



**PRODUCT INFORMATION &
MANUAL**

**Tissue Preservation Solution -
HRP/AP assays
*NBP3-12178***

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

Tissue Preservation Solution - HRP/AP assays

Catalog Number: NBP3-12178

Effective Date: 3/7/2017

Intended Use

For In Vitro Diagnostic Use.

Product Description

Tissue Preservation Solution - HRP/AP assays is a preservation solution that acts as a sealant to improve tissue morphology and stain over an extended period of time. When applied to the stained section, specimens can be permanently mounted without chromogens fading. Additional benefits of the Tissue Preservation Solution - HRP/AP assays include: saves hands on time when mounting slides, avoids the use of toxic chemicals such as xylene and eliminates accumulation of hazardous waste.

Summary and Explanation

This product is designed for tissue specimens stained with peroxidase, alkaline phosphatase-based systems and various chromogens. After treating with Tissue Preservation Solution - HRP/AP assays, water solution chromogens can be mounted in Permanent Mounting Media.

Format

Clear viscous, ready to use solution.

Storage and Handling

Store at room temperature. Do not use after expiration date printed on label.

Preparation of Working Solutions

Tissue Preservation Solution - HRP/AP assays is ready to use and does not require additional preparation.

Protocol Recommendations

1. After counterstaining, rinse slide with water.
2. Position the slide at an angle on a stack of paper towels to allow excess fluid to run off.
3. Lay slides flat and pipette or add a few drops of Tissue Preservation Solution - HRP/AP assay over the tissue, making sure you cover the tissue section.
4. Let tissue section soak for about 5 minutes.
5. Position the slide at an angle on a stack of paper towels to allow excess fluid to run off.
6. Let the fluid dry completely without further rinsing for about a minimum of 20 minutes or longer if needed. It is important that the slide is completely dry before moving on to the next step.
7. Mount and coverslip slide using any conventional mounting method.

Performance Characteristics

The protocols for a specific application can vary. These include, but are not limited to: fixation, heat-retrieval method, incubation times, and tissue section thickness and detection kit used. Due to the superior sensitivity of these unique reagents, the recommended incubation times and titers listed are not applicable to other detection systems, as results may vary. Ultimately, it is the responsibility of the investigator to determine optimal conditions. These products are tools that can be used for interpretation of morphological findings in conjunction with other diagnostic tests and pertinent clinical data by a qualified pathologist.

Precautions

1. Wear disposable gloves when handling reagents.
2. Specimens, before and after fixation, and all materials exposed to them should be handled as if capable of transmitting infection and disposed of with proper precautions. Never pipette reagents by mouth and avoid contacting the skin and mucous membranes with reagents and specimens. If reagents or specimens come in contact with sensitive areas, wash with copious amounts of water.
3. Microbial contamination of reagents may result in an increase in nonspecific staining.
4. Incubation times or temperatures other than those specified may give erroneous results. The user must validate any such change.
5. Do not use reagent after the expiration date printed on the label.
6. The MSDS is available upon request.
7. Consult OSHA, federal, state or local regulations for disposal of any toxic substances.