

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

ab150680

[68 References](#) [2 Images](#)

Overview

Product name	Periodic Acid Schiff (PAS) Stain Kit (Mucin Stain)
Sample type	Tissue sections
Assay time	0h 45m
Product overview	<p>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.</p>

- PAS staining protocol summary:
- deparaffinize sections (if necessary) and hydrate
 - immerse slide in PAS solution for 5-10 min
 - rinse 4 times with water
 - immerse slide in Schiff's solution for 15-30 min and rinse
 - stain slide with Hematoxylin for 2-3 min and rinse
 - incubate in Bluing reagent for 30 s and rinse
 - incubate in Light Green Solution for 2 min and rinse
 - dehydrate with alcohol, 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

Fungal Organisms	Magenta
PAS Positive Material	Magenta
Nuclei	Black/Blue

Other Tissue Components

Green/Blue

Control Tissue: Kidney, Intestine, Liver.

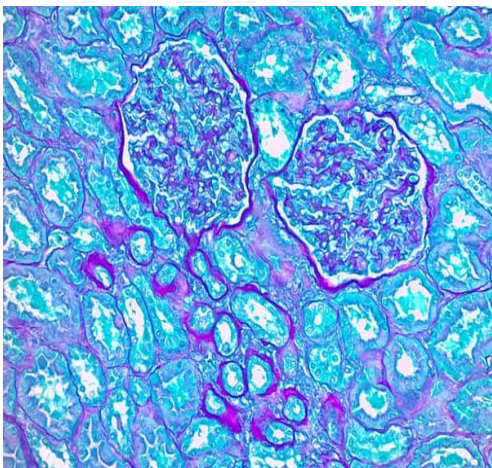
Properties

Storage instructions

Store at +4°C. Please refer to protocols.

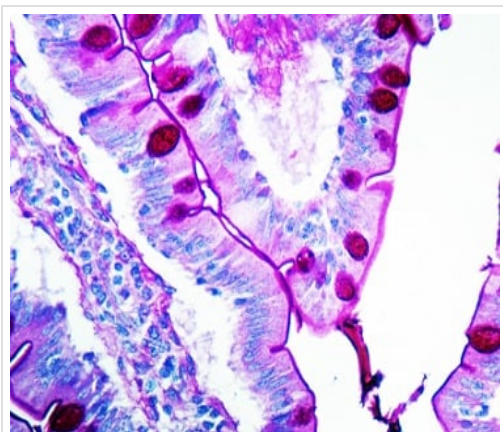
Components	100 tests
Bluing Reagent	1 x 125ml
Hematoxylin (Modified Mayer's Solution)	1 x 125ml
Light Green Solution	1 x 125ml
Periodic acid solution	1 x 250ml
Schiff's solution	1 x 250ml

Images



ab150680 Periodic Acid Schiff (PAS) Stain Kit (Mucin Stain) staining formalin-fixed-paraffin embedded human kidney. Showing fungal organisms (magenta), PAS Positive Material (Magenta), Nuclei (Black/Blue) and other tissue components (Green/Blue).

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



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

Staining of Intestine using ab150680 - Periodic Acid Schiff (PAS) Stain Kit. Showing fungal organisms (magenta), PAS Positive Material (Magenta), Nuclei (Black/Blue) and other tissue components (Green/Blue).

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