

## ab245880– H&E Stain Kit (Hematoxylin and Eosin)

For use in histology and cytology applications

**View Datasheet:** [www.abcam.com/ab245880](http://www.abcam.com/ab245880) [use [www.abcam.cn/ab245880](http://www.abcam.cn/ab245880) for China, or [www.abcam.co.jp/ab245880](http://www.abcam.co.jp/ab245880) for Japan]

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

### Overview

H&E Stain Kit (Hematoxylin and Eosin) ab245880 is intended for use in histology and cytology applications. Included in this kit is a newly formulated Eosin that provides the benefits of a traditional alcoholic formula with significant improvements in usability. Advantages include lower evaporation rate, better color patterns, reduced tendency to spill over container, hands, and countertops, and improved surface tension to remain on tissue section. Our Hematoxylin produces crisp, intense blue nuclei providing optimal contrast to the Eosin stained cytoplasm.

### Staining Interpretation:

Cytoplasm: Light pink  
Collagen: Pink  
Muscle: Pink/Rose  
Erythrocytes: Pink/Red  
Nuclei: Blue

### Control Tissue:

Any well fixed paraffin embedded or frozen tissue section. Cell smear.

### Precautions

**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](http://www.abcam.com/assaykitguidelines)

For typical data produced using the assay, please see the assay kit datasheet on our website.

### Materials Supplied:

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)
Hematoxylin, Mayer's (Lillie's Modification)	500 mL	RT	RT
Bluing Reagent	500 mL	RT	RT
Eosin Y Solution (Modified Alcoholic)	500 mL	RT	RT

### 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. Apply adequate Hematoxylin, Mayer's (Lillie's Modification) to completely cover tissue section and incubate for 5 mins.
3. Rinse slide in two changes of distilled water to remove excess stain.
4. Apply adequate Bluing Reagent to completely cover tissue section and incubate for 10-15 secs.
5. Rinse slide in two changes of distilled water.
6. Dip slide in absolute alcohol and blot excess off.
7. Apply adequate Eosin Y Solution (Modified Alcoholic) to completely cover tissue section to excess and incubate for 2-3 mins.
8. Rinse slide using absolute alcohol.
9. Dehydrate slide in three changes of absolute alcohol.
10. Clear slide and mount in synthetic resin

### Technical Support

Copyright © 2020 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](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)