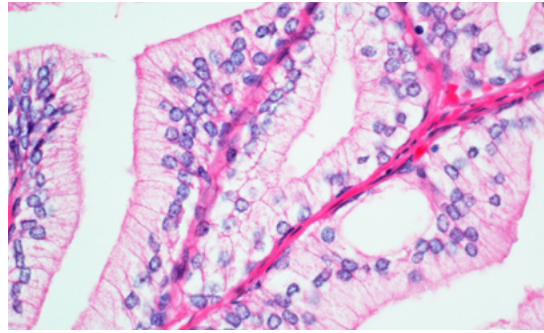


HEMATOXYLIN & EOSIN STAIN KIT

Cat. No.: H-3502

Storage: 15-25 °C



Prostate tissue section stained with Hematoxylin & Eosin Stain Kit.

DESCRIPTION

The Hematoxylin and Eosin Stain Kit is intended for use in histology and cytology applications. Included in this kit is a modified Eosin that provides the benefits of a traditional alcoholic formulation with significant improvements in usability. Advantages include lower evaporation rate, better color patterns, and improved surface tension to remain on tissue section. Our Hematoxylin produces crisp, intense blue nuclei providing optimal contrast to the Eosin stained cytoplasm.

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

Control Tissues: Small Intestine, Appendix, Colon, Prostate

COMPONENTS

<u>Product Name</u>	<u>Volume</u>
Hematoxylin, Mayer's (Modified)	30 ml
Bluing Reagent	30 ml
Eosin Y Solution (Modified)	30 ml

NOTES:

- a) Intensity should be optimized for each tissue type, fixative used, and presentation (frozen or paraffin-embedded)
- b) Use standard laboratory precautions to avoid contact with skin and eyes.
- c) Harmful if swallowed. Do not take internally.
- d) These reagents are for research use only.
- e) Follow all Federal, State and Local regulations for disposal.

PROCEDURE

1. Deparaffinize sections, if required, and hydrate to distilled water.
2. Apply adequate Hematoxylin to completely cover tissue section and incubate for 5 minutes.
3. Rinse slide in 2 changes of distilled water (15 seconds each) to remove excess stain.
4. Apply adequate Bluing Reagent to completely cover tissue section and incubate for 10-15 seconds.
5. Rinse for slide in 2 changes of distilled water (15 seconds each).
6. Dip slide in 100% ethanol (10 seconds) and blot excess off.
7. Apply adequate Eosin Y Solution to completely cover tissue section and incubate for 2-3 minutes.
8. Rinse slide using 100% ethanol (10 seconds).
9. Dehydrate slide in 3 changes of 100% ethanol (1-2 minutes each).
10. Clear and coverslip.

Related Reagents

<u>Product Name</u>	<u>Unit Size</u>	<u>Cat. No.</u>
Vector Hematoxylin	500 ml	H-3401
Vector Hematoxylin QS	100 ml	H-3404
Vector Methyl Green	500 ml	H-3402
Vector Nuclear Fast Red	500 ml	H-3403
Alcian Blue (pH 2.5) Stain Kit	1 Kit	H-3501
VectaMount [®] Mounting Medium	60 ml	H-5000
ImmPrint [™] Histology Pen	5-pen set	H-6100