



---

**VECTOR®**  
**HEMATOXYLIN QS**  
**NUCLEAR COUNTERSTAIN**  
**(Modified Mayer's Formula)**  
Catalog Number H-3404

This nuclear counterstain was designed and developed to provide quick, convenient, single step staining and to complement immunocytochemical procedures. Nuclei in stained sections will be a blue-violet color with crisp nuclear detail.

**INSTRUCTIONS FOR USE:**

1. Rinse slides in tap water.
2. Immerse slides in Vector® Hematoxylin QS counterstain or apply counterstain directly to slide, completely covering the tissue section.
3. Incubate sections with Vector® Hematoxylin QS for 5-45 seconds.
4. Rinse sections with running tap water until rinse water is colorless or dip for approximately 10 seconds in water.
5. Mount sections in aqueous or permanent mounting media as per standard protocols.

**NOTES:**

Counterstain intensity should be optimized for each tissue type, antigen unmasking protocol, and immunocytochemical staining intensity desired.

Generally, hematoxylin solutions produce a more intense nuclear stain with age. Staining times can be adjusted to achieve the desired nuclear intensity.

See reverse side for substrate compatibility.

Use ordinary precautions to avoid contact with skin and eyes.

This product is for research use only.