



## **POLYVIEW<sup>®</sup> IHC reagent (mouse-HRP), kit**

Catalog #: ADI-950-112



# Product Manual

## **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.



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



Please contact Enzo Life Sciences Technical Support if necessary.

|                                     |   |
|-------------------------------------|---|
| Intended Use.....                   | 2 |
| Kit Contents .....                  | 2 |
| Storage .....                       | 2 |
| Stability .....                     | 2 |
| Format .....                        | 3 |
| Other Materials Needed .....        | 3 |
| Precautions .....                   | 3 |
| Recommended Staining Protocol ..... | 4 |
| Contact Information.....            | 6 |

## **INTENDED USE**

The POLYVIEW<sup>®</sup> IHC reagent (mouse-HRP), kit is a non-biotin one-step detection reagent suitable for demonstrating antigens in formalin-fixed paraffin-embedded tissues and frozen sections. The POLYVIEW<sup>®</sup> detection reagent may also be used with blood smears, cytosmears, and cell preparations.

POLYVIEW<sup>®</sup> detection kits have been developed by directly labeling anti-mouse immunoglobulins with enzymes using a proprietary tandem hyperlabelling technology. This ensures consistent and reproducible immunodetection of mouse antibodies against nuclear, cytoplasmic and membrane antigens in different types of tissues. The single step POLYVIEW<sup>®</sup> Detection Reagent enables faster staining procedures than traditional two-step methods using biotin and avidin/streptavidin conjugates, with significantly lower background.

The POLYVIEW<sup>®</sup> IHC reagent (mouse-HRP), kit is suitable for use with all mouse antibodies, both monoclonal and polyclonal. The reagents can be used for manual staining or with automated staining instruments and are well suited for multiplex immunohistochemical staining assays.

## **KIT CONTENTS**

Anti-mouse HRP detection reagent (10 mL)

## **STORAGE**

Store at 2°-8°C. Do not freeze.

## **STABILITY**

12-24 months (see expiration date on reagent bottles).

## **FORMAT**

Mouse/Rabbit POLYVIEW<sup>®</sup> HRP ready-to-use reagent is formulated without azide or thimerosal preservatives.

## **OTHER MATERIALS NEEDED**

1. Xylene or dewaxing reagents
2. Absolute ethanol
3. Distilled or deionized water
4. IHC wash buffer (Prod. no. ADI-950-235)
5. IHC background blocker (Prod. no. ADI-950-230, optional)
6. IHC diluent (primary Ab) (Prod. no. ADI-950-244)
7. Counterstain
8. HIGHDEF<sup>®</sup> IHC mount (Prod. no. ADI-950-261)
9. Chromogens HIGHDEF<sup>®</sup> IHC chromogen substrate (DAB, stable) (Prod. no. ADI-950-212) or HIGHDEF<sup>®</sup> yellow IHC chromogen (HRP) (Prod. no. ADI-950-170)
10. IHC tissue primer (Prod. no. ADI-950-234) or Peroxide Block

## **PRECAUTIONS**

1. Wear appropriate personal 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. Interpretation of the results is the sole responsibility of the user.

## RECOMMENDED STAINING PROTOCOL

1. Paraffin-embedded tissue sections must be deparaffinized with xylene or dewaxing agent and rehydrated with a graded series of ethanol and water washes before staining. Follow the standard dewaxing and rehydration protocol used in your lab.
2. The investigator needs to optimize the dilution and incubation times for primary antibodies.
3. Each immunostaining run should include known positive and negative controls to assure proper functioning of the staining system and aid in valid interpretation of the results.

### **Typical controls:**

Positive Control: A tissue known to contain the desired antigen, which has yielded positive staining in the past.

Negative Controls:

#### *Reagent Controls*

- A. Substitute normal non-immune serum from the same host animal as the primary antibody (e.g. if using mouse monoclonal primary antibodies, use mouse non-immune serum).
- B. Substitute matching host species isotype control for primary antibody
- C. Use antigen-adsorbed primary antibody (i.e. antibody reagent which has been adsorbed with the target antigen to remove specific antibody)

*Tissue control* – A tissue known to not contain the desired antigen.

4. Consult the primary antibody supplier for recommended for antigen recovery treatments. Perform epitope recovery pretreatments before starting the staining procedure.
5. Once the slide treatment has been started, DO NOT let tissues or specimens dry. This can cause undesirable background or artifacts.
6. After the Primary Antibody application wash slides with IHC wash buffer (Prod. no. ADI-950-235).
7. Apply POLYVIEW<sup>®</sup> reagent (Make sure tissue sections/smears are completely covered with the reagent) and incubate for 20 min at room temperature.
8. Wash slides with IHC wash buffer then follow the further steps with Chromogen.



# 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)