

H&E Staining Kit (Hematoxylin and Eosin) ab245880

★★★★★ [3 Abreviews](#) [129 References](#) [3 Images](#)

Overview

Product name H&E Staining Kit (Hematoxylin and Eosin)

Sample type Tissue sections

Product overview H&E Staining 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. The Hematoxylin produces crisp, intense blue nuclei providing optimal contrast to the Eosin stained cytoplasm.

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.

Notes The **Hematoxylin solution ab220365** and **Eosin Y Solution ab246824** included in this kit are also available as separate products.

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](#).

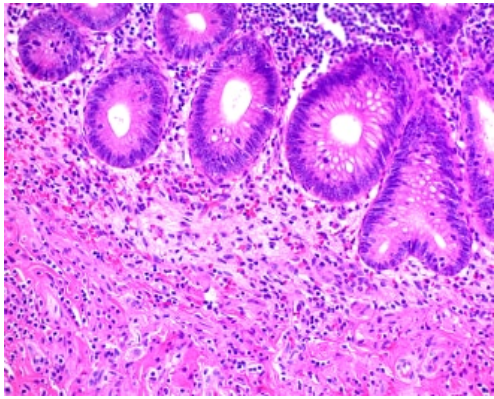
Properties

Storage instructions Store at room temperature. Please refer to protocols.

Components	1 kit
Bluing Reagent	1 x 500ml
Eosin Y Solution	1 x 500ml

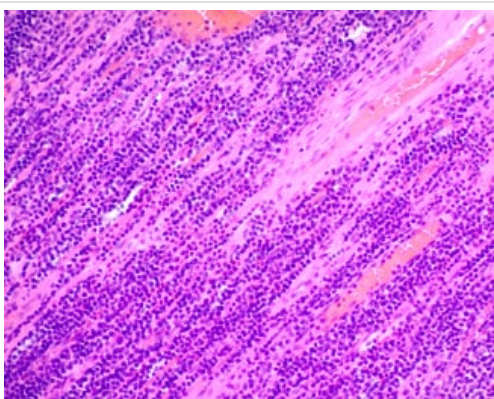
Components	1 kit
Hematoxylin (Modified Mayer's Solution)	1 x 500ml

Images



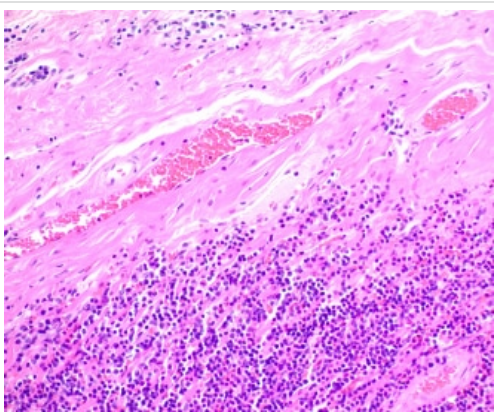
H&E Staining Kit (Hematoxylin and Eosin) ab245880

ab245880 H&E Staining Kit used to perform Hematoxylin and Eosin staining of formalin-fixed-paraffin embedded tissue sections.



H&E Staining Kit (Hematoxylin and Eosin) ab245880

ab245880 H&E Staining Kit used to perform Hematoxylin and Eosin staining of formalin-fixed-paraffin embedded tissue sections.



H&E Staining Kit (Hematoxylin and Eosin) ab245880

ab245880 H&E Staining Kit used to perform Hematoxylin and Eosin staining of formalin-fixed-paraffin embedded tissue sections.

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