



## VitroView™ H&E Stain Kit

Cat. No. VB-3000

### Introduction

Hematoxylin and Eosin is the most widely used and important general purpose histological stain combination. This specially formulated kit provides clear and sharp blue nuclear and red cytoplasmic staining with very low background and with no need to perform differentiation steps. Mayer's Hematoxylin solution is also good for immunohistochemistry counterstaining when DAB or AEC substrates are used.

### Kit Components

1. Mayer's Hematoxylin Solution - 250 ml
2. Eosin Solution - 250 ml

### Storage

Room temperature.

### Protocol

#### H&E Staining of paraffin sections:

1. Deparaffinize in xylene I for 6 minutes and II for 6 minutes.
2. Rehydrate
  - a. Ethanol 100% (2 minutes)
  - b. Ethanol 100% (2 minutes)
  - c. Ethanol 95% (2 minutes)
  - d. Ethanol 95% (2 minutes)
  - e. Ethanol 70% (2 minutes)
3. Rinse in DI water (2 minutes).
4. Place in Mayer's Hematoxylin Solution for 5 minutes.
5. Rinse in tap water for 5 min with 3 changes.
6. Place in 95% Ethanol for 1 minute (DO NOT RINSE).
7. Place in Eosin for 1 min 30sec.
8. Dehydrate with 2 changes of 95% Ethanol and 2 changes of 100% Ethanol (2 minute per change).
9. Clear with 3 changes of xylene (5 minute per change)
10. Mount coverslip onto glass slide with Permount or some other suitable organic mounting medium.

#### H&E Staining of frozen sections:

1. Frozen sections are fixed in acetone at -20 °C for 3 min or 80% Methanol at 4°C for 5 min
2. PBS 1X 10 min.
3. Rinse in DI water (2 minutes).
4. Follow step 3-12 above.

#### For Immunohistochemistry (IHC) Counterstaining:

Prepare a 1:10 dilution with DI Water and use for IHC counterstain for 1 minute.

**Precautions:** Handle with care. Avoid contact with eyes, skin and clothing. Do not ingest. Wear gloves.