

Version 2 Last updated 24 January 2019

ab245886

PAS Stain Kit (Fungus, Periodic Acid Schiff Stain)

For the histological visualization of fungal organisms in tissue sections

View PAS Stain Kit datasheet:

<https://www.abcam.com/ab245886>

[use www.abcam.cn/ab245886 for China, or www.abcam.co.jp/ab245886 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 PAS Stain Kit (Fungus, Periodic Acid Schiff Stain) is intended for use in histological demonstration of fungal organisms in tissue sections. The PAS reaction is also useful in the demonstration of lymphocytes and mucopolysaccharides. The staining patterns of the lymphocytes are helpful in making therapeutic decisions in established cases of lymphocytic leukemia.

Staining Interpretation:

Fungal Organisms:	Magenta
PAS Positive Material:	Magenta
Other Tissue Components :	Green/Blue

Control Tissue:

Any fungal infected tissue, Kidney, intestine and liver tissues.

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	Storage temperature (before prep)	Storage temperature (after prep)
Periodic Acid Solution	250 mL	2-8° C	2-8° C
Schiff's Solution	250 mL	2-8° C	2-8° C
Light Green Solution	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 to distilled water.
- 1.2 If sections are Zenker-fixed, remove mercuric chloride crystals using iodine and clear with sodium thiosulfate. Rinse in running
- 1.3 tap water.
- 1.4 Immerse slide in Periodic Acid Solution for 5 minutes (10 minutes for Kidney, skin and diastase digested liver sections).
- 1.5 Rinse slide in 4 changes of distilled water.
- 1.6 Immerse slide in Schiff's Solution for 15 minutes (30 minutes for Kidney, skin and diastase digested liver sections).
- 1.7 Rinse slide in hot running tap water.
- 1.8 Rinse slide in distilled water.
- 1.9 Stain slide in Light Green Solution (LGA125) for 2 minutes.
- 1.10 Rinse slide using absolute alcohol.
- 1.11 Dehydrate in 2 changes of absolute alcohol, clear, and mount in synthetic resin.

ΔNote: A crystal precipitate may be seen when staining with small volumes of Schiff's solution on horizontal slides. This precipitate can be removed by rinsing vigorously in warm tap water for 5 minutes or by reapplying Periodic Acid Solution to the tissue and agitating the slide for 30-60 seconds. These modifications should be performed before counterstaining.

5. Notes

Technical Support

Copyright © 2019 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

Austria

wissenschaftlicherdienst@abcam.com | 019-288-259

France

supportscientifique@abcam.com | 01.46.94.62.96

Germany

wissenschaftlicherdienst@abcam.com | 030-896-779-154

Spain

soportecientifico@abcam.com | 91-114-65-60

Switzerland

technical@abcam.com

Deutsch: 043-501-64-24 | Français: 061-500-05-30

UK, EU and ROW

technical@abcam.com | +44(0)1223-696000

Canada

ca.technical@abcam.com | 877-749-8807

US and Latin America

us.technical@abcam.com | 888-772-2226

Asia Pacific

hk.technical@abcam.com | (852) 2603-6823

China

cn.technical@abcam.com | 400 921 0189 | +86 21 2070 0500

Japan

technical@abcam.co.jp | +81-(0)3-6231-0940

Singapore

sg.technical@abcam.com | 800 188-5244

Australia

au.technical@abcam.com | +61-(0)3-8652-1450

New Zealand

nz.technical@abc.com | +64-(0)9-909-7829