

## **AMPIXTRACT<sup>®</sup> Paraffin Tissue Section DNA Isolation Kit**

Catalog #: *ENZ-GEN501*

*ENZ-GEN501-0050 – for 50 samples*

*ENZ-GEN501-0100 – for 100 samples*

For the latest product information, including support documentation, visit  
us online:

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

Technical Support (US): 800-942-0430  
Technical Support (EU): +41 61 926 8989

## **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 use only. Not for use in diagnostic procedures. 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.  
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**

## TABLE OF CONTENTS



Carefully note the handling and storage conditions of each kit component.

INTRODUCTION .....	2
MATERIALS SUPPLIED .....	2
ADDITIONAL MATERIALS NEEDED .....	3
STORAGE & STABILITY .....	3
SAFETY WARNINGS & PRECAUTIONS .....	4
PRINCIPLE OF THE ASSAY .....	4
PROCEDURE .....	5
CONTACT INFORMATION .....	6

## INTRODUCTION

Retrospective studies with DNA on archival tissue samples would provide significant information for disease-related molecular processes. However, isolating high-quality genomic DNA from formalin-fixed, paraffin-embedded tissue can be difficult because only minimal amounts of intact DNA may be present in the sample. The AMPIXTRACT® Paraffin Tissue Section DNA Isolation Kit uses a unique procedure and composition to efficiently isolate DNA from paraffin archives. The kit has the following features:

- The fastest procedure currently available, which can be finished within 2 hours, depending on sample types, with consistent isolation conditions.
- High efficiency of DNA isolation from tissue sections containing tiny amounts of DNA (as low as 1 ng).
- No requirement for pre-deparaffinization.
- Use of non-toxic reagents and no phenol chloroform.

## MATERIALS SUPPLIED

Components	50 samples (ENZ-GEN501-0050)	100 samples (ENZ-GEN501-0100)
PS1 (DNA Digestion Solution)	0.5 ml	1 ml
PS2 (DNA Digestion Powder)	1 vial	1 vial
PS3 (DNA Isolation Buffer)	10 ml	20 ml
PS4 (DNA Binding Buffer)	12 ml	22 ml
PS5 (DNA Elution Solution)	1 ml	2 ml
F-Spin Column	50	100
F-Collection Tube	50	100
Product Manual	1	1

## ADDITIONAL MATERIALS NEEDED

The following solutions are required but not provided:

- 90% Ethanol
- 70% Ethanol



Storage temp

## STORAGE & STABILITY

The AMPIXTRACT® Paraffin Tissue Section DNA Isolation Kit can be stored at room temperature (20-22°C) for 6 months from shipping date, with the exception of PS2. Upon receipt, PS2 should be stored at -20°C or stored at 4°C as soon as it is dissolved in PS1 (stable for up to 6 months).



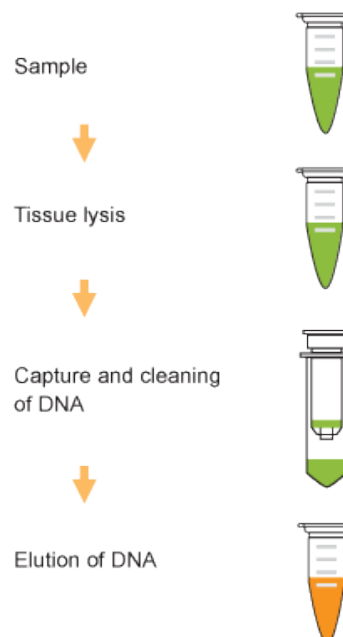
Important/ Warning

## SAFETY WARNINGS & PRECAUTIONS

1. Wear appropriate personnel protective apparel. Avoid contact with clothes and exposed skin. In case of accidental skin exposure, flush with water immediately. Consult a physician if required.
2. Use a safety pipetting device for all pipetting. Never pipet by mouth.
3. Interpretation of the results is the sole responsibility of the user.

## PRINCIPLE OF THE ASSAY

The AMPIXTRACT<sup>®</sup> Paraffin Tissue Section DNA Isolation Kit simply applies our proprietary DNA isolation buffer to the samples. After treatment with DNA digestion buffer, the DNA is easily recovered in 8-20  $\mu$ l by our specially designed Fast-Spin Column. DNA is ready for down-stream application.



Schematic Procedure for Using the AMPIXTRACT<sup>®</sup> Paraffin Tissue Section DNA Isolation Kit

## PROCEDURE

Note: Always cap spin columns before placing them in the microcentrifuge.

**Before starting, prepare the following required solutions (not included):**

**90% Ethanol and 70% Ethanol**

1. Add 0.5 ml (for P-1009-1) or 1 ml (for P-1009-2) of **PS1** to **PS2**. Vortex until **PS1/PS2 solution** is clear.
2. Remove the tissue area you need (1-40 mm<sup>2</sup> of 10 µm thick tissue) from the slide with a scalpel. Transfer it to a 1.5 ml vial and add 100 µl of **PS3** (for tissue area > 20 mm<sup>2</sup>, add 200 µl of **PS3**).
3. Incubate the sample at 95°C for 10-15 minutes. Vortex for 5-10 seconds and leave it at room temperature for 1 minute.
4. Add 5 µl of **PS1/PS2 solution** to each 100 µl of the sample solution. Mix and incubate this mixed solution at 65°C for 60-90 minutes or until paraffin tissue is completely lysed (it is usually less than 2 hours). Vortex the sample for 5-10 seconds every 30 minutes. If tissue is not properly in this mixed solution after vortexing, spin it down into the solution.
5. Vortex the mixture for 5 seconds after incubation. Add 200 µl of **PS4** to the mixture and centrifuge at 12,000 rpm for 1 minute. The paraffin layer should now be on the top and the clear solution on the bottom. Place a spin column into a 2 ml collection tube. Carefully penetrate through the paraffin layer and transfer the clear solution to the column. Spin for 30 seconds at 12,000 rpm. Discard the flowthrough. Replace the column to the collection tube (Note: maximum volume of the column is 600 µl.)
6. Add 200 µl of 70% ethanol to the column and spin at 12,000 rpm for 20 seconds. Add 200 µl of 90% ethanol to the column and spin at 12,000 rpm for 20 seconds. Discard the flowthrough and replace the column to the collection tube.
7. Add additional 200 µl of 90% ethanol to the column and spin at 12,000 rpm for 40 seconds.
8. Place the column in a new 1.5 ml vial. Add 8-18 µl of **PS5** directly to the column filter, and centrifuge at 12,000 rpm for 20 seconds to elute DNA.

DNA is now ready for use or storage at -20°C.



# 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**

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)