

Version 2 Last updated 11 September 2019

ab245881 Orcein Stain Kit (Hepatitis B and Elastic Fibers)

For use in histology procedures for the visualization of Hepatitis B surface Antigen (HBsAg), elastic fibers, and copper associated proteins.

View Orcein Stain Kit (Hepatitis B and Elastic Fibers) datasheet:

www.abcam.com/ab245881

[use www.abcam.cn/ab245881 for China, or www.abcam.co.jp/ab245881 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 Orcein Stain Kit (Hepatitis B and Elastic Fibers) is intended for use in histology procedures for the visualization of Hepatitis B surface Antigen (HBsAg), elastic fibers, and copper associated proteins.

Staining Interpretation:

HBsAg:	Dark Brown/Purple
HBsAg:	Dark Brown/Purple
Copper Assoc. Proteins:	Dark Purple
Background:	Light Reddish/Purple

Control Tissue:

Known hepatitis positive liver. Lung for elastic fiber. Any well fixed tissue cut 3-5 microns.

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)
Potassium Permanganate Sol. (5%)	30 mL	RT	RT
Sulfuric Acid Solution (3%)	30 mL	RT	RT
Oxalic Acid Solution (2%)	125 mL	RT	RT
Orcein Solution	125 mL	RT	RT
Differentating 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 Prepare the Oxidizing Solution immediately prior to beginning procedure by combining:

Reagent	Volume
Distilled water	50 mL
Potassium Permanganate Solution (5%)	5 mL
Sulfuric Acid Solution (3%)	3 mL

Mix thoroughly.

- 1.2 Deparaffinize sections if necessary and hydrate in distilled water.
- 1.3 Incubate slide in freshly prepared Oxidizing Solution for 10 mins.
- 1.4 Rinse slide briefly in running tap water followed by one dip in distilled water.
- 1.5 Incubate slide in Oxalic Acid Solution (2%) for 10 mins.
Δ Note: Section should be colorless following this step.
- 1.6 Rinse slide for 1 min in running tap water followed by two dips in distilled water.
- 1.7 Apply adequate Orcein Solution to cover tissue and incubate slide in Orcein Solution for 2 hrs.
Δ Note: Check periodically and apply additional stain as needed to avoid drying.
- 1.8 Rinse slide in Alcohol, Reagent (70%).
- 1.9 Apply Differentiating Solution for 3-5 secs.
- 1.10 Dip slide in Alcohol, Reagent (70%) and check slide microscopically for proper differentiation.
Δ Note: Repeat step 1.9 if necessary.
- 1.11 Dehydrate quickly in three changes of absolute alcohol.
- 1.12 Clear, and mount in synthetic resin.

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