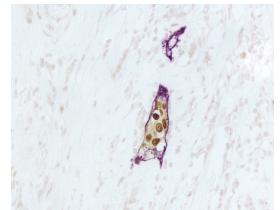
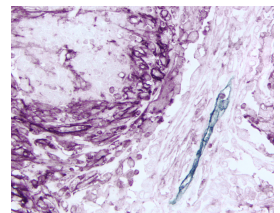
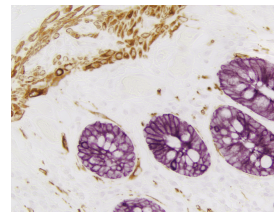
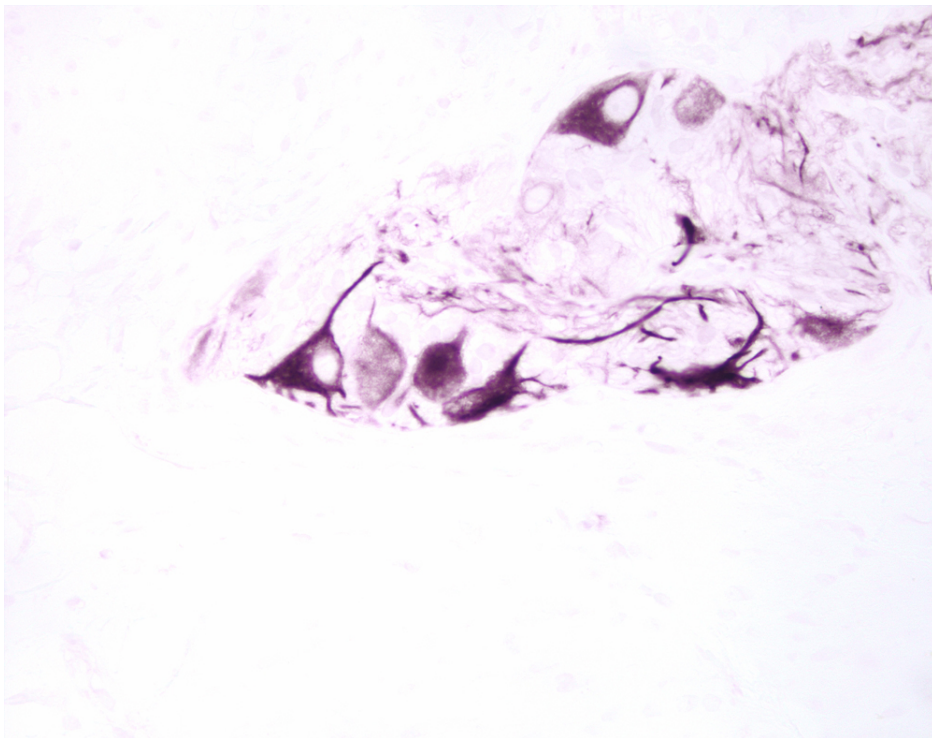


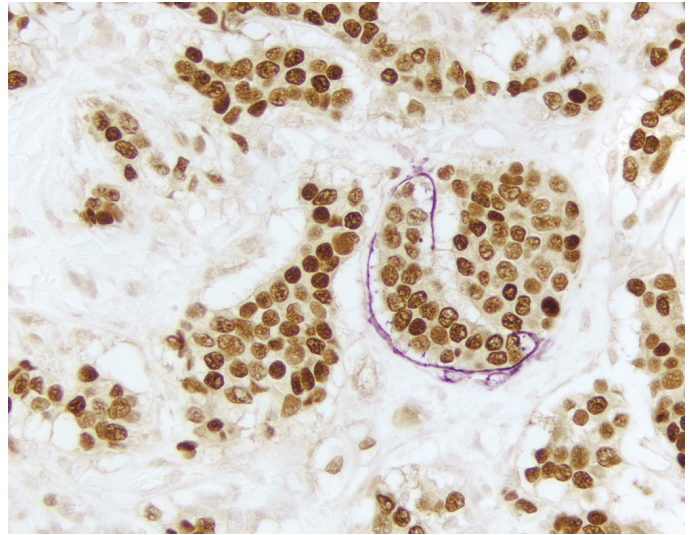
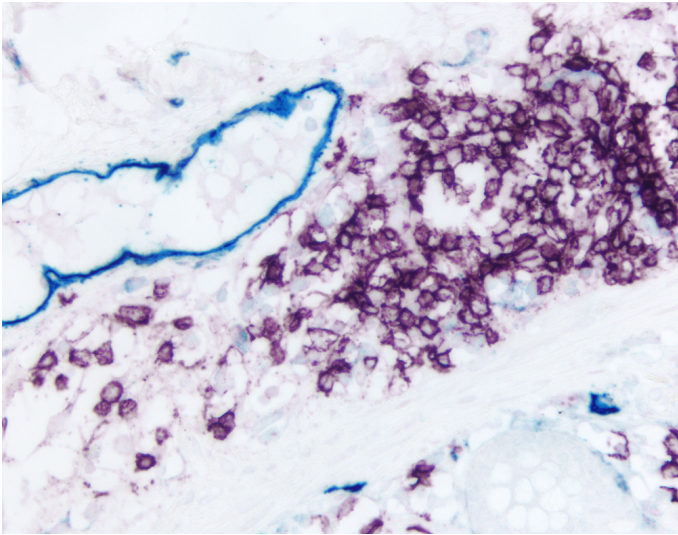


Vector® VIP Substrate Kit, Peroxidase (HRP)

SK-4600

[Product Images](#)





Short Description

The Vector VIP HRP substrate produces an intense, violet (purple) reaction product, and can be used as an alternative to DAB 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
- Stock solutions supplied in convenient dropper bottles promoting ease of handling
- No wait times for mixing and dissolving powders or tablets
- Sufficient reagents to produce 300 ml of working solution

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	Purple
Microscopy	Brightfield, Darkfield, Electron, Spectral Imaging
Additional Substrate Properties	Contrast with Tissue Pigments, Multiple Labeling
Mounting	Non-Aqueous (Permanent)

