

AccuPrime[™] Taq DNA Polymerase System

Cat. No. 12339-016 Cat. No. 12339-024 Size: 200 Reactions Size: 1,000 Reactions

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

Description

The AccuPrime[™] *Taq* DNA Polymerase System provides qualified reagents for the amplification of nucleic acid templates by polymerase chain reaction (PCR). AccuPrime[™] *Taq* DNA polymerase contains anti-*Taq* DNA polymerase antibodies. The 10X AccuPrime[™] buffers contain thermostable AccuPrime[™] protein, Mg⁺⁺, and deoxyribonucleotide triphosphates at concentrations sufficient to allow amplification during PCR. Two individual buffer systems (10X AccuPrime[™] PCR Buffer I and II) are provided for amplification of specific types of templates.* Reagents sufficient for 200 or 1,000 amplification reactions of 25 µl each are provided.

Anti-*Taq* DNA polymerase antibodies inhibit polymerase activity, providing an automatic "hot start" (1,2) and permitting room temperature set-up. The thermostable AccuPrimeTM protein enhances specific primer-template hybridization during every cycle of PCR. Antibody/AccuPrimeTM protein-mediated amplification dramatically improves PCR specificity, improves the fidelity of *Taq* by two fold, and provides the most robust PCR for multiplex PCR and suboptimal primer sets.

| | Amount | |
|--|-------------|----------------------|
| <u>Component</u> | 200 rxn kit | <u>1,000 rxn kit</u> |
| AccuPrime [™] <i>Taq</i> DNA Polymerase | 100 µl | 500 µl |
| 10X AccuPrime [™] PCR Buffer I* | 500 µl | 2×1.25 ml |
| 10X AccuPrime [™] PCR Buffer II* | 500 µl | 2×1.25 ml |
| 50 mM Magnesium Chloride | 500 µl | 500 µl |

*10X AccuPrime[™] PCR Buffer I is designed for small genomic DNA amplicons (≤200 bp), cDNA, or plasmids. 10X AccuPrime[™] PCR Buffer II is designed for genomic DNA (200 bp–4 kb).

Storage Buffer

20 mM Tris-HCl (pH 8.0), 0.1 mM EDTA, 1 mM DTT, stabilizers, 50% (v/v) glycerol

10X AccuPrime[™] PCR Buffer I and II

200 mM Tris-HCl (pH 8.4), 500 mM KCl, 15 mM MgCl₂, 2 mM dGTP, 2 mM dATP, 2 mM dTTP, 2 mM dCTP, thermostable AccuPrime[™] protein, 10% glycerol

Product Qualification

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

Intended Use

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

Part No. 12339.pps

MAN0001076

Rev. date: 2 Jun 2010

PCR Precautions

Since PCR is a powerful technique capable of amplifying trace amounts of DNA, all appropriate precautions should be taken to avoid cross-contamination. Ideally, amplification reactions should be assembled in a DNA-free environment. Use of aerosol-resistant barrier tips is recommended. Take care to avoid contamination with the primers or template DNA used in individual reactions. PCR products should be analyzed in an area separate from the reaction assembly area.

General Protocol

The following procedure is suggested as a general guideline when using AccuPrimeTM *Taq* DNA Polymerase in any PCR amplification. Optimal reaction conditions (incubation times and temperatures; the amounts of AccuPrimeTM *Taq* DNA Polymerase, primers, MgCl₂, and template DNA) will vary and may need to be optimized. Reaction sizes may be altered to suit user preference, as shown in the tables below.

