



PRODUCT INFORMATION & MANUAL

BLUE HRP Staining Kit (Colorimetric) *NBP3-12182*

For research use only.
Not for diagnostic or therapeutic procedures.

www.novusbio.com - P: 303.730.1950 - P: 888.506.6887 - F: 303.730.1966 - technical@novusbio.com

Novus kits are guaranteed for 6 months from date of receipt

BLUE HRP

NBP3-12182

Effective Date: 3/31/2020

Intended Use

For In Vitro Diagnostic Use

Summary and Explanation

BLUE HRP is a substrate-chromogen system designed to be used for either IHC or ISH when utilizing horseradish peroxidase. BLUE HRP can be permanently mounted to produce a strong azure blue color that can be easily distinguished from other stains.

Principles of the Procedures

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

Reagents Provided

Kit Contents	30 mL	110 mL
BLUE HRP Substrate Buffer	30 mL	110 mL
BLUE HRP Chromogen	1 mL	3 mL
Empty Mixing Bottle	1	1

Prepare the Following Solutions Before Use

1. Aliquot 1mL of BLUE HRP Substrate Buffer in a mixing bottle.
2. Add one drop (~20µl) of concentrated BLUE HRP Chromogen solution.
3. Replace tip, mix, and allow solution to reach room temperature before using.
4. The BLUE HRP chromogen-substrate working solution is light sensitive and should be kept away from light as much as possible.
5. Working solution is stable for up to 24 hour, for optimal results prepare fresh reagent.

Materials Required But Not Provided

Some of the reagents and materials required for IHC are not provided.

Storage and Handling

Store at 2°-8°C away from light. Do not use product after the expiration date printed on vial. If reagents are stored under conditions other than those specified here, they must be verified by the user. Diluted reagents should be used promptly.

Staining Procedure

- 1) Once sections have been incubated with peroxidase, wash sections with wash buffer then follow protocol of choice:

***** bring the solutions to Room Temperature before mixing**

- a. **Pre-Mix Working Solution: (Automation)** BLUE HRP has a working solution stability of up to 24 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. Incubate for 5-10 min.
 - b. **On Board Mixing: (Automation)** 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. Incubate for 5-10 min.
 - c. **Manual Use:** Mix substrate-chromogen and buffer in a 1:50 ratio and apply directly to slide. Incubate for 5-10 min.
- 2) Counterstain with Hematoxylin or other counterstain. Wash with DI water followed by immuno wash buffer.
 - 3) Mounting: Slides should be air dried (do not dehydrate in alcohol or xylene). After rinsing off counterstain in DI H₂O, leave slides on benchtop for at least 20 minutes to air dry, and then permanently mount.

Recommendation:

For best color preservation and long term slide storage, we recommend to use Tissue Preservation Solution - HRP/AP assays (**NBP3-12178**) after counterstaining.

Precautions

- 4) Consult local and/or state authorities with regard to recommended method of disposal.
- 5) Materials of human or animal origin should be handled as biohazardous materials and disposed of with proper precautions.
- 6) Avoid microbial contamination of reagents. Contamination could produce erroneous results.
- 7) This reagent may cause irritation. Avoid contact with eyes and mucous membranes.
- 8) If reagent contacts these areas, rinse with copious amounts of water.
- 9) Do not ingest or inhale any reagents.