

Version 3b Last updated 21 April 2026

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

For use in histology and cytology applications

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

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.

## 2. 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.

| <b>Item</b>                                     | <b>Quantity<br/>(30mL<br/>pack<br/>size)</b> | <b>Quantity<br/>(1 kit/<br/>500mL<br/>pack<br/>size)</b> | <b>Storage<br/>temp<br/>(before<br/>prep)</b> | <b>Storage<br/>temp<br/>(after<br/>prep)</b> |
|---|--|--|---|--|
| Hematoxylin, Mayer's<br>(Lillie's Modification) | 30 mL  | 500 mL   | RT  | RT   |
| Bluing Reagent                                  | 30 mL  | 500 mL   | RT  | RT   |
| Eosin Y Solution<br>(Modified Alcoholic)        | 30 mL  | 500 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](http://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 in distilled water.
- 1.2** Apply adequate Hematoxylin, Mayer's (Lillie's Modification) to completely cover tissue section and incubate for 5 mins.
- 1.3** Rinse slide in two changes of distilled water to remove excess stain.
- 1.4** Apply adequate Bluing Reagent to completely cover tissue section and incubate for 10-15 secs.
- 1.5** Rinse slide in two changes of distilled water.
- 1.6** Dip slide in absolute alcohol and blot excess off.
- 1.7** Apply adequate Eosin Y Solution (Modified Alcoholic) to completely cover tissue section to excess and incubate for 2-3 mins.
- 1.8** Rinse slide using absolute alcohol.
- 1.9** Dehydrate slide in three changes of absolute alcohol.
- 1.10** Clear slide and mount in synthetic resin.

## 5. Notes

## **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)