



## **EpiXtract™ Nuclear Protein Isolation Kit II (Nucleic acid-free)**

Catalog #: ENZ-45015

100 Assays



## **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.  
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**

## TABLE OF CONTENTS



Please read  
entire booklet  
before  
proceeding with  
the assay.

<b>Introduction.....</b>	<b>2</b>
<b>Safety Warnings and Precautions.....</b>	<b>2</b>
<b>Reagents Provided and Storage .....</b>	<b>3</b>
<b>Additional Materials Required.....</b>	<b>3</b>
<b>Methods and Procedure.....</b>	<b>4</b>
Monolayer and adherent cells preparation .....	4
Suspension cells preparation.....	4
Tissue sample preparation .....	5
Nuclear extraction.....	5
<b>Contact Information.....</b>	<b>6</b>

## INTRODUCTION

The EpiXtract™ Nuclear Protein Isolation Kit II (nucleic acid-free) uses a selective and easy to use method for extracting nuclear proteins that can be used for a variety of downstream applications. Applications include Western blot analysis, protein-DNA binding assays, nuclear enzyme assays, and any other procedures that require optimization or the use isolated nucleic acid-free nuclear proteins. The EpiXtract™ Nuclear Protein Isolation Kit II (nucleic acid-free) is compatible with mammalian adherent or suspension cells and tissue samples.

Using the EpiXtract™ Nuclear Protein Isolation Kit II, the extraction procedure can be completed within 1 hour. Under optimal extraction conditions a total yield can be up to 100µg.

## SAFETY WARNINGS & PRECAUTIONS

### **FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.**

Observe good laboratory practices. Gloves, lab coat, and protective eyewear should always be worn. Never pipet by mouth. Do not eat, drink or smoke in the laboratory areas. All blood components and biological materials should be treated as potentially hazardous and handled as such. They should be disposed of in accordance with established safety procedures.

## REAGENTS PROVIDED AND STORAGE

The EpiXtract™ Nuclear Protein Isolation Kit II is shipped on ice.

Upon receipt: (1) Store **Extraction Pre-Cleaner** and **Extraction Cleaner** store at  $-20^{\circ}\text{C}$  in aliquots; and (2) Store all other components at  $4^{\circ}\text{C}$ .

Reagent	Volume	Storage Temperature
10x Pre-Extraction Buffer	10 mL	$4^{\circ}\text{C}$
Extraction Buffer	10 mL	$4^{\circ}\text{C}$
Extraction Pre-Cleaner	1 mL	$-20^{\circ}\text{C}$
Extraction Cleaner	100 $\mu\text{L}$	$-20^{\circ}\text{C}$
1000x DTT Solution	100 $\mu\text{L}$	$4^{\circ}\text{C}$
1000x Protease Inhibitor Cocktail	100 $\mu\text{L}$	$4^{\circ}\text{C}$

## ADDITIONAL MATERIALS REQUIRED

- Adjustable pipette
- Aerosol resistant pipette tips
- Distilled water
- Conical tubes
- microcentrifuge vials
- tissue homogenizer
- Vortex
- Microcentrifuge
- Hemacytometer
- Freshly made PBS solution
- UV Spectrometer
- Ice

