaction times. The reaction protocol for iProof polymerase was compared to the recommended protocols for two competing polymerases. Each protocol was designed to amplify 1, 8, and 15 kb products in 30 cycles. iProof polymerase reactions used a two-step protocol with combined annealing and extension, while the other reactions used three-step protocols with the minimum recommended extension times. Overall reaction times include temperature ramping times.

Ordering Information

Catalog #	Description
172-5300	iProof High-Fidelity DNA Polymerase, 2 U/ μ l, 20 U, includes 5x reaction buffers, MgCl ₂ solution, DMSO
172-5301	iProof High-Fidelity DNA Polymerase, 2 U/ μ l, 100 U
172-5302	iProof High-Fidelity DNA Polymerase, 2 U/ μ l, 500 U
172-5310	iProof HF Master Mix, 100 x 50 μ l reactions, includes 2x master mix, DMSO (for highest fidelity with most templates)
172-5311	iProof HF Master Mix, 500 x 50 μ l reactions
172-5320	iProof GC Master Mix, 100 x 50 μ l reactions, includes 2x master mix, DMSO (for GC-rich templates)
172-5321	iProof GC Master Mix, 500 x 50 μ l reactions
172-5330	iProof High-Fidelity PCR Kit, 2 U/ μ l, 50 U, includes 5x reaction buffers, MgCl ₂ solution, DMSO, dNTPs, λ DNA, 1.3 and 10 kb primers, DNA standard
172-5331	iProof High-Fidelity PCR Kit, 2 U/ μ l, 200 U
172-5391	5x iProof HF Buffer
172-5392	5x iProof GC Buffer
172-5393	5x iProof HPLC HF Buffer, detergent-free
172-5394	5x iProof HPLC GC Buffer, detergent-free
170-8870	iTaq™ DNA Polymerase, 5 U/ μ l, includes 250 U polymerase, 1.25 ml of 10x PCR buffer, 1.25 ml of 50 mM MgCl ₂ solution
170-8882	MgCl ₂ Solution, 50 mM
170-8874	dNTP Mix, 200 μ l premixed solution, contains 10 mM each dNTP (dATP, dCTP, dGTP, dTTP)
170-8890	iScript™ cDNA Synthesis Kit, 25 x 20 μ l reactions, includes 5x iScript reaction mix, iScript reverse transcriptase, nuclease-free water
170-8891	iScript cDNA Synthesis Kit, 100 x 20 μ l reactions
170-8896	iScript™ Select cDNA Synthesis Kit, 25 x 20 μ l reactions, includes 5x iScript Select reaction mix, iScript reverse transcriptase, oligo(dT) mix, random primer mix, gene-specific primer (GPS) enhancer solution, nuclease-free water
170-8897	iScript Select cDNA Synthesis Kit, 100 x 20 μ l reactions

Thermal Cyclers

185-1096	C1000™ Thermal Cycler With 96-Well Fast Reaction Module, includes thermal cycler chassis, 96-well fast reaction module, USB flash drive, power cord, reagent and consumable samples, instructions
185-1048	C1000 Thermal Cycler With Dual 48/48 Fast Reaction Module, includes thermal cycler chassis, dual 48/48 fast reaction module, USB flash drive, power cord, reagent and consumable samples, instructions
185-1384	C1000 Thermal Cycler With 384-Well Reaction Module, includes thermal cycler chassis, 384-well reaction module, USB flash drive, power cord, reagent and consumable samples, instructions
185-2096	S1000™ Thermal Cycler With 96-Well Fast Reaction Module, includes thermal cycler chassis, 96-well fast reaction module, power cord, reagent and consumable samples, instructions
185-2048	S1000 Thermal Cycler With Dual 48/48 Fast Reaction Module, includes thermal cycler chassis, dual 48/48 fast reaction module, power cord, reagent and consumable samples, instructions
185-2384	S1000 Thermal Cycler With 384-Well Reaction Module, includes thermal cycler chassis, 384-well reaction module, power cord, reagent and consumable samples, instructions

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.

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 applied fields other than Human In Vitro Diagnostics under one or more of U.S. Patents Nos. 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

**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