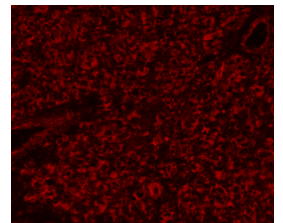
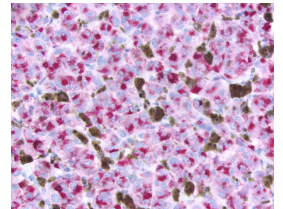