## METHODS AND PROCEDURES

### Monolayer or adherent cells preparation

1. Grow cells to 70-80% confluency on a culture plate or flask. Cell density should not exceed  $2-5 \times 10^6$  cells for a 100 mm plate. Remove the growth medium and wash cells with PBS twice and then remove PBS.
2. Add 1mL of fresh PBS per 20 cm<sup>2</sup> area (e.g., add 3 mL of PBS to a 100 mm plate) and scrape cells into a 15 mL conical tube. Alternatively, cells can be detached with trypsin/EDTA, collected into a 15 mL conical tube and counted with a hemacytometer.
3. Centrifuge the cells for 5 min at 1000 rpm and discard the supernatant.
4. Dilute the **10x Pre-Extraction Buffer** with distilled water at a 1/10 ratio (e.g., 1 mL of **10x Pre-Extraction Buffer** and 9 mL of distilled water) Add **1000x DTT Solution** and **1000x Protease Inhibitor Cocktail** to ice cold diluted (**1x**) **Pre-Extraction Buffer** at a 1/1000 ratio. Resuspend cell pellet in 100  $\mu$ L of this mix per  $10^6$  cells. Transfer to a microcentrifuge vial.
5. Incubate on ice for 10 minutes. Vortex vigorously for 10 seconds and centrifuge the preparation for 1 minute at 12,000rpm.
6. Carefully remove the cytoplasmic extract from the nuclear pellet.
7. Proceed to **Section D, Nuclear Extraction**.

### Suspension cells preparation

1. Grow cells to a  $2 \times 10^6$ /ml. Collect the cells in a 15 mL conical tube.
2. Centrifuge the cells for 5 minutes at 1000rpm and discard the supernatant. Wash cells with PBS once by centrifugation for 5 minutes at 1000rpm. Discard the supernatant.
3. Dilute the **10x Pre-Extraction Buffer** with distilled water at a 1/10 ratio (e.g., 1 mL of **10x Pre-Extraction Buffer** and 9mL of distilled water). Add **1000x DTT Solution** and **1000x Protease Inhibitor Cocktail** to ice cold diluted (**1x**) **Pre-Extraction Buffer** at a 1/1000 ratio. Resuspend cell pellet in 100 $\mu$ L of this mix per  $10^6$  cells. Transfer to a microcentrifuge vial.

4. Incubate on ice for 10 minutes. Vortex vigorously for 10 seconds and centrifuge the preparation for 1 minute at 12,000rpm.
5. Carefully remove the cytoplasmic extract from the nuclear pellet.
6. Proceed to **Section D, Nuclear Extraction**.

#### Tissue sample preparation

1. Weigh tissue and cut it into small pieces. Place the cut pieces in a clean homogenizer
2. Dilute the **10x Pre-Extraction Buffer** with distilled water at a 1/10 ratio (e.g., 1 mL of **10x Pre-Extraction Buffer** and 9mL of distilled water). Add 5 mL of diluted (1x) **Pre-Extraction Buffer** containing 5  $\mu$ L of **DTT Solution** per gram of tissue, and homogenize tissue pieces (50-60 strokes).
3. Incubate on ice for 15 minutes and centrifuge for 10 minutes at 12,000rpm at 4°C. Remove the supernatant.
4. Proceed to **Section D, Nuclear Extraction**.

#### Nuclear extraction

1. Add **DTT Solution** and **Protease Inhibitor Cocktail** to **Extraction Buffer** at a 1/1000 ratio, followed by adding **Extraction Pre-Cleaner** to **Extraction Buffer** at a 1/10 ratio. Add 2 volumes of **Extraction Buffer** mix to the nuclear pellet (about 10 $\mu$ L **Extraction Buffer** mix per 10<sup>6</sup> cells or per 2 mg of tissue).
2. Incubate the extract on ice for 15 minutes. Vortex (for 5 seconds) every 3 minutes. The extract (especially tissue extract) can be further sonicated for 3 x 10 seconds to increase nuclear protein extraction.
3. Centrifuge the suspension for 10 minutes at 14,000rpm at 4°C. Transfer the supernatant into a new microcentrifuge vial.
4. Add **Extraction Cleaner** to the supernatant at a 1/100 ratio (e.g., add 10  $\mu$ L of **Extraction Cleaner** to 990  $\mu$ L of the supernatant) and incubate for 15-20 minutes at room temperature.
5. Centrifuge the suspension for 1 minute at 14,000rpm at 4°C and transfer the supernatant into a new microcentrifuge vial.
6. Measure the protein concentration of the nuclear extract.
7. Use immediately or aliquot and freeze the supernatant at -80°C until further use. Avoid freeze/thaw cycle.



# 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)