

## **AMPIXTRACT™ General Tissue Section DNA Isolation Kit**

Catalog #: ENZ-GEN500

*ENZ-GEN500-0050 - for 50 samples*

*ENZ-GEN500-0100 – 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



Please read entire  
booklet before  
proceeding with the  
assay.

INTRODUCTION .....	2
BACKGROUND .....	2
MATERIALS SUPPLIED.....	3
ADDITIONAL MATERIALS REQUIRED .....	4
STORAGE & STABILITY .....	4
SAFETY WARNINGS & PRECAUTIONS.....	5
PRINCIPLE OF THE ASSAY.....	5
PROCEDURE.....	6
CONTACT INFORMATION .....	10

## INTRODUCTION

The AMPIXTRACT™ kits are very suitable for isolating tiny amounts of DNA from microdissection samples, fresh tissue sections, formalin-fixed and paraffin-embedded tissues, plasma, serum, body fluids, etc. The quality of extracted DNA from formalin-fixed and paraffin-embedded tissues may be affected by the quality of the embedded tissue.

The AMPIXTRACT™ kits allow isolation of DNA size from 50 bp to 20 kb; DNA quantity from 1 ng to 2 µg, optimal at between 10 ng and 1 µg.

## BACKGROUND

The *AMPIXTRACT™* General Tissue Section DNA Isolation Kit is designed for isolating DNA from tissue sections. The kit uses a unique procedure and composition to efficiently isolate DNA in any targeted microscopic tissue area on a slide. The kit has the following features:

- The fastest procedure 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).
- Use of non-toxic reagents and no phenol chloroform.

## MATERIALS SUPPLIED

Components	50 samples (ENZ-GEN500-0050)	100 samples (ENZ-GEN500-0100)
<b>S1</b> (DNA Digestion Solution)	0.3 ml	0.6 ml
<b>S2</b> (DNA Digestion Powder)	1 vial	1 vial
<b>S3</b> (DNA Isolation Buffer)	6 ml	11 ml
<b>S4</b> (DNA Binding Buffer)	12 ml	22 ml
F-Spin Column	50	100
F-Collection Tube	50	100
Product Manual	1	1

## **ADDITIONAL MATERIALS REQUIRED**

The following solutions are required but not provided

- 90% ethanol
- 70% ethanol



Storage temp

## **STORAGE & STABILITY**

Upon receipt: (1) S2 should be stored at  $-20^{\circ}\text{C}$ , or stored at  $4^{\circ}\text{C}$  as soon as it is dissolved in S1 (up to 6 months); (2) Store all other components at room temperature. The kit can be stable for up to 6 months from the shipment date when stored properly.



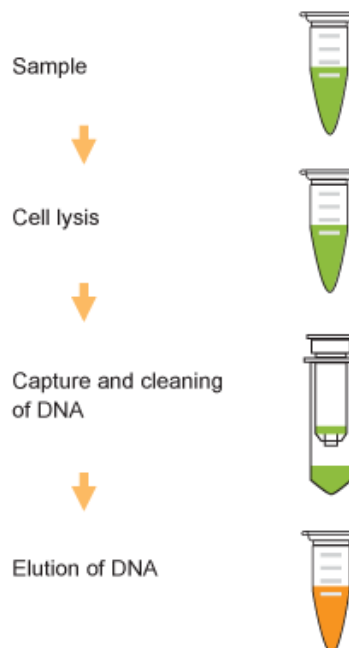
Important/ Warning

## 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™ General Tissue Section DNA Isolation Kit simply applies our proprietary DNA isolation buffer to a selected microscopic tissue area. The area is removed and transferred into a tube. 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™ General 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 0.3 ml of **S1** to **S2** in order to create the **S1/S2 solution**. Vortex until solution is clear. Spin the solution down to the bottom.
2. Treat the tissue with the DNA Isolation Buffer:

**For microdissection samples**, directly collect the sample into a vial containing 100  $\mu$ l of **S3**, followed by adding 5  $\mu$ l of **S1/S2 solution**. Mix well and incubate at 65°C for 60-90 minutes.

**For tissues from fresh sections**, add 0.5  $\mu$ l of **S3** to 1 mm<sup>2</sup> (about 500-1000 cells) of tissue area and immediately remove the tissue area you need (1- 20 mm<sup>2</sup>) from the slide with a scalpel. Transfer it to a 1.5 ml vial containing 100  $\mu$ l of **S3**, followed by adding 5  $\mu$ l of **S1/S2 solution**. Mix well and incubate this mixed solution at 65°C for 60-90 minutes or until the tissue is completely lysed (usually it is less than 2 hours). Vortex the sample for 5 seconds every 30 minutes. If tissue is not properly in this mixed solution after vortexing, spin it down into the solution.

**For paraffin samples**, remove the paraffin first with deparaffin reagents according to the manufacturer's instructions or according to the following procedures:

- 1) Drop the slide into 100% of *xylene* at room temperature for 10 min. Repeat once with new *xylene*.
- 2) Drop the slide in 100% of *ethanol*, 95% and 70% for 5 minutes each. Air dry the slide. Add 0.5  $\mu$ l of **S3** to 1 mm<sup>2</sup> of tissue area and immediately remove the tissue area you need (1-20 mm<sup>2</sup>) from the slide. Transfer it to a vial containing 100  $\mu$ l of **S3**, followed by adding 5  $\mu$ l of **S1/S2 solution**. Mix well and incubate this mixed solution at 65°C for 60-90 minutes, or until tissue is completely lysed (it is usually less than 2 hours). Vortex the sample for 5 seconds, every 30 minutes. If tissue is not properly in this mixed solution after vortexing, spin it down into the solution.



3. Place a spin column into a 2 ml collection tube. Vortex the mixture for 5 seconds after incubation. Add 200  $\mu$ l of **S4** to the mixture and transfer it to the column. Centrifuge at 12,000 rpm for 30 seconds. Discard the flowthrough. Replace the column to the collection tube (*Note: maximum volume of the column is 600  $\mu$ l.*)
4. Add 200  $\mu$ l of 70% ethanol to the column and centrifuge at 12,000 rpm for 20 seconds. Add 200  $\mu$ l of 90% ethanol to the column and centrifuge at 12,000 rpm for 20 seconds. Discard the flowthrough and replace the column to the collection tube.
5. Add an additional 200  $\mu$ l of 90% ethanol to the column and centrifuge at 12,000 rpm for 40 seconds.
6. Place the column in a new 1.5 ml vial. Add 8-18  $\mu$ l of **S5** directly to the column filter and centrifuge at 12,000 rpm for 20 seconds to elute DNA.



# Product Manual

## NOTES

---



# 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**

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)**