

## AccuPrime™ GC-Rich DNA Polymerase

**Cat. No:**

**12337-016**

**12337-024**

**Size:**

**200 reactions**

**1000 reactions**

**Conc: 2 U/μl**

**Store at -20°C** in a non-frost-free freezer

### Description

AccuPrime™ GC-Rich DNA Polymerase is a robust enzyme formulation designed for high-specificity, high-yield PCR amplification of difficult GC-rich templates (>65% GC content). This extremely thermostable DNA polymerase, from the archaebacterium *Pyrolobus fumarius*, retains full activity after incubation at 95°C for 4 hours and has five-fold better processivity than *Taq* DNA polymerase.

The enzyme is supplied with two separate 5X AccuPrime™ GC-Rich Buffer mixtures (A and B) containing thermostable AccuPrime™ proteins, MgSO<sub>4</sub>, and dNTPs. Thermostable AccuPrime™ proteins enhance primer-template hybridization during every cycle of PCR, greatly increasing the specificity and robustness of the reaction. Buffer A is optimized for GC-rich genomic DNA targets, while Buffer B is optimized for non-GC-rich genomic DNA, cDNA, and plasmids.

Sufficient reagents are provided for 200 or 1000 amplification reactions of 25 μl each (at 1 unit of enzyme per reaction).

<b>Component</b>	<b>Kit Size</b>	
	<b>200 Rxns</b>	<b>1000 Rxns</b>
AccuPrime™ GC-Rich DNA Polymerase	100 μl	500 μl
5X AccuPrime™ GC-Rich Buffer A	1 ml	5 ml
5X AccuPrime™ GC-Rich Buffer B	1 ml	5 ml
50-mM MgSO <sub>4</sub>	1 ml	1 ml

### Unit Definition

One unit of enzyme is the amount of enzyme required to incorporate 10 nmoles of dNTPs into acid insoluble material in 30 minutes at 74°C.

Part no. 12337.pps

MAN0001075

Rev. Date: 11 Jun 2010

### Enzyme Storage Buffer

2 U/ $\mu$ l in 50-mM Tris-HCl (pH 8.0), 100-mM KCl, 1-mM Dithiothreitol (DTT), 0.1-mM EDTA, 50% Glycerol, and 0.1% Triton<sup>®</sup> X-100

### 5X AccuPrime™ GC-Rich Buffer

Buffer A and B differ in their concentration of MgSO<sub>4</sub> and enhancers. Key components are:

300-mM Tris-HCl (pH 9.2), MgSO<sub>4</sub> at 10 mM (Buffer A) or 7.5 mM (Buffer B), 150-mM NaCl, 1-mM dGTP, 1-mM dATP, 1-mM dTTP, 1-mM dCTP, thermostable AccuPrime™ proteins, and enhancers

### Product Qualification

The Certificate of Analysis (CofA) provides detailed quality control information for each product. The CofA is available on our website at [www.invitrogen.com/cofa](http://www.invitrogen.com/cofa), and is searchable by product lot number, which is printed on each box.

### Recommendations and Guidelines

**Template:** Use 5–100 ng genomic DNA or plasmid DNA, or 10–100 ng cDNA or bacteriophage lambda DNA

**Primers:** Use  $\geq$ 50 ng each primer per 25- $\mu$ l reaction. A T<sub>m</sub> of 65–70°C is optimal for most applications. Primer design is one of the most important factors in successful PCR. We recommend using the OligoPerfect™ Designer, available at [www.invitrogen.com/oligos](http://www.invitrogen.com/oligos).

**Buffers:** In general, we recommend using Buffer A for GC-rich genomic DNA targets and Buffer B for non-GC-rich genomic DNA, cDNA, and plasmids. Also use Buffer B if you find that Buffer A is inhibitory with your genomic targets.

**Magnesium:** MgSO<sub>4</sub> is included in Buffer A at a final concentration of 2 mM and Buffer B at 1.5 mM. For some targets, more Mg<sup>2+</sup> may be required; use the 50-mM MgSO<sub>4</sub> provided in the kit to prepare a titration from 2 mM to 4 mM (final concentration) in 0.25-mM increments.

**Reaction:** Take appropriate precautions to avoid cross-contamination of DNA between reactions. Ideally, amplification reactions should be assembled in a DNA-free environment. Use of aerosol-resistant barrier tips is recommended.

## Protocol

The following protocol is recommended as a starting point. Optimal reaction conditions (incubation times and temperatures; concentrations of enzyme, primers, and template) may vary. After preparation of the samples, transfer them immediately to a preheated thermal cycler and start the amplification program.

1. Add components in the following order to each reaction vessel. Prepare a master mix for multiple reactions to enable accurate pipetting.

DNA template (see previous page)	x $\mu\text{l}$
Sense primer (10 $\mu\text{M}$ )	0.5 $\mu\text{l}$
Anti-sense primer (10 $\mu\text{M}$ )	0.5 $\mu\text{l}$
5X Buffer A or B	5 $\mu\text{l}$
AccuPrime™ GC-Rich DNA Polymerase (2 U/ $\mu\text{l}$ )*	0.5 $\mu\text{l}$
Sterile water	to 25 $\mu\text{l}$

\*Up to 2 U of enzyme (1  $\mu\text{l}$ ) may be added for difficult templates.

