



PRODUCT INFORMATION & MANUAL

NBP3-12176

Purple Hematoxylin Solution

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

Purple Hematoxylin Solution

Catalog Number:
NBP3-12176

Effective Date: 07/28/2020

Intended Use

For In Vitro Diagnostic Use.

Product Description

Hematoxylin is a histological staining reagent suitable for visualization of nuclei in tissue sections.

Summary and Explanation

Purple Hematoxylin Solution is intended for use in the histologic demonstration of nuclear staining. This staining technique is used to make the critical distinction between a normal nucleus and an abnormal one.

Format

Ready To Use

Volume/UOM

20mL/ 100mL

Storage and Handling

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

Preparation of Working Solutions

Purple Hematoxylin Solution is ready to use and does not require any preparation.

Protocol Recommendations

Counterstain

1. Rinse slides with deionized or distilled water.
2. Place slides into Purple Hematoxylin Solution for 30-60 seconds.
3. Rinse slides in deionized or distilled water until all free hematoxylin has been removed.
4. Bluing the slides in an alkaline solution is not required and is not recommended.

Performance Characteristics

The protocols for a specific application can vary. These include, but are not limited to: fixation, heat-retrieval method, incubation times, 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.