

Version 4b Last updated 20 May 2020

ab150662

Alcian Blue, pH 2.5 (Mucin Stain)

For the histological visualization of sulfated and carboxylated acid mucopolysaccharides and sialomucins (glycoproteins).

[View kit datasheet: www.abcam.com/ab150662](http://www.abcam.com/ab150662)

(use www.abcam.cn/ab150662 for China, or www.abcam.co.jp/ab150662 for Japan)

This product is for research use only and is not intended for diagnostic use.

Table of Contents

1. Overview	1
2. Materials Supplied and Storage	2
3. Materials Required, Not Supplied	3
4. General guidelines, precautions, and troubleshooting	4
5. Staining Protocol	5
6. FAQs / Troubleshooting	6
7. Typical Data	7
8. Notes	8

1. Overview

The Alcian Blue, pH 2.5 (Mucin Stain) (ab150662) is intended for use in the histological visualization of sulfated and carboxylated acid mucopolysaccharides and sulfated and carboxylated sialomucins (glycoproteins).

Staining Interpretation:

Acidic Sulfated Mucosubstances	Blue
Hyaluronic Acid	Blue
Sialomucins	Blue
Nuclei	Red
Background	Pink

Control Tissue:

Small intestine, Colon, Appendix.

2. Materials Supplied and Storage

Store kit at room temperature immediately on receipt and check below for storage for individual components. Kit can be stored for 1 year from receipt, if components have not been reconstituted.

Keep away from open flame and refer to the safety datasheet.

ΔNote: This Kit originally used Safranin O Solution but this has now been replaced with Nuclear Fast Red Solution.

Item	Quantity	Storage temperature (before prep)
Alcian Blue, pH 2.5	250 mL	RT
Nuclear Fast Red Solution	250 mL	RT
Acetic Acid Solution	500 mL	RT

3. Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully perform this assay:

- Graded alcohol solutions
- Distilled water

4. General guidelines, precautions, and troubleshooting

Please observe safe laboratory practice and consult the safety datasheet.

For general guidelines, precautions, limitations on the use of our assay kits and general assay troubleshooting tips, particularly for first time users, please consult our guide:

www.abcam.com/assaykitguidelines

For typical data produced using the assay, please see the assay kit datasheet on our website.

5. Staining Protocol

- Equilibrate all materials and prepared reagents to room temperature just prior to use and gently agitate.

- 5.1** Deparaffinize sections if necessary and hydrate in distilled water.
- 5.2** Incubate slide in Acetic Acid Solution for 3 minutes.
- 5.3** Incubate slide in Alcian Blue (pH 2.5) solution for 30 minutes at room temperature or 15 minutes at 37° C.
- 5.4** If desired, rinse slide briefly in Acetic acid solution to remove excess Alcian Blue.
- 5.5** Rinse for 2 minutes in running tap water followed by 2 changes of distilled water.
- 5.6** Stain slide in Nuclear Fast Red Solution for 5 minutes.
- 5.7** Rinse for 2 minutes in running tap water followed by 2 changes of distilled water.
- 5.8** Dehydrate through graded alcohols.
- 5.9** Clear and mount in synthetic resin.

6. FAQs / Troubleshooting

General troubleshooting points are found at www.abcam.com/assaykitguidelines.

7. Notes

Technical Support

Copyright © 2020 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

www.abcam.com/contactus

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)