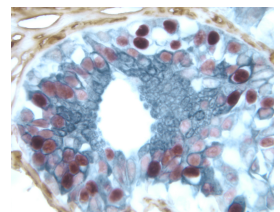
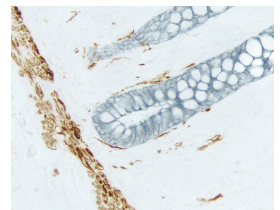
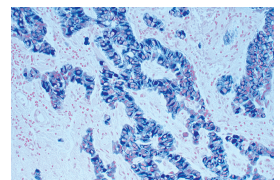
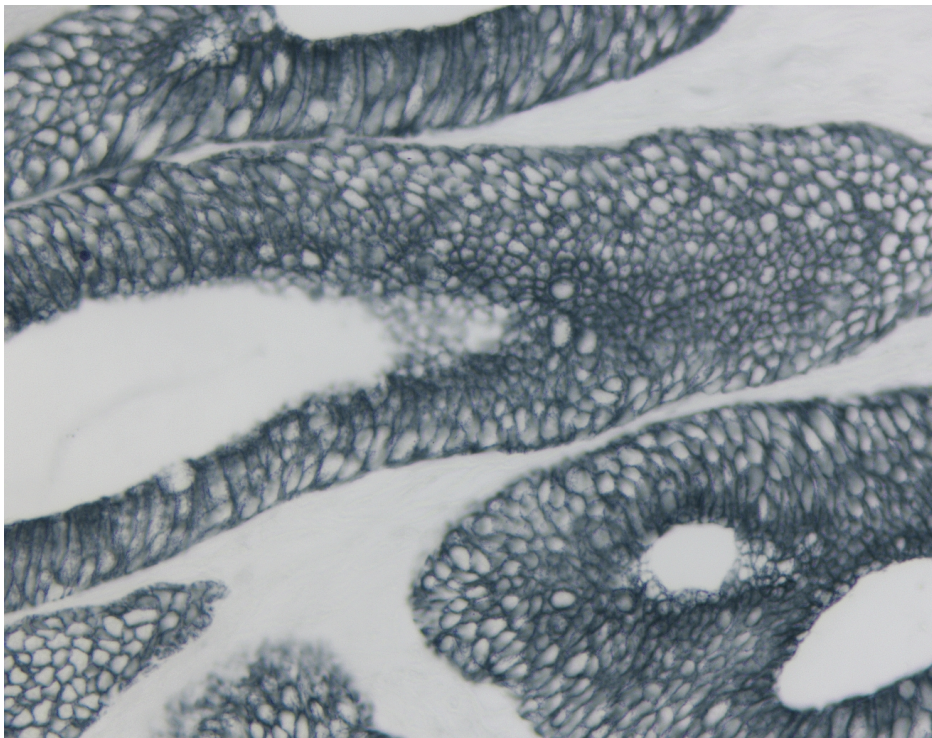


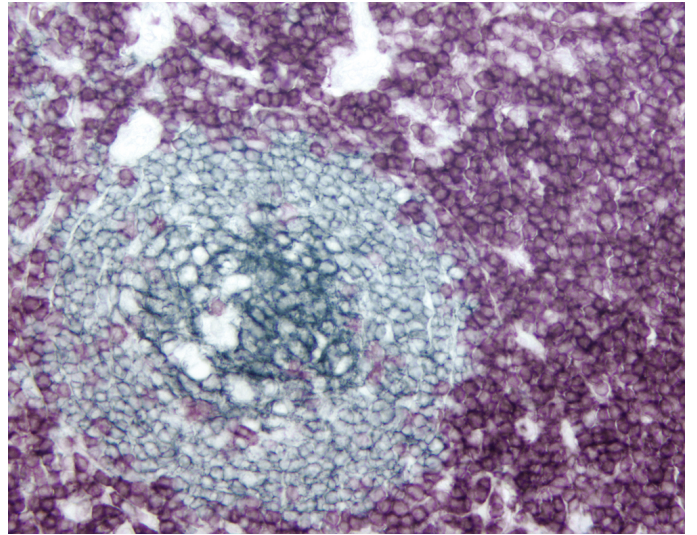
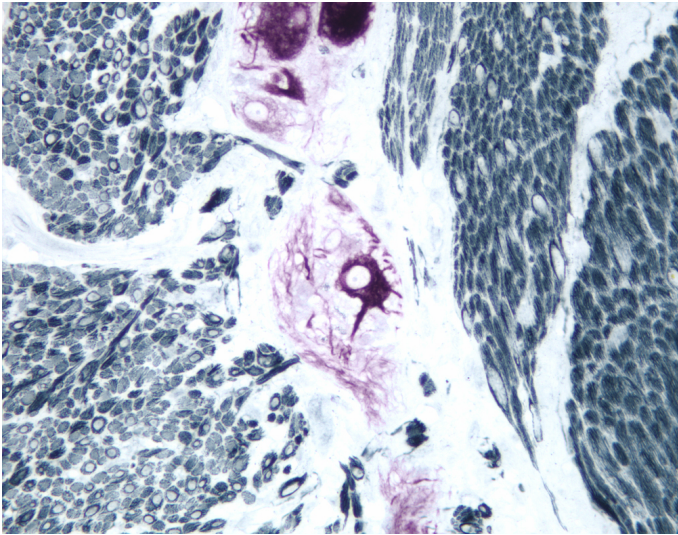


Vector® SG Substrate Kit, Peroxidase (HRP)

SK-4700

Product Images





Short Description

The Vector SG HRP substrate produces a blue-gray reaction product that can be used as a single label or as a second color for multiple antigen labeling.

Features:

- Greater sensitivity than conventional substrates
- Consistent and reliable
- Ideal for IHC/ICC/ISH and blots
- Permanent mounting
- No wait times for mixing and dissolving powders or tablets

Additional Information

Unit Size	1 kit
Applications	Immunohistochemistry / Immunocytochemistry, In situ hybridization, Blotting Applications, Elispot
Detection Enzyme(s)	Peroxidase
Substrate Reaction Product	Precipitating Reaction Product
Chromogen Color	Blue-Gray
Microscopy	Brightfield, Darkfield, Electron, Spectral Imaging
Additional Substrate Properties	Contrast with Tissue Pigments, Multiple Labeling
Mounting	Aqueous (Hardening), Non-Aqueous (Permanent)

