# AccuPrime<sup>™</sup> *Pfx* SuperMix

Catalog Number 12344-040

Doc. Part No. 12344040.pps Pub. No. MAN0001080 Rev. A.0



**WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from **thermofisher.com/support**.

## **Product description**

The Invitrogen<sup>™</sup> AccuPrime<sup>™</sup> Pfx SuperMix provides qualified reagents for the high-fidelity PCR amplification of DNA templates. It includes recombinant DNA polymerase from *Thermococcus* species KOD, anti-KOD antibodies, thermostable AccuPrime<sup>™</sup> proteins, MgSO<sub>4</sub>, dNTPs, and stabilizers in a convenient and highly optimized SuperMix formulation for ease of reaction setup.

The AccuPrime™ Pfx DNA Polymerase possesses a proofreading 3′ to 5′ exonuclease activity that provides higher fidelity than Pfu DNA polymerase. This highly processive enzyme is provided in an antibody-bound form that is inactive at room temperatures. The enzyme regains activity after the initial denaturation step at 94°C in PCR cycling, providing an automatic "hot start". The hot start increases specificity, sensitivity, and yield, while allowing room-temperature assembly. The thermostable AccuPrime™ proteins enhance specific primer-template hybridization during every cycle of PCR. The high specificity, fidelity, and yield offered by AccuPrime™ Pfx SuperMix make it ideal for demanding PCR applications such as site-directed mutagenesis and PCR expression cloning.

The AccuPrime<sup>™</sup> Pfx SuperMix is suitable for targets up to 15 kb in length. It is supplied at 1.1X concentration to allow ~10% of the final reaction volume to be used for the addition of primer and template solutions.

## Contents and storage

Contents	Number of reactions	Amount	Storage
AccuPrime <sup>™</sup> <i>Pfx</i> SuperMix <sup>[1]</sup>	200	4 × 1.125 mL	-20°C in a non-frost-free freezer

<sup>[1]</sup> The AccuPrime" Pfx SuperMix contains 22 U/mL of Thermococcus species KOD thermostable polymerase complexed with anti-KOD antibodies, 66 mM Tris-SO<sub>4</sub> (pH 8.4), 30.8 mM [NH4]<sub>2</sub>SO<sub>4</sub>, 11 mM KCl, 1.1 mM MgSO<sub>4</sub>, 330 µM dNTPs, AccuPrime" proteins, and stabilizers.

**Note: Unit (U) definition:** One unit of AccuPrime<sup>™</sup> *Pfx* DNA Polymerase incorporates 10 nmol of deoxyribonucleotide into acid-insoluble material in 30 minutes at 74°C.

## Procedural guidelines

- Take appropriate precautions to avoid cross-contamination.
- For multiple reactions, prepare a master mix of AccuPrime™ Pfx SuperMix and the component(s) common to all reactions.
- Use an annealing temperature that is 5–10°C lower than the T<sub>m</sub> of the primers. If needed, gradually increase the annealing temperature by 2–3°C for higher specificity.
- If the PCR efficiency is not optimal, repeat the reaction with different primer concentrations from 100–500 nM, in 100 nM increments.

#### Perform the PCR

1. Add the following components in any order to each reaction tube:

Component	Amount for one 25-µL reaction
AccuPrime™ <i>Pfx</i> SuperMix	22.5 μL
Forward and reverse primers	200 nM final concentration of each is recommended
Template DNA solution	10 pg to 200 ng

**Note:** A standard 25- $\mu$ L PCR reaction includes a combined primer and template volume of 2.5  $\mu$ L. We have observed no decrease in product yield if the amount of primer and template solution is between 0.5  $\mu$ L and 7.5  $\mu$ L.

2. Mix the tube contents; if needed, cover with mineral or silicone oil.

Note: The oil is not needed in thermal cyclers equipped with a heated lid.

- 3. Cap the tubes, then load the tubes in a thermal cycler.
- 4. Use the following PCR program as a starting point for your template and primers:

C1	Time				
Step	Temperature	Time			
Initial denaturation	95°C	5 minutes			
35 cycles of:					
Denature	95°C	15 seconds			
Anneal	55-65°C	30 seconds			
Extend	68°C	1 minute per kb			

 Maintain the reactions at 4°C after cycling. Samples can be stored at -20°C until use.

## Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.

The information in this guide is subject to change without notice.

#### DISCLAIMER

TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

### Revision history: Pub. No. MAN0001080

Revision	Date	Description
A.0	5 May 2016	Format, style, and legal updates
_	7 June 2010	Baseline for this revision history

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

Corporate entity: Life Technologies Corporation | Carlsbad, CA 92008 USA | Toll Free in USA 1 800 955 6288

©2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.