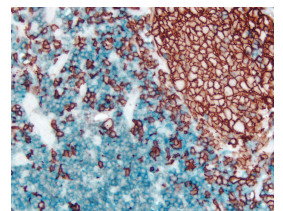
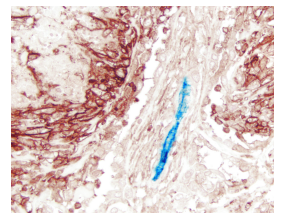
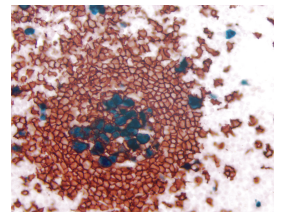
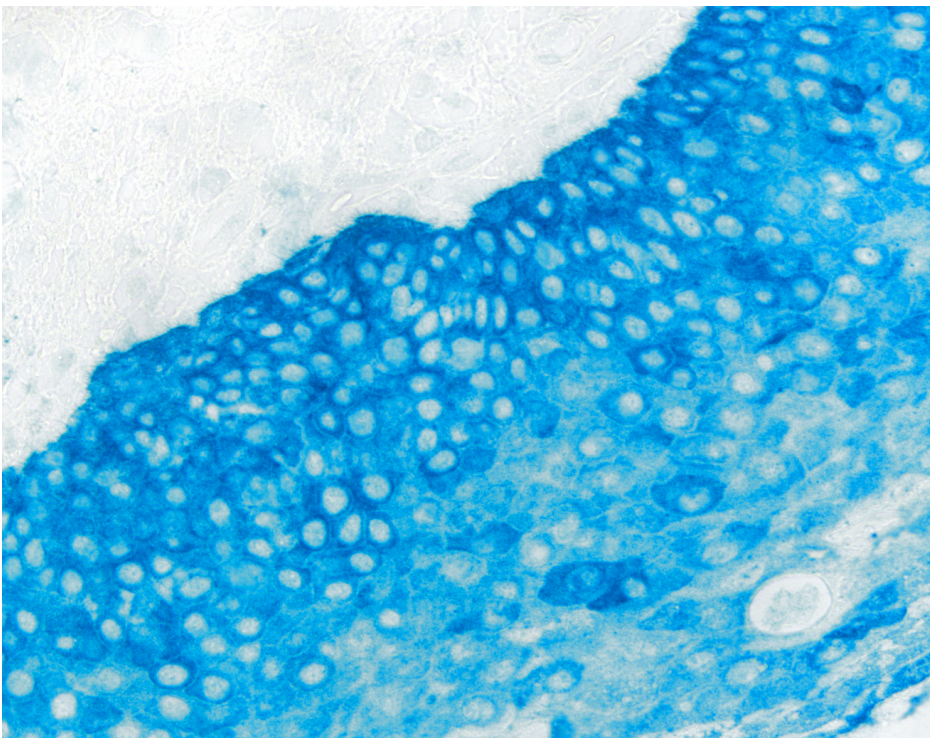


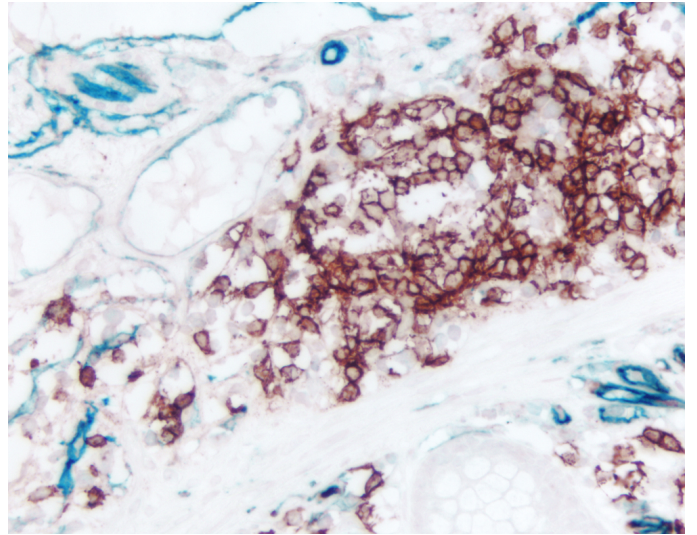
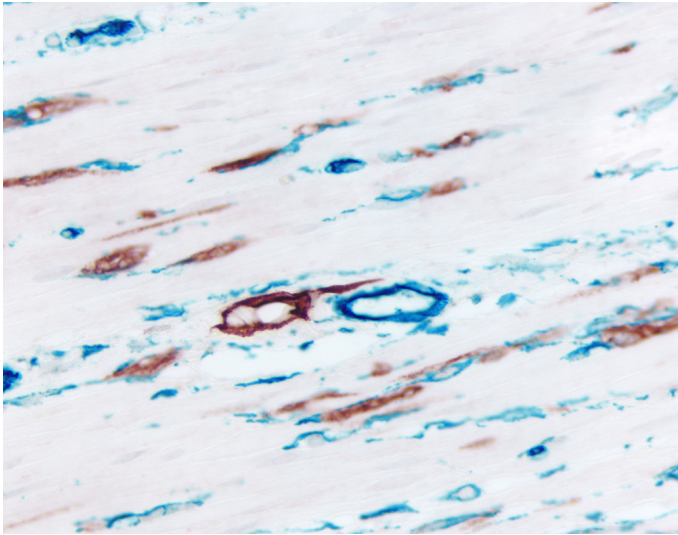


Vector® Blue Substrate Kit, Alkaline Phosphatase (AP)

SK-5300

Product Images





Short Description

The Vector Blue AP Substrate Kit produces a blue, heat stable reaction product. This chromogen can be used singly or in combination with other alkaline phosphatase or peroxidase substrates for multiple label applications.

Features:

- Greater sensitivity than conventional substrates
- Consistent and reliable
- Ideal for IHC/ICC/ISH and blots
- Fluorescent
- Heat Stable
- 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 200 ml of working solution
- Excitation: broad, peak at ~500 nm*
- Emission: broad, peak at ~680 nm*

*Histochem J. (1998) 8:577

Additional Information

Unit Size	1 kit
Applications	Immunohistochemistry / Immunocytochemistry, Immunofluorescence, In situ hybridization, Blotting Applications, Elispot
Detection Enzyme(s)	Alkaline Phosphatase
Substrate Reaction Product	Precipitating Reaction Product
Chromogen Color	Blue
Color of Fluorescence	Far Red
Microscopy	Brightfield, Fluorescence, Spectral Imaging
Additional Substrate Properties	Contrast with Tissue Pigments, Fluorescent, Heat Resistant, Multiple Labeling
Mounting	Aqueous (Hardening), Aqueous (Non-Hardening), Non-Aqueous (Permanent)
Safety Title	Reproductive Harm

