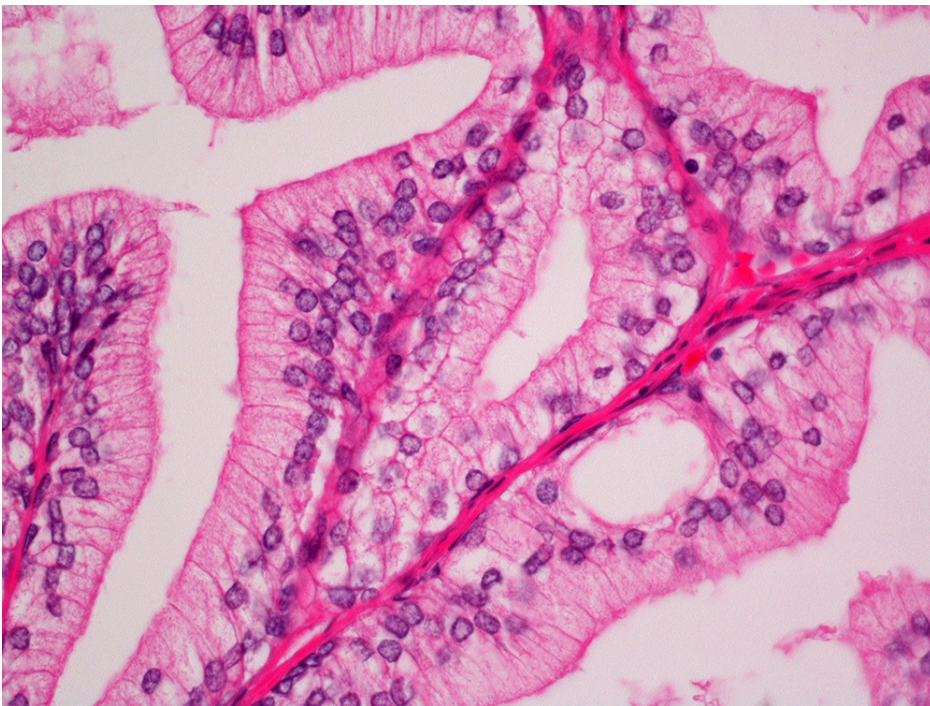




# Hematoxylin and Eosin Stain Kit

## H-3502

### Product Images



## Short 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.

## Additional Information

---

Unit Size	1 Kit
Applications	Histology
Microscopy	Brightfield
Cellular Stain	Nucleus, Cytoplasm
Color	Blue, Pink
Format	Ready-to-Use

