



AMPIGENE[®] 1-Step RT-PCR Kit

Catalog #: ENZ-KIT105

ENZ-KIT105-0050 for 50 rxns

ENZ-KIT105-0200 for 200 rxns

ENZ-KIT105-1000 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

Enzo Life Sciences' AMPIGENE[®] 1-Step RT-PCR Kit uses the latest developments in reverse transcriptase technology and buffer chemistry for efficient cDNA synthesis and PCR in a single tube.

Our modified MMLV reverse transcriptase (RTase) is both thermostable and extremely active. The enzyme is blended with RNase inhibitor preventing degradation of RNA by contaminating RNase. The RTase is not inhibited by ribosomal and transfer RNAs, total RNA is an ideal substrate.

AMPIGENE[®] 1-Step RT-PCR Kit uses proprietary small molecular inhibitor technology that prevents formation of primer-dimers to improve reaction sensitivity and specificity.

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.

Component	50 reactions	200 reactions	1000 reactions
2x 1-Step Mix	1 x 1.25ml	4 x 1.25ml	20 x 1.25ml
20x RTase with RNase inhibitor	1 x 125µl	4 x 125µl	20 x 125µl

SHIPPING AND STORAGE



Protect from prolonged exposure to light.

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

2x 1-Step Mix: The 2x mix contains HS Taq DNA Polymerase, 6mM MgCl₂, 2mM 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.

20x RTase: The 20x RTase also contains RNase inhibitor. It is essential to use the correct volume per reaction. Using the incorrect volume will result in loss of sensitivity.

Template: Use 1pg to 1µg total RNA per reaction, use minimum 0.01pg mRNA 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.

Reverse Transcription: We recommend incubating with a temperature of 45°C for 10 minutes for the majority of applications. Where regions of interest contain high secondary structure incubation temperatures up to 55°C may be used. For amplicons above 1kb the incubation time should be increased to 20 minutes.

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 3kb.

REACTION SETUP

1. Before starting, briefly vortex 2x 1-Step Mix
2. Prepare a master mix based on following table; we recommend also setting up a no-RTase control:

Reagent	50µl reaction	Final concentration	Notes
2x 1-Step Mix	25µl	1x	
Forward primer (10µM)	2.0µl	400nM	See above for optimal primer design
Reverse primer (10µM)	2.0µl	400nM	
20x RTase	2.5µl	1x	Correct volume is critical, do not reduce
Template RNA	1pg to 1µg total RNA >0.01pg mRNA	variable	
PCR grade dH ₂ O	Up to 50µl final volume		

3. Program the instrument using following conditions:

Cycles	Temperature	Time	Notes
1	45°C to 55°C	10min	Reverse transcription, 45°C is recommended for most applications, 55°C should be used only when amplicon contains regions of high secondary structure
1	95°C	2min	Polymerase activation, 2 minutes
40	95°C	10 seconds	Denaturation
	60°C to 65°C	10 seconds	Anneal
	72°C	30-60 seconds	15 seconds per kb



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