

# Alcian Blue Stain Kit (pH 1.0, Mucin Stain) ab150661

[2 References](#)   [2 Images](#)

Overview

Product name

Alcian Blue Stain Kit (pH 1.0, Mucin Stain)

Product overview

The Alcian Blue (pH 1.0) (Mucin Stain) is intended for use in the histological visualization of strongly sulfated mucosubstances.

The component Acetic Acid Solution has been replaced with Hydrochloric Acid Solution in this kit.

Alcian Blue staining protocol summary:

- deparaffinize sections if necessary and hydrate
- prepare Working Rinse solution
- apply Working Rinse solution to tissue for 30 sec
- incubate in Alcian Blue pH 1.0 for 30 min
- rinse in Working Rinse Solution
- blot and air dry slide
- stain in nuclear fast red for 1-2 min
- rinse in water and allow to dry completely
- clear and mount

Other products for staining tissue sections

Find more kits and reagents in the [special stains guide](#), or products for antigen retrieval, blocking, signal amplification, visualization, counterstaining, and mounting in the [IHC kits and reagents guide](#).

Notes

Staining Interpretation

Strongly Sulfated Mucosubstances	Blue
Nuclei	Red
Background	Pink

**Control Tissue:** Small Intestine, Appendix, Colon

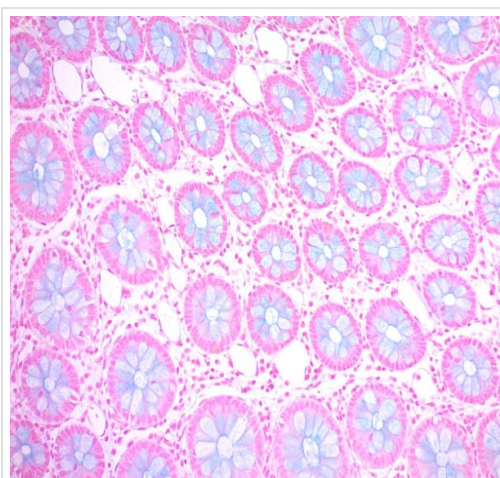
Previously called Alcian Blue pH 1.0 Stain Kit (Mucin Stain)

## Storage instructions

Store at room temperature. Please refer to protocols.

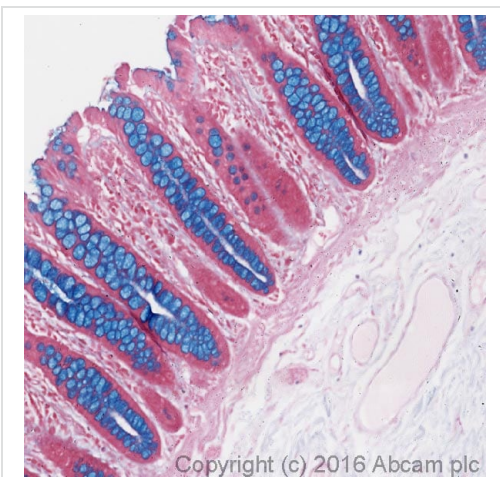
Components	100 tests
1M Hydrochloric Acid Solution	1 x 500ml
Alcian Blue pH 1	1 x 250ml
Nuclear Fast Red Solution (Enhanced)	1 x 250ml

## Images



ab150661 - Alcian Blue pH 1.0 Stain Kit (Mucin Stain)

ab150661 Alcian Blue pH 1 Stain Kit (Mucin Stain) staining formalin-fixed-paraffin embedded human GI tract.



ab150661 - Alcian Blue pH 1.0 Stain Kit (Mucin Stain)

ab150661 (Alcian Blue pH 1.0) staining strongly sulphated mucins (blue) in formalin fixed paraffin embedded human colon, counterstained with safranin o (pink).

\*Tissue obtained from the Human Research Tissue Bank, supported by the NIHR Cambridge Biomedical Research Centre.

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

### **Our Abpromise to you: Quality guaranteed and expert technical support**

---

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
  
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

### **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors