



# Product Manual

## **HIGHDEF<sup>®</sup> blue IHC chromogen (HRP)**

Catalog #: ADI-950-151

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

## INTENDED USE

Substrate/chromogen in conjunction with horseradish peroxidase (HRP)-based immunostaining or *in situ* hybridization (ISH) systems. With at least 2 hour working stability, this substrate/chromogen system can be used in any automation system as well as manual use.

## INTRODUCTION

HIGHDEF<sup>®</sup> blue IHC chromogen (HRP) is a substrate-chromogen system designed to be used for either immunohistochemistry (IHC) or ISH when utilizing horseradish peroxidase. HIGHDEF<sup>®</sup> blue IHC chromogen (HRP) produces a sharply contrasting ebony blue color that can easily be distinguished from other stains.

## KIT CONTENTS

| Description  |
|--|
| HIGHDEF <sup>®</sup> blue IHC chromogen (HRP) Substrate Buffer |
| HIGHDEF <sup>®</sup> blue IHC chromogen (HRP) Chromogen        |
| Empty Mixing Bottle  |

## STORAGE

Store at 2-8°C away from light. Do not use beyond the expiration date stated on the label.

## WORKING SOLUTION

Aliquot 1mL of HIGHDEF<sup>®</sup> blue IHC chromogen (HRP) substrate Buffer in a mixing bottle. Add one drop (~20 µL) of concentrated HIGHDEF<sup>®</sup> blue IHC chromogen (HRP) chromogen solution. Replace tip, mix, and allow solution to reach room temperature before using.

**Note:** *The HIGHDEF<sup>®</sup> blue IHC chromogen (HRP) chromogen-substrate working solution is light sensitive and should be kept away from light as much as possible. Working solution is stable for at least 2 hours, for optimal results prepare fresh reagent.*

## PROCEDURE

| Protocol   | STAINING PROCEDURE:  | INCUBATION TIMES                            |
|--|--|---|
| <p>Following peroxidase incubation, wash tissue sections with wash buffer, then follow protocol of choice:</p> |  |   |
| <p><b>Pre-Mix Working Solution: (Automation)</b></p>   | <p>HIGHDEF<sup>®</sup> blue IHC chromogen (HRP) has a working solution stability of at least 2 hours and can be loaded directly onto instrument as a single solution. Reduce exposure to light to achieve optimal staining. Working solution is applied directly to slide.</p> | <p><b>Working Solution: 10 - 20 min</b></p> |
| <p><b>On Board Mixing (Automation)</b></p>   | <p>Instruments that have on-board mixing capability can load the chromogen and substrate-buffer components independently. Working solution is made mixing reagents 1:50 in on-board mixing station before application to slide.</p>  | <p><b>Working Solution: 10 - 20 min</b></p> |
| <p><b>Manual Use</b></p>   | <p>Mix substrate-chromogen and buffer in a 1:50 ratio and apply directly to slide.</p>   | <p><b>Working Solution: 10 - 20 min</b></p> |

## COUNTERSTAIN

Counterstain with Hematoxylin or other counterstain for good contrast. Wash with distilled or de-ionized H<sub>2</sub>O followed by immuno wash buffer.

## MOUNTING

Slides should be air dried (do not dehydrate in alcohol or xylene), After rinsing off counterstain in distilled or de-ionized H<sub>2</sub>O, leave slides on benchtop for at least 20 minutes to air dry, then permanently mount. (e.g., using HIGHDEF<sup>®</sup> IHC mount, Prod. no. ADI-950-261).



# 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 NONINFRINGEMENT 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.

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

***Enabling Discovery in Life Science®***

### Global Headquarters

#### Enzo Life Sciences Inc.

10 Executive Blvd  
Farmingdale, NY 11735  
(p) 1-800-942-0430  
(f) 1-631-694-7501  
(e) [info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

#### Enzo Life Sciences (ELS) AG

Industriestrasse 17, Postfach  
CH-4415 Lause / Switzerland  
(p) +41/0 61 926 89 89  
(f) +41/0 61 926 89 79  
(e) [info-ch@enzolifesciences.com](mailto:info-ch@enzolifesciences.com)

Please visit our website at [www.enzolifesciences.com](http://www.enzolifesciences.com) for additional contact information.