

Platinum[®] *Taq* DNA Polymerase High Fidelity

Cat. No. 11304-050

Size: 50 reactions

Conc: 5 U/μl

Store at -20°C (non-frost-free freezer)

Kit Contents

This sample kit includes sufficient reagents for 50 amplification reactions of 50 μl each. Ordering information for standard kit sizes is provided on page 4.

Component

Amount

Platinum[®] *Taq* DNA Polymerase High Fidelity (5 U/μl)

10 μl

10X High Fidelity PCR Buffer

1.25 ml

50-mM Magnesium Sulfate

1 ml

Description

Platinum[®] *Taq* DNA Polymerase High Fidelity is an enzyme mixture composed of recombinant *Taq* DNA polymerase, *Pyrococcus species* GB-D polymerase, and Platinum[®] *Taq* Antibody (1,2). *Pyrococcus species* GB-D polymerase possesses a proofreading ability by virtue of its 3' to 5' exonuclease activity (3). Mixture of the proofreading enzyme with *Taq* DNA polymerase increases fidelity approximately six times over that of *Taq* DNA polymerase alone and allows amplification of simple and complex DNA templates over a large range of target sizes. Targets 12–20 kb can be amplified with some optimization. Targets over 20 kb require additional optimization. The enzyme mixture is provided with an optimized buffer that improves enzyme fidelity and amplification of difficult templates.

The anti-*Taq* DNA polymerase antibody complexes with and inhibits polymerase activity at room temperature. Activity is restored after the initial denaturation step in PCR cycling at 94°C, providing an automatic “hot start” for increased specificity, sensitivity, and yield (4,5).

Platinum[®] *Taq* DNA Polymerase High Fidelity is supplied at the same 5 unit per μl concentration as Platinum[®] *Taq* DNA Polymerase. No modification to PCR reactions or protocols are necessary. Like regular *Taq*, Platinum[®] *Taq* DNA Polymerase High Fidelity has a nontemplate-dependent terminal transferase activity that adds a single deoxyadenosine (A) to the 3' ends of PCR products.

Platinum[®] *Taq* DNA Polymerase High Fidelity Storage Buffer

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

10X High Fidelity PCR Buffer

600 mM Tris-SO₄ (pH 8.9), 180 mM Ammonium Sulfate

Unit Definition

One unit of Platinum[®] *Taq* DNA Polymerase High Fidelity incorporates 10 nmol of deoxyribonucleotide into acid-preceptible material in 30 minutes at 74°C.

Quality Control

The Certificate of Analysis provides detailed quality control information for this product. Certificates of Analysis are available on our website. Go to www.invitrogen.com/support and search for the Certificate of Analysis by product lot number, which is printed on the box.

Notes

- Genomic DNA and cDNA: Use a primer concentration of 0.2 μM
- Plasmid and lambda DNA: Increase the primer concentration to 0.4 μM
- Do not perform the initial denaturation for more than 30 seconds if the target is greater than 12 kb.
- For longer genomic DNA targets (>15 kb), we recommend using 2–2.5 U of Platinum[™] *Taq* DNA Polymerase High Fidelity and increasing the extension time as specified (1 min per kb).
- We recommend an extension temperature of 68°C
- A concentration of 2 mM MgSO₄ is sufficient for most targets. For further optimization, prepare a titration from 2 mM to 4 mM in 0.25-mM increments.

Protocol

The following procedure is suggested as a guideline and starting point. The reaction size may be altered to suit user preference. Due to the “hot-start” capability of the polymerase, the reaction can be set up at room temperature.

1. Program the thermal cycler as follows:

Initial denaturation: 94°C for 30 seconds to 2 minutes (see **Notes**, page 2)

25–35 cycles of:

Denature: 94°C for 15–30 seconds

Anneal: 55°C for 15–30 seconds

Extend: 68°C for 1 minute per kb of PCR product

2. Add the following components to a DNase/RNase-free microcentrifuge tube. For multiple reactions, prepare a master mix of common components to minimize reagent loss and enable accurate pipetting.

