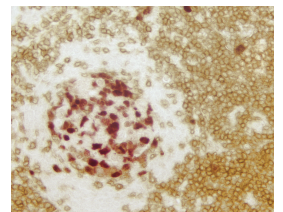
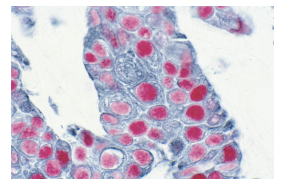
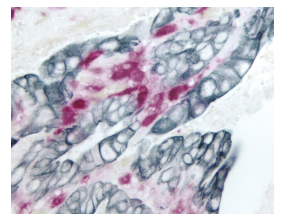
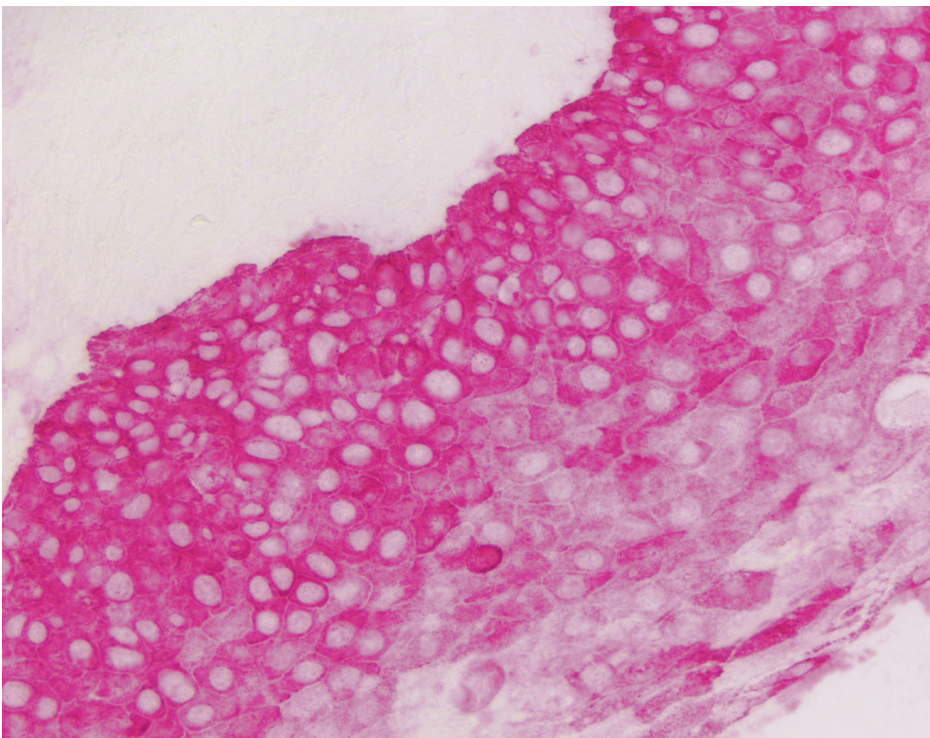


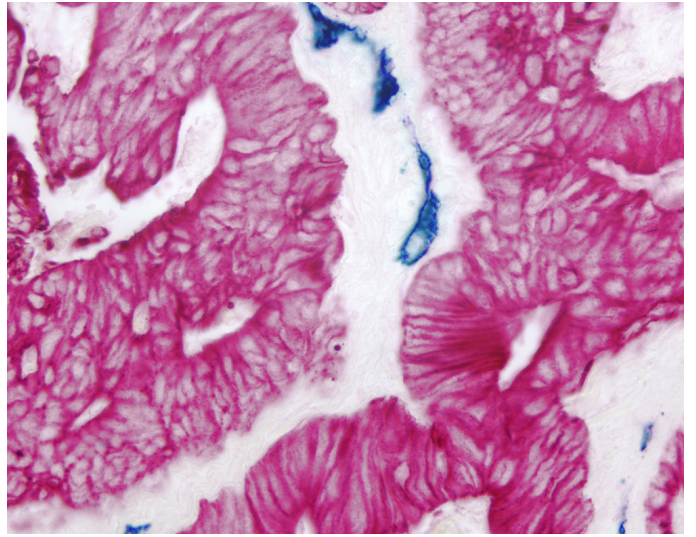
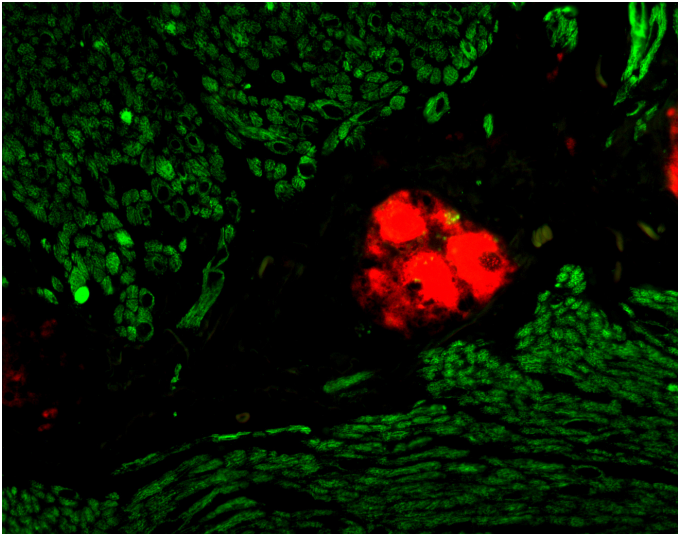


Vector® Red Substrate Kit, Alkaline Phosphatase (AP)

SK-5100

[Product Images](#)





Short Description

Vector® Red AP Substrate Kit produces a magenta heat stable reaction product that is also fluorescent. Sections stained with the Vector® Red substrate can be dehydrated, cleared, and permanently mounted.

Features:

- Greater sensitivity than conventional substrates
- Consistent and reliable
- Ideal for IHC/ICC/ISH and blots
- Ideal for single and multiple labeling
- Stock solutions supplied in convenient dropper bottles promoting ease of handling
- No wait times for mixing and dissolving powders or tablets
- One year expiry date
- Sufficient reagents to produce 200 ml of working solution
- Excitation: broad excitation spectrum, approx. 365 nm to 560 nm
- Emission: Above 560 nm
- Color: Red

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	Magenta
Color of Fluorescence	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)

