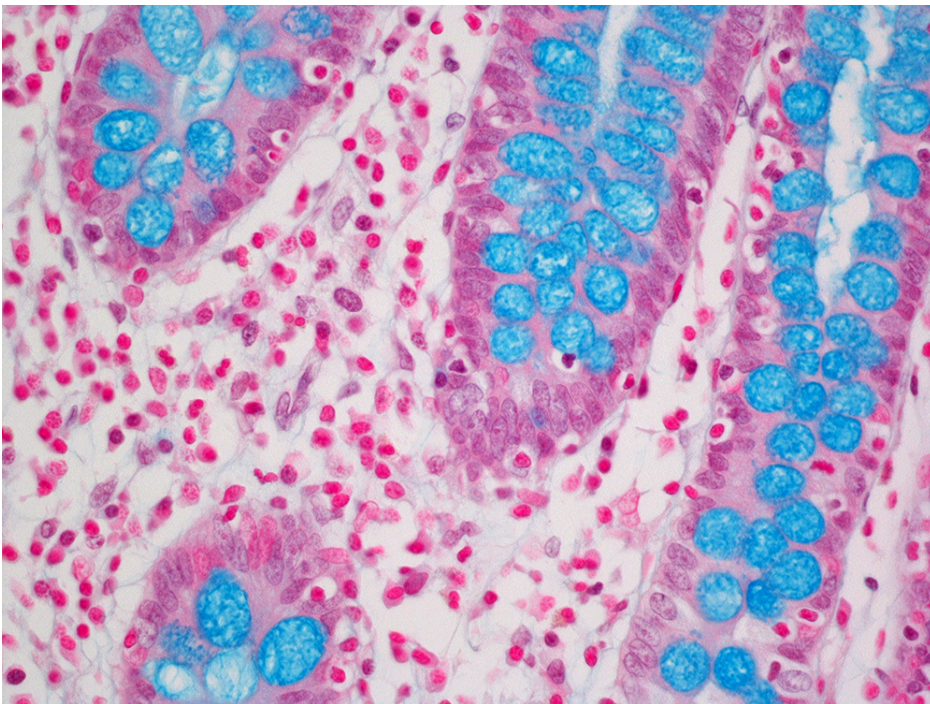




Alcian Blue (pH 2.5) Stain Kit

H-3501

Product Images



Short 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

Additional Information

Unit Size	1 Kit
Applications	Histology
Microscopy	Brightfield
Cellular Stain	Nucleus, Glycoproteins
Color	Blue, Red
Format	Ready-to-Use

