



Product Manual

AMPIGENE[®] HS Taq DNA Polymerase

Catalog #: ENZ-PRT101

ENZ-PRT101-0500 for 500 units



Product Manual

USE FOR RESEARCH PURPOSES ONLY

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NON-INFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

TRADEMARKS AND PATENTS

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending. Enzo is a trademark of Enzo Life Sciences, Inc.

FOR RESEARCH USE ONLY.

**THE PRODUCT MAY BE USED ONLY FOR *IN VITRO*
RESEARCH PURPOSES.**



Please read entire booklet before proceeding with the assay.



Carefully note the handling and storage conditions.



Please contact Enzo Life Sciences Technical Support if necessary.

TABLE OF CONTENTS

| | |
|--------------------------------|---|
| Description | 2 |
| Shipping and Storage | 2 |
| Important Considerations | 3 |
| Reaction Setup | 4 |
| Contact Information | 6 |

DESCRIPTION

AmipiGene™ HS Taq DNA Polymerase uses the latest developments in polymerase technology and buffer chemistry to enhance PCR speed, yield and specificity. The enzyme uses advanced hot-start technology for superior sensitivity. The enzyme and buffer system allow for superior PCR performance on complex templates such as mammalian genomic DNA. Due to enhanced efficiency and specificity the enzyme is perfectly suited to difficult PCR.

AMPIGENE[®] HS Taq is a robust enzyme for all your everyday PCR applications including genotyping, multiplex PCR, screening, library construction, colony PCR and PCR direct from blood and urine. AMPIGENE[®] HS Taq DNA Polymerase can perform consistently well on a broad range of templates (including both GC and AT rich).

AMPIGENE[®] HS Taq DNA Polymerase has an error rate of approximately 1 error per 2.0×10^5 nucleotides incorporated. PCR products generated with AMPIGENE[®] HS Taq DNA Polymerase are A-tailed and may be cloned into TA cloning vectors.

| Component | 500 units |
|---|-----------|
| AMPIGENE [®] HS Taq DNA polymerase (5u/μl) | 2 x 50μl |
| 5x AMPIGENE [®] reaction buffer | 4 x 1ml |

High-throughput screening has resulted in a buffer system that allows efficient amplification from GC-rich and AT-rich templates, under fast and standard cycling conditions.

AMPIGENE[®] HS Taq Mix is particularly resistant to PCR inhibitors. The mix is suitable for direct PCR from unprocessed samples including bacterial culture, bacterial colonies, blood and urine.



Protect from prolonged exposure to light.

SHIPPING AND STORAGE

On arrival the kit should be stored at -20°C. Avoid prolonged exposure to light. If stored correctly the kit will retain full activity for 12 months. The kit can be stored at 4°C for 1 month. The kit can go through 30 freeze/thaw cycles with no loss of activity.

IMPORTANT CONSIDERATIONS

5x AMPIGENE[®] Reaction buffer: The 5x reaction buffer contains 15mM MgCl₂, 5mM dNTPs, enhancers and stabilizers. It is not recommended to add further PCR enhancers or MgCl₂ to the reaction. The buffer composition has been optimized to maximize PCR success rates.

Template: For eukaryotic DNA use between 5ng and 500ng per reaction, for cDNA use below 100ng per reaction.

Primers: Primers should have a predicted melting temperature of around 60°C, using default Primer 3 settings. The final primer concentration in the reaction should be between 0.2μM and 0.6μM.

Annealing: We recommend performing a temperature gradient to experimentally determine the optimal annealing temperature. Alternatively, we recommend a 55°C annealing temperature then increase in 2°C increments if non-specific products are present.

Extension: Optimal extension is achieved at 72°C. The optimal extension time is dependent on amplicon length and complexity of template. 15 seconds per kilobase (kb) is recommended for amplification from eukaryotic DNA for amplicons between 1kb and 6kb. For shorter amplicons a 1 second extension is sufficient.

Multiplex PCR: When first performing multiplex PCR it is recommended to run an annealing temperature gradient from 55°C to 65°C. The annealing temperature that results in the best specificity should be used in subsequent experiments. Fast cycling conditions should not be used for multiplex PCR. Initially, we recommend a 90 second extension time. This time may be further extended to increase yield.

Colony PCR: From bacterial colonies use a sterile tip to pick a colony and resuspend into a 50μl reaction as described below. From liquid culture add 5μl of overnight culture to the final mix. Increase initial denaturation time to 10 minutes.

Direct blood/urine PCR: Add 2μl mammalian blood or urine to a 50μl reaction as described below.

REACTION SETUP

1. Prepare a master mix based on the following table:

| Reagent | 50 μ l reaction | Final concentration | Notes |
|---|--------------------------------|---------------------|---------------------------------------|
| 5x AMPIGENE [®] Reaction buffer | 10.0 μ l | 1x | |
| Forward primer (10 μ M) | 2.0 μ l | 400nM | See above for optimal primer design |
| Reverse primer (10 μ M) | 2.0 μ l | 400nM | |
| Template DNA | <100ng cDNA, <500ng genomic | variable | See above for template considerations |
| AMPIGENE [®] HS Taq DNA polymerase (5u/ μ l) | 0.25 μ l - 1.0 μ l | | |
| PCR grade dH ₂ O | Up to 50 μ l final volume | | |

2. Cycle using conditions based on the following table:

| Cycles | Temperature | Time | Notes |
|--------|--------------|--------------|---|
| 1 | 95°C | 1-2min | Initial denaturation and enzyme activation. For colony PCR increase denaturation time to 10 minutes |
| 40 | 95°C | 15 seconds | Denaturation |
| | 55°C to 65°C | 15 seconds | Anneal |
| | 72°C | 1-90 seconds | Extension (15 seconds per kb). For multiplex PCR use 90 seconds. |



Product Manual

NOTES



Product Manual

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Toll-Free: 1.800.942.0430
Phone: 631.694.7070
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE/ASIA

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen
Switzerland
Phone: +41/0 61 926 89 89
Fax: +41/0 61 926 89 79
info-ch@enzolifesciences.com

For local distributors and detailed product information visit us online:
www.enzolifesciences.com