2. Cap/seal the reaction vessels and flick with your finger for several seconds to mix.
3. Program the thermal cycler as follows. Note that the annealing temperature will vary depending on the  $T_m$  of your primers. The optimal annealing temperature is typically 5°C below the  $T_m$  of the primers.

Step	Temp (GC-rich template)	Time	Cycle
Denaturation	95°C	3 min	1
Denaturation	95°C	30 sec	25–30
Annealing	55–65°C (5°C < $T_m$ )	30 sec	
Extension	72°C	1 min/kb	
Final Extension	72°C	10 min	1

4. Maintain the reaction at 4°C after cycling. The samples can be stored at –20°C until use. Analyze 5–10  $\mu\text{l}$  of sample by agarose gel electrophoresis.

## Purchaser Notification

### Limited Use Label License No. 1: Thermostable Polymerases

Use of this product is covered by one or more of the following US patents and corresponding patent claims outside the U.S.: 5,789,224, 5,618,711, 6,127,155. The purchase of this product includes a limited, non-transferable immunity from suit under the foregoing patent claims for using only this amount of product for the purchaser's own internal research. No right under any other patent claim, no right to perform any patented method, and no right to perform commercial services of any kind, including without limitation reporting the results of purchaser's activities for a fee or other commercial consideration, is conveyed expressly, by implication, or by estoppel. This product is for research use only. Diagnostic uses under Roche patents require a separate license from Roche. 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.

### Limited Use Label License No. 5: Invitrogen Technology

The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) not to transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. For products that are subject to multiple limited use label licenses, the terms of the most restrictive limited use label license shall control. Life Technologies Corporation will not assert a claim against the buyer of infringement of patents owned or controlled by Life Technologies Corporation which cover this product based upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Life Technologies is willing to accept return of the product with a full refund. For information about purchasing a license to use this product or the technology embedded in it for any use other than for research use please contact Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008 ; Phone (760) 603-7200 or e-mail: outlicensing@lifetech.com.

### Limited Use Label License No. 33: AccuPrime™ Reagent

This product is sold under licensing arrangements with Stratagene. The purchase price of this product includes limited, nontransferable rights under U.S. Patents Nos. 5,449,603; 5,605,824; 5,646,019; and 5,773,257 and foreign equivalents owned by Stratagene to use only this amount of the product to practice the claims in said patents solely for activities of the purchaser within the field of research. Further information on purchasing licenses under the above patents may be obtained by contacting the Director of Business Development, Stratagene, 11011 North Torrey Pines Road, La Jolla, California 92037.

### Limited Use Label License No. 82: ThermalAce™ Enzyme

Life Technologies has an exclusive license under US Patent No. 5,948,666, pending US patent applications and corresponding foreign patents and patent applications owned and licensed by Diversa Corporation to sell the ThermalAce™ enzyme to scientists for internal research purposes only, under the terms described below. Use of the ThermalAce™ enzyme by commercial entities for purposes other than research requires the user to obtain a commercial license as detailed below. The ThermalAce™ enzyme is for research use only and may not be used in clinical diagnostics without a separate license. Life Technologies Corporation ("Life Technologies") grants you a non-exclusive license to use the ThermalAce™ enzyme for research purposes only. The enzyme is being transferred to you in furtherance of, and reliance on, such license. You may not use the ThermalAce™ enzyme for commercial purposes other than research without a license for such purpose from Diversa Corporation, San Diego, California. Commercial purposes include: any use of ThermalAce™ in the manufacture of a commercial product; any sale or transfer for consideration of ThermalAce™ enzyme; any clinical diagnostic or therapeutic application employing ThermalAce™ enzyme. Access to the ThermalAce™ enzyme must be limited solely to those officers, employees and students of your institution who need access to perform the above-described internal research. You must inform each such officer, employee and student of the provisions of this license agreement and require them to agree to be bound by the provisions of this license agreement. You may not distribute the enzyme, even to those within your own institution. You may not assign, sub-license, rent, lease or otherwise transfer this license agreement or any of the rights or obligation there under, except as expressly permitted by Life Technologies and Diversa Corporation. This license agreement is effective until terminated. You may terminate it at any time by destroying all the ThermalAce™ enzyme in your control. It will also terminate automatically if you fail to comply with the terms and conditions of this license agreement. You shall, upon termination of the license agreement, destroy all ThermalAce™ enzyme in your control. For information on commercial use, please contact: Licensing Coordinator, Diversa Corporation, 4955 Directors Place, San Diego, CA 92121. Tel: (858) 526-5128, Fax: (858) 526-5628.

©2010 Life Technologies Corporation. All rights reserved.

For research use only. Not intended for any animal or human therapeutic or diagnostic use.

The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.