<u>Component</u>	<u>Volume</u>	<u>Final Concentration</u>
10X High Fidelity PCR Buffer	5 µl	1X
10 mM dNTP mixture	1 µl	0.2 mM each
50 mM MgSO ₄	2 µl	2 mM
Primer mix (10 µM each)	1–2 µl	0.2–0.4 µM each (see Notes)
Template DNA	≥1 µl	(as required)
Platinum [®] <i>Taq</i> High Fidelity	0.2 µl	1.0 unit*
Autoclaved, distilled water	to 50 µl	Not applicable

*1.0 unit is sufficient for amplifying most targets. In some cases, more enzyme may be required (up to 2.5 units).

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 the amplification products by agarose gel electrophoresis. We recommend using E-Gel[®] 1.2% gels and TrackIt[™] 100 bp or 1kb Plus DNA ladders (see **Additional Products** on page 4).

References

1. Innis, M.A., Myambo, K.B., Gelfand, D.H. and Brow, M.A.D. (1988) *Proc. Natl. Acad. Sci. USA* 85, 9436.
2. Barnes, W.M. (1994) *Proc. Natl. Acad. Sci. USA* 91, 2216.
3. Tindall, K.R. and Kunkel, T.A. (1988) *Biochemistry* 27, 6008.
4. Chou, Q., Russel, M., Birch, D., Raymond, J., Bloch, W. (1992) *Nucl. Acids Res.*, 20, 1717.
5. Sharkey, D.J., Scalice, E.R., Christy, K.G., Atwood, S.M., Daiss, J.L. (1994) *BioTechnology*, 12, 506.

Additional Products

<u>Product</u>	<u>Amount</u>	<u>Catalog no.</u>
Platinum [®] <i>Taq</i> DNA Polymerase High Fidelity	100 rxns	11304-011
	500 rxns	11304-029
	5,000 rxns	11304-102
10 mM dNTP Mix, PCR Grade	100 µl	18427-013
10 mM dNTP Mix, PCR Grade	1 ml	18427-088
E-Gel [®] 1.2% Starter Pak	6 gels plus PowerBase [™]	G6000-01
E-Gel [®] 1.2% 18-Pak	18 gels	G5018-01
TrackIt [™] 100 bp DNA Ladder	100 applications	10488-058
TrackIt [™] 1kb Plus DNA Ladder	100 applications	10488-085

Limited Use Label License No. 1: Thermostable Polymerases

A license under the non-U.S. counterparts of U.S. Patents Nos. 4,683,202, 4,683,195 and 4,965,188 owned by F. Hoffmann-La Roche Ltd, for use in research and development, has an up-front fee component and a running-royalty component. The purchase price of this product includes a limited, nontransferable license under the running-royalty component to use only this amount of product to practice the Polymerase Chain Reaction (PCR) and related processes described in said patents solely for the research and development activities of the purchaser when this product is used in conjunction with a thermal cycler whose use is covered by the up-front fee component. Rights to the up-front fee component must be obtained by the end user in order to have a complete license to use this product in the PCR process. Rights under the up-front fee component may be purchased from Applied Biosystems or obtained by purchasing an Authorized Thermal Cycler. This product is also an authorized reagent and may also be used under service sublicenses from Applied Biosystems under the foregoing patents. No right to perform or offer commercial services of any kind using PCR, including without limitation reporting the results of purchaser's activities for a fee or other commercial consideration, is hereby granted by implication or estoppel. Further information on purchasing licenses to practice the PCR process 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. 117: Thermostable DNA Polymerase Blend

This product is the subject of U.S. Patent No. 5,436,149 and corresponding foreign patents licensed by Takara Corporation to Invitrogen. This product may used only for research purposes and may not be resold or used for therapeutic purposes.

Limited Use Label License No. 14: Direct Inhibition by Anti-Polymerase Antibodies

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