

ab150680 - Periodic Acid Schiff (PAS) Stain Kit (Mucin Stain)

For the histological identification of lymphocytes and mucopolysaccharides.

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

Overview

The Periodic Acid Schiff (PAS) Stain Kit (Mucin Stain) is intended for use in histological demonstration of lymphocytes and mucopolysaccharides.

The staining pattern of the lymphocytes are helpful in making therapeutic decisions in established cases of lymphocytic leukemia. The PAS reaction in tissue sections is useful for the demonstration of mucopolysaccharides. PAS staining may also be used for the demonstration of fungal organisms in tissue sections.

Staining Interpretation:

Fungal Organisms	Magenta
Pas positive Material	Magenta
Nuclei	Black/Blue
Other Tissue Components	Green/Blue

Control Tissue: Kidney, intestine and liver.

Materials Supplied and Storage

Store kit at Room temperature immediately on receipt and check below for storage for individual components.

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

Item	Quantity (30 mL pack size)	Quantity (100 test/ 125 mL pack size)	Storage temperature (before prep)
Hematoxylin (Modified Mayer's)	30 mL	125 ml	RT
Periodic Acid Solution	30 mL	250 ml	4°C
Schiff's Solution	30 mL	250 ml	4°C
Bluing Reagent	30 mL	125 ml	RT
Light Green Solution	30 mL	125 ml	RT

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

Assay Procedure

- Equilibrate all materials and prepared reagents to room temperature just prior to use and gently agitate.
 1. Deparaffinize sections if necessary and hydrate in distilled water.
 2. If sections are Zenker-fixed, remove mercuric chloride crystals using iodine and clear with sodium thiosulfate. Rinse in running tap water.
 3. Immerse slide in Periodic Acid Solution for 5 minutes (10 minutes for kidney, skin and diastase digested liver sections).
 4. Rinse slide in four changes of distilled water.
 5. Immerse slide in Schiff's Solution for 15 minutes (30 minutes for kidney, skin and diastase digested liver sections).
 6. Rinse slide in hot running tap water.
 7. Rinse slide in distilled water.
 8. Stain slide in Hematoxylin (Modified Mayer's) for 2-3 minutes.
 9. Rinse slide in running tap water for 2-3 minutes.
 10. Apply Bluing Reagent for 30 seconds.
 11. Rinse in distilled water.
 12. Apply Light Green Solution for 2 minutes (optional).
 13. Rinse quickly in distilled water.
 14. Dehydrate through graded alcohols.
 15. Clear slide and mount in synthetic resin.

Technical Support

Copyright © 2026 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:

<https://www.abcam.com/en-us/contact-us>

<https://www.abcam.cn/contact-us> (China)

<https://www.abcam.co.jp/contact-us> (Japan)