

ab245876

Alcian Blue PAS Stain Kit

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

View Alcian Blue PAS Stain Kit datasheet:

<http://www.abcam.com/ab245876>

[use <http://www.abcam.cn/ab245876> for China, or <http://www.abcam.co.jp/ab245876> 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. General guidelines, precautions, and troubleshooting	3
4. Assay Procedure	4
5. Notes	5

1. Overview

The Alcian Blue PAS Stain Kit is intended for use in the simultaneous histological visualization of sulfated and carboxylated acid mucopolysaccharides, sulfated and carboxylated sialomucins (glycoproteins) and neutral mucins.

Staining Interpretation:

Acidic Sulfated Mucosubstances:	Blue
Hyaluronic Acid:	Blue
Sialomucins:	Blue
Neutral Mucins:	Magenta
Mixtures of Acidic and Neutral Mucins:	Blue - Mauve depending on dominant entity

Control Tissue:

Small or large 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.

Item	Quantity (30 mL pack size)	Quantity (1 kit/ 250 mL pack size)	Storage temperature (before prep)	Storage temperature (after prep)
Acetic Acid Solution (3%)	30 mL	2 x 125 mL	RT	RT
Alcian Blue Solution, pH 2.5	30 mL	250 mL	RT	RT
Periodic Acid Solution (1%)	30 mL	250 mL	4°C	4°C
Schiff's Solution	30 mL	250 mL	4°C	4°C
Hematoxylin (modified Mayer's solution)	30 mL	2 x 125 mL	RT	RT

3. 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.

4. Assay Procedure

- Equilibrate all materials and prepared reagents to room temperature just prior to use and gently agitate.
- 1.1 Deparaffinize sections if necessary and hydrate in distilled water.
- 1.2 Apply Acetic Acid Solution (3%) to tissue section for 2 minutes.
- 1.3 Remove excess Acetic Acid Solution (3%) and without rinsing apply Alcian Blue (pH 2.5) solution for 15-20 minutes.
- 1.4 Rinse for 2 minutes in running tap water followed by 2 changes of distilled water.
- 1.5 Apply Periodic Acid Solution to tissue section for 5 minutes.
- 1.6 Rinse slide in 2 changes of distilled water.
- 1.7 Apply Schiff's Solution to tissue section for 10-20 minutes.
- 1.8 Rinse for 2 minutes in warm running tap water followed by 2 changes of distilled water.
- 1.9 Apply Hematoxylin (modified Mayer's solution) to tissue section for 2 minutes.
- 1.10 Rinse for 2 minutes in running tap water followed by 2 changes of distilled water.
- 1.11 Dehydrate through graded alcohols.
- 1.12 Clear, and mount in synthetic resin.

5. Notes

Technical Support

Copyright © 2025 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)