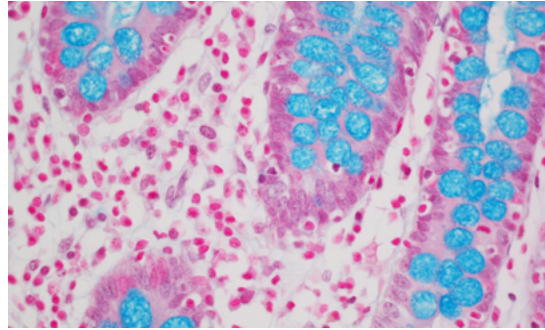


## ALCIAN BLUE (pH 2.5) STAIN KIT

**Cat. No.:** H-3501

**Storage:** 15-25 °C



Colon tissue section stained with Alcian Blue (pH 2.5) Stain Kit.

### DESCRIPTION

The Alcian Blue (pH 2.5) Stain Kit is intended for use in the histological visualization of sulfated and carboxylated acid mucopolysaccharides and sulfated and carboxylated sialomucins (glycoproteins).

Acidic Sulfated Mucosubstances:	Blue
Hyaluronic Acid:	Blue
Sialomucins:	Blue
Nuclei:	Red

Control Tissue: Small Intestine, Appendix, Colon

### COMPONENTS

<u>Product Name</u>	<u>Volume</u>
Alcian Blue Solution (pH 2.5)	30 ml
Nuclear Fast Red (Enhanced)	30 ml
Acetic Acid Solution (3%)	60 ml (30 ml x 2)

### NOTES:

- Intensity should be optimized for each tissue type, fixative used, and presentation (frozen or paraffin-embedded)
- Use standard laboratory precautions to avoid contact with skin and eyes.
- Harmful if swallowed. Do not take internally.
- These reagents are for research use only.
- Follow all Federal, State and Local regulations for disposal.

### PROCEDURE

- Deparaffinize sections, if required, and hydrate to distilled water.
- Apply Acetic Acid Solution onto tissue section for 3 minutes.
- Tip off excess and apply Alcian Blue Solution for 30 minutes at room temperature or 15 minutes at 37 °C.
- Rinse slide briefly (10-30 seconds) in Acetic Acid Solution to remove excess Alcian Blue.
- Rinse for 2 minutes in running tap water followed by 2 changes of distilled water (1-2 minutes each).
- Apply Nuclear Fast Red Solution for 5 minutes.
- Rinse for 2 minutes in running tap water followed by 2 changes of distilled water (1-2 minutes each).
- Dehydrate through graded ethanols.
- Clear and coverslip.

### Related Reagents

<u>Product Name</u>	<u>Unit Size</u>	<u>Cat. No.</u>
Vector® Hematoxylin	500 ml	H-3401
Vector Hematoxylin QS	100 ml	H-3404
Vector Methyl Green	500 ml	H-3402
Vector Nuclear Fast Red	500 ml	H-3403
Hematoxylin & Eosin Stain Kit	1 Kit	H-3502
VectaMount® Mounting Medium	60 ml	H-5000
ImmPrint™ Histology Pen	5-pen set	H-6100