

Pfx50™ DNA Polymerase

Cat. no. 12355-012
12355-036

Size: 100 Reactions
500 Reactions

Conc. 5 U/μl

Store at -20°C

Description

Pfx50™ DNA Polymerase is a fusion enzyme consisting of recombinant DNA polymerase from the archaean *Thermococcus zilligii* fused to an accessory protein. The highly thermostable polymerase possesses a proofreading 3'→5' exonuclease activity, while the accessory protein stabilizes primer-template complexes in PCR.

Pfx50™ DNA Polymerase offers 50 times better fidelity than *Taq* DNA polymerase, coupled with high specificity and an extremely fast elongation rate (as fast as 15 seconds per kb). In addition, the fusion enzyme has an intrinsic hot-start capability for room-temperature reaction assembly.

10X *Pfx50™* PCR buffer contains MgSO₄ at a final 1X concentration of 1.2 mM. A tube of 50-mM MgSO₄ is provided for further optimization. Reagents are provided for 100, 250 or 500 PCRs of 50 μl each.

Note:

- 10X *Pfx50™* PCR buffer contains BSA; store at -20°C.
- *Pfx50™* DNA Polymerase produces blunt-end PCR products, which can be used with Directional TOPO® Cloning and Zero Blunt® TOPO® Cloning technologies.

Component

Pfx50™ DNA Polymerase (5 U/μl)
10X *Pfx50™* PCR Mix
50-mM Magnesium Sulfate

100 Rxn Kit

100 μl
1.3 ml
1 ml

500 Rxn Kit

500 μl
2 × 1.3 ml
1 ml

Part. no. 12355.pps

Rev. date: 1 Sep 2006

***Pfx50*[™] DNA Polymerase Storage Buffer**

20 mM Tris-HCl (pH 8.0), 40 mM KCl, 0.1 mM EDTA, 1 mM DTT, stabilizers, and 50% (v/v) glycerol

Unit Definition

One unit of *Pfx50*[™] DNA Polymerase incorporates 10 nmol of deoxyribonucleotide into acid-insoluble material in 30 min at 74°C.

General Recommendations and Guidelines for PCR

General PCR parameters and troubleshooting information are documented in Innis, et al (Innis et al., 1990).

Template: *Pfx50*[™] DNA Polymerase is suitable for amplifying targets up to 4 kb from the following templates:

<u>Template</u>	<u>Amount</u>
Genomic DNA	1–200 ng
Plasmid DNA	1–100 pg
cDNA	3–5 µl from 10 ng to 1 µg starting total RNA

Amplification of longer targets (up to 7 kb) may be possible, but may require more template and longer elongation times.

Primers: Use 0.3 µM per primer as a general starting point. For larger amounts of template (e.g., 200 ng genomic DNA), increasing the concentration up to 0.5 µM per primer may improve yield.

Annealing Temperature: The annealing temperature is slightly higher than with typical PCR. **The optimal annealing temperature should be ~2°C lower than the T_m of the primers used. A range of 60–68°C is recommended.**

MgSO₄: MgSO₄ is included in the 10X *Pfx50*[™] PCR Mix at a final 1X working concentration of 1.2 mM, which is sufficient for most templates. For further optimization, add 0.1 µl to 1.0 µl of 50-mM MgSO₄.

Extension Time: As little as 15 seconds per kb may be used; 30 seconds per kb is suitable for most targets. Use up to 60 seconds per kb for maximum yield.

PCR Protocol

The following procedure is suggested as a starting point when using *Pfx50*TM DNA Polymerase in any PCR amplification.

1. Program the thermal cycler as follows (see the note on annealing temperature on page 2):

Initial denaturation: 94°C for 2 minutes

35 cycles of:

Denaturation: 94°C for 15 seconds

Annealing: 60–68°C (T_m of primers minus 2°C) for 10–30 seconds

Extension: 68°C for 30–60 seconds per kb of PCR product

Final extension: 68°C for 5 minutes

2. Add the following components to an autoclaved microcentrifuge tube at room temperature (for multiple reactions, prepare a Master Mix of common components to enable accurate pipetting):

<u>Component</u>	<u>Volume</u>	<u>Final Conc.</u>
10X <i>Pfx50</i> TM PCR Mix	5 µl	1X
10 mM dNTP Mix	1.5 µl	0.3 mM each
Primer mix (10 µM each)	1.5 µl	0.3 µM each
Template DNA (see page 2 for amounts)	≥1 µl	As required
<i>Pfx50</i> TM DNA Polymerase (5 U/µl)	1 µl	5 units
Autoclaved, distilled water	to 50 µl	

3. Cap the tube, tap gently to mix, and centrifuge briefly to collect the contents.
4. Place the tube in the thermal cycler and run the program from Step 1. After cycling, maintain the reaction at 4°C. Samples can be stored at –20°C until use.
5. Analyze products using E-Gel[®] Pre-Cast agarose gels or standard agarose gel electrophoresis. Visualize by staining with SYBR SafeTM DNA gel stain or ethidium bromide.

Quality Control

*Pfx50*TM DNA Polymerase is tested in a PCR functional assay, a double-strand-endonuclease assay, and a 5'-exonuclease assay.

References

- Innis, M. A., Gelfand, D. H., Sninsky, J. J., and White, T. S. (eds) (1990) *PCR Protocols: A Guide to Methods and Applications*, Academic Press, San Diego, CA

Additional Products

<u>Product</u>	<u>Amount</u>	<u>Catalog no.</u>
10 mM dNTP Mix, PCR Grade	100 µl	18427-013
10 mM dNTP Mix, PCR Grade	1 ml	18427-088
Invitrogen custom oligonucleotides	visit www.invitrogen.com	
E-Gel [®] 1.2% Starter Pak	6 gels plus PowerBase [™]	G6000-01
E-Gel [®] 1.2% 18-Pak	18 gels	G5018-01
TrackIt [™] 1 Kb Plus DNA Ladder	100 applications	10488-085
TrackIt [™] 100 bp DNA Ladder	100 applications	10488-058
SYBR Safe [™] DNA gel stain (10,000X)	400 µl	S33102

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. Invitrogen Corporation will not assert a claim against the buyer of infringement of patents owned or controlled by Invitrogen 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, Invitrogen is willing to accept return of the product with a full refund. For information on purchasing a license to this product for purposes other than research, contact Licensing Department, Invitrogen Corporation, 1600 Faraday Avenue, Carlsbad, California 92008. Phone (760) 603-7200. Fax (760) 602-6500. Email: outlicensing@invitrogen.com.

©2005–2006 Invitrogen Corporation. All rights reserved.

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