1. Add the following components to a sterile thin walled 0.25-ml or 0.5-ml PCR tube either at room temperature or on ice:

| Component | 10-µl Reaction | 25-µl Reaction | 50-µl Reaction |
|---|----------------|----------------|----------------|
| 10X AccuPrime [™] PCR Buffer I | 1 µl | 2.5 μl | 5 µl |
| Primer Mix (10 µM each) | 0.2 μl | 0.5 µl | 1 µl |
| Template DNA | 10 pg–200 ng | 10 pg–200 ng | 10 pg–200 ng |
| AccuPrime [™] <i>Taq</i> DNA Polymerase | 0.25 μl | 0.5 μl | 1 µl |
| Autoclaved distilled water | To 10 μl | To 25 μl | To 50 μl |

For Small Genomic DNA (<200 bp), Plasmids, or cDNA:

For Genomic DNA (200 bp-4 kb):

| Component | 10-µl Reaction | 25-µl Reaction | 50-µl Reaction |
|--|----------------|----------------|----------------|
| 10X AccuPrime [™] PCR Buffer II | 1 µl | 2.5 µl | 5 µl |
| Primer Mix (10 μM each) | 0.2 μl | 0.5 µl | 1 µl |
| Template DNA | 1–200 ng | 1–200 ng | 1–200 ng |
| AccuPrime™ <i>Taq</i> DNA Polymerase | 0.25 μl | 0.5 μl | 1 µl |
| Autoclaved distilled water | To 10 μl | To 25 μl | To 50 μl |

Note: Prepare a master mix for multiple reactions to minimize reagent loss and variation due to pipetting.

- 2. Mix contents of the tubes and overlay with 50 µl of mineral or silicone oil, if necessary.
- 3. Cap the tubes and centrifuge briefly to collect the contents.
- 4. Incubate tubes in a thermal cycler at 94°C for 2 min to completely denature the template and activate the enzyme.
- 5. Perform 25-35 cycles of PCR amplification as follows:

Denature: 94°C for 15-30 s

Anneal: 55°C-60°C for 15-30 s

Extend: 68°C for 1 min per kb

- 6. Maintain the reaction at 4°C after cycling. The samples can be stored at -20°C until use.
- 7. Analyze the amplification products by agarose gel electrophoresis and visualize by ethidium bromide staining. Use appropriate molecular weight standards.

Multiplex PCR Protocol

The following specialized procedure is suggested as a guideline and as a starting point when using AccuPrime[™] *Taq* DNA Polymerase in multiplex PCR amplification. Optimal reaction conditions (incubation times and temperatures; the amounts of AccuPrime[™] *Taq* DNA Polymerase, primers, MgCl₂, and template DNA) will vary and may need to be optimized. Reaction sizes may be altered to suit user preference, as shown in the tables below.

1. Add the following components to a sterile, thin-walled 0.25-ml or 0.5-ml PCR tube either at room temperature or on ice:

| Component | Amount |
|--|-------------------------|
| 10X AccuPrime [™] PCR Buffer I (for genomic DNA ≤200 bp, cDNA, or plasmids) <i>or</i> 10X AccuPrime [™] PCR Buffer II (for genomic DNA 200 bp–4 kb) | 5 µl |
| Primer mix (10 μM each) | 1 μl each (0.2 μM each) |
| Template DNA | 100-200 ng |
| AccuPrime [™] Taq DNA Polymerase* | 1-2.5 μl |
| Autoclaved, distilled water | to 50 μl |

*For primer mixes up to 5 sets, $1 \mu l$ of enzyme is sufficient.

If desired, a master mix can be prepared for multiple reactions, to minimize reagent loss and to enable accurate pipetting.

2. Continue with steps 2-7 of the General Protocol.

References

- 1. Chou, Q., Russel, M., Birch, D., Raymond, J., Bloch, W. (1992) Nucl. Acids Res. 20, 1717.
- 2. Sharkey, D.J., Scalice, E.R., Christy, K.G., Atwood, S.M., Daiss, J.L. (1994) *BioTechnology* 12, 506.
- 3. Westfall, B., Sitaraman, K., Solus, J., Hughes, J., Rashtchian, A. (1997) *Focus*[®] 19, 46.

Limited 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 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. 14: Direct Inhibition by Anti-Polymerase Antibodies

Licensed to Life Technologies Corporation, under U.S. Patent Nos. 5,338,671; 5,587,287, and foreign equivalents for use in research only.

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.

©2010 Life Technologies Corporation. All rights reserved.

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