

## Platinum<sup>®</sup> *Taq* DNA Polymerase

**Cat. No. 10966-050**

**Size: 50 reactions**

**Conc: 5 U/μl**

**Store at -20°C in a 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.

<u>Component</u>	<u>Amount</u>
Platinum <sup>®</sup> <i>Taq</i> DNA Polymerase (5 U/μl)	10 μl
10X PCR Buffer, Minus Mg	1.25 ml
50-mM Magnesium Chloride	1 ml

### Description

Platinum<sup>®</sup> *Taq* DNA Polymerase is recombinant *Taq* DNA polymerase complexed with a proprietary antibody that blocks polymerase activity at ambient temperatures. Activity is restored after the denaturation step in PCR cycling at 94°C, thereby providing an automatic “hot start” for *Taq* DNA polymerase in PCR (1,2,3). Hot starts in PCR provide increased sensitivity, specificity, and yield, while allowing assembly of reactions at room temperature. The use of this antibody helps reduce PCR optimization requirements, reaction set-up time and effort, handling of reaction components, and contamination risk, thereby improving PCR results.

Platinum<sup>®</sup> *Taq* DNA Polymerase is supplied at the same 5 unit per μl concentration as *Taq* DNA Polymerase (Cat. No. 10342-053). No modification to PCR reactions or protocols are necessary. This enzyme formulation can also be used in larger volume cocktail mixes without difficulty.

### Storage Buffer

20 mM Tris-HCl (pH 8.0), 40 mM NaCl, 2 mM Sodium Phosphate, 0.1 mM EDTA, 1 mM DTT, stabilizers, 50% (v/v) glycerol

Part No: 100004289

Rev. date: 26 Jun 2008

## 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](http://www.invitrogen.com/support) and search for the Certificate of Analysis by product lot number, which is printed on the box.

## Unit Definition

One unit of Platinum<sup>®</sup> *Taq* DNA Polymerase incorporates 10 nmol of deoxyribonucleotide into acid-precipitable material in 30 min at 74°C.

## Guidelines and Recommendations

- Since PCR is a powerful technique capable of amplifying trace amounts of DNA, all appropriate precautions should be taken to avoid cross-contamination.
- If the PCR efficiency is not optimal, repeat the reaction with different primer concentrations from 100 to 500 nM (final concentration), in 100-nM increments.
- A concentration of 1.5 mM MgCl<sub>2</sub> is sufficient for most targets. For further optimization, prepare a titration from 1.5 mM to 3 mM in 0.25-mM increments.
- For longer genomic DNA targets, we recommend using 2–2.5 U of Platinum<sup>®</sup> *Taq* DNA Polymerase and increasing the extension time as specified (1 min per kb).

## Protocol

The procedure on the following page is suggested as a guideline and starting point when using Platinum<sup>®</sup> *Taq* DNA Polymerase in any PCR amplification. Optimal reaction conditions (incubation times and temperatures, concentration of Platinum<sup>®</sup> *Taq* DNA Polymerase, primers, MgCl<sub>2</sub>, and template DNA) vary and need to be optimized. Reaction size may be altered to suit user preferences.

**Protocol, continued**

1. Add the following components to a sterile 0.5-ml microcentrifuge tube:

<u>Components</u>	<u>Volume</u>	<u>Final Concentration</u>
10X PCR Buffer, Minus Mg	5 $\mu$ l	1X
10 mM dNTP mixture	1 $\mu$ l	0.2 mM each
50 mM MgCl <sub>2</sub>	1.5 $\mu$ l	1.5 mM
Primer mix (10 $\mu$ M each)	1 $\mu$ l	0.2 $\mu$ M each
Template DNA	$\geq$ 1 $\mu$ l	(as required)
Platinum <sup>®</sup> <i>Taq</i> DNA Polymerase	0.2 $\mu$ l	1.0 unit*
Autoclaved, distilled water	to 50 $\mu$ l	Not applicable

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

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

2. Mix contents of the tubes and overlay with 50  $\mu$ l 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 30 s to 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 30 s
Anneal	55°C for 30 s
Extend	72°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 products by agarose gel electrophoresis and visualize by ethidium bromide staining. Use appropriate molecular weight standards.

## Additional Products

The following products are available on our website at [www.invitrogen.com](http://www.invitrogen.com) or by contacting technical support.

	<u>Cat. No.</u>	<u>Amount</u>
Platinum® <i>Taq</i> DNA Polymerase	10966-018	100 Reactions
	10966-026	250 Reactions
	10966-034	500 Reactions
	10966-083	5,000 Reactions

## References

1. Chou, Q., *et al.* (1992) *Nucl. Acids Res.* 20, 1717.
2. Sharkey, D.J., *et al.* (1994) *BioTechnology* 12, 506.
3. Westfall, B.A., *et al.* (1997) *Focus*® 19.3, 46.

### 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 US: 5,079,352, 5,789,224, 5,618,711, 6,127,155 and claims outside the US corresponding to US Patent No. 4,889,818. 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 (such as the patented 5' Nuclease Process claims in US Patents Nos. 5,210,015 and 5,487,972), 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. 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.

©2003–2008 Invitrogen Corporation. All rights reserved.

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

## Headquarters

5791 Van Allen Way | Carlsbad, CA 92008 USA  
Phone +1 760 603 7200 | Toll Free in USA 800 955 6288

## For support visit

[www.lifetechnologies.com/support](http://www.lifetechnologies.com/support) or email [techsupport@lifetech.com](mailto:techsupport@lifetech.com)

[www.lifetechnologies.com](http://www.lifetechnologies.com)

