



PRODUCT INFORMATION & MANUAL

Tissue Glue Kit ***NBP3-12174***

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 Glue

Catalog Number: NBP3-12174

Effective Date: 8/21/2020

Intended Use

For In Vitro Diagnostic Use. This product is intended for qualitative immunohistochemistry with normal and neoplastic formalin-fixed, paraffin-embedded tissue sections, to be viewed by light microscopy. Clinical interpretation of staining results should be accompanied by histological studies with proper controls. Patients' clinical histories and other relevant diagnostic tests should be utilized by a qualified person(s) when evaluating and interpreting results.

Product Description

Tissue Glue uses a two-step process to adhere tissues onto a microscope slide. Tissue Glue is particularly useful for maintaining strong attachment of the tissue to the slide during Heat-induced Epitope Retrieval (HIER) in preparation for immunohistochemical staining. The first reagent is applied to the tissue to form attachment points on the slide and on the tissue. The second reagent contains a glue-like substance that covalently bind to the attachment points, thereby firmly attaching the tissue to the slide. Although Tissue-Glue works with all slides, positively charged slides are recommended in order to achieve the optimal results.

Summary and Explanation

This product is designed for use with both frozen tissue sections and paraffin-embedded tissue sections for adhering the tissues firmly onto a glass microscope slide.

Format

Reagent 1 and Reagent 2 are clear solutions that are provided as ready-to-use reagents and do not require any further dilutions.

Volume/UOM

50mL + 50mL

Storage and Handling

Store at 4 C. Do not use after the expiration date printed on the label.

Preparation of Working Solutions

The reagents are provided ready-to-use and no further preparation is required.

Protocol

1. This procedure will work for both frozen tissue sections and paraffin-embedded tissue sections.
2. Sections should be mounted on positively-charged glass slides for best results.
3. For paraffin-embedded sections, the slides should be deparaffinized and rehydrated using standard procedures for deparaffinization.
4. Place slides into a water bath for five minutes

5. Remove slides from water bath and carefully remove excess water by wiping around the tissue with an absorbent wipe.

6. Lay the slides flat onto a level surface and pipette 200 ul of Reagent 1 onto the slides, making sure that the tissue is completely covered with reagent.

7. Incubate slides for 10 minutes at room temperature.

8. Drain off Reagent 1 by tipping the slides onto an absorbent towel and allowing reagent to run off the slides into the towel. Do not rinse slides.

9. Lay the slides flat onto a level surface and pipette 200 ul of Reagent 2 onto the slides, making sure that the tissue is completely covered with reagent.

10. Incubate slides for 10 minutes at room temperature.

11. Drain off Reagent 2 by tipping the slides onto an absorbent towel and allow reagent to run off of the slides into the towel. Do not rinse slides.

12. Lay the slides flat onto a level surface and allow slides to completely dry. Drying time can be shortened by drying slides in a 37 C incubator.

13. The polymers contained in these reagents prevent drying artifacts, and dried slides can be stored for several days before staining.

14. At the time of staining, place the slides into an appropriate antigen retrieval solution. The tissue glue will be removed from the slides during the antigen retrieval process.

15. After antigen retrieval commence immunohistochemical staining.

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. 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. Protective gloves should be worn and work should be performed in a ventilated area.
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.