



## **AMPIGENE<sup>®</sup> HS Taq Mix**

Catalog #: ENZ-NUC101

ENZ-NUC101-0040 for 40 rxns

ENZ-NUC101-0200 for 200 rxns

ENZ-NUC101-1000 for 1000 rxns



## **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

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

AMPIGENE<sup>®</sup> HS Taq Mix has an error rate of approximately 1 error per  $2.0 \times 10^5$  nucleotides incorporated. PCR products generated with AMPIGENE<sup>®</sup> HS Taq DNA Polymerase are A-tailed and may be cloned into TA cloning vectors.

Component	40 reactions	200 reactions	1000 reactions
2x AMPIGENE <sup>®</sup> HS Taq Mix	1 x 1ml	5 x 1ml	25x 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<sup>®</sup> 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.

## SHIPPING AND STORAGE

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



Protect from prolonged exposure to light.

## IMPORTANT CONSIDERATIONS

**2x AMPIGENE<sup>®</sup> HS Taq Mix:** The 2x mix contains AMPIGENE<sup>®</sup> HS Taq DNA Polymerase, 6mM MgCl<sub>2</sub>, 2mM dNTPs, enhancers and stabilizers. It is not recommended to add further PCR enhancers or MgCl<sub>2</sub> 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. 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	50µl reaction	Final concentration	Notes
2x AMPIGENE <sup>®</sup> HS Taq Mix	25.0µl	1x	
Forward primer (10µM)	2.0µl	400nM	See above for optimal primer design
Reverse primer (10µM)	2.0µl	400nM	
Template DNA	<100ng cDNA, <500ng genomic	variable	See above for template considerations
PCR grade dH <sub>2</sub> O	Up to 50µl final volume		

2. Cycle using conditions based on the following table:

Cycles	Temperature	Time	Notes
1	95°C	1min to 2min	Initial denaturation and enzyme activation. For colony PCR increase denaturation time to 10 minutes
40	95°C 55°C to 65°C 72°C	15 seconds 15 seconds 1-90 seconds	Denaturation Anneal 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](http://www.enzolifesciences.com)