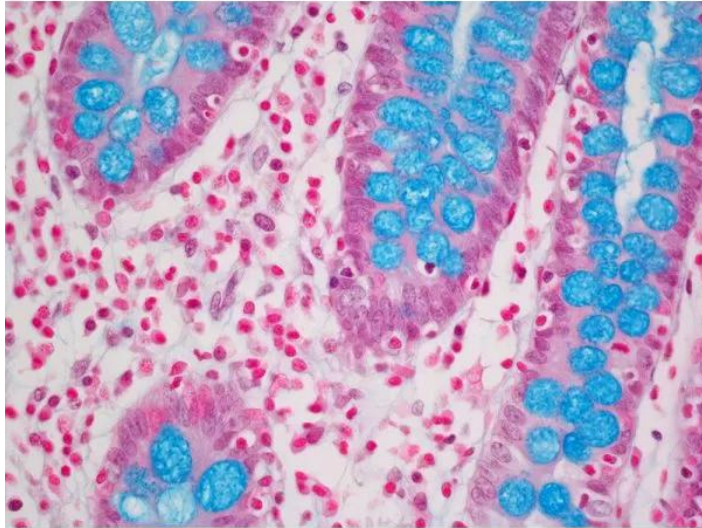




ALCIAN BLUE (PH 2.5) STAIN KIT

SKU: H-3501



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

SPECIFICATIONS

Format

Ready-to-Use

Microscopy

Brightfield

For research use only. Not intended for therapeutic or diagnostic use in animals or humans.



Unit Size	1 Kit
Applications	Histology
Cellular Stain	Nucleus, Glycoproteins

KIT COMPONENTS

- Alcian Blue Solution (pH 2.5), 30 ml
- Nuclear Fast Red (Enhanced), 30 ml
- Acetic Acid Solution (3%), 60 ml (30 ml x 2)

TECHNICAL INFORMATION

Components:

- 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.

CITATIONS



Powered by Bioz © 2023 See more details on Bioz

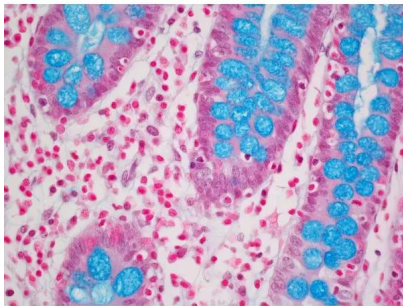
For research use only. Not intended for therapeutic or diagnostic use in animals or humans.



DOCUMENTS

- [User Guide](#)
- [Safety Data Sheet](#)
- [Download CoA](#)
- [Datasheet](#)

GALLERY IMAGES



For research use only. Not intended for therapeutic or diagnostic use in animals or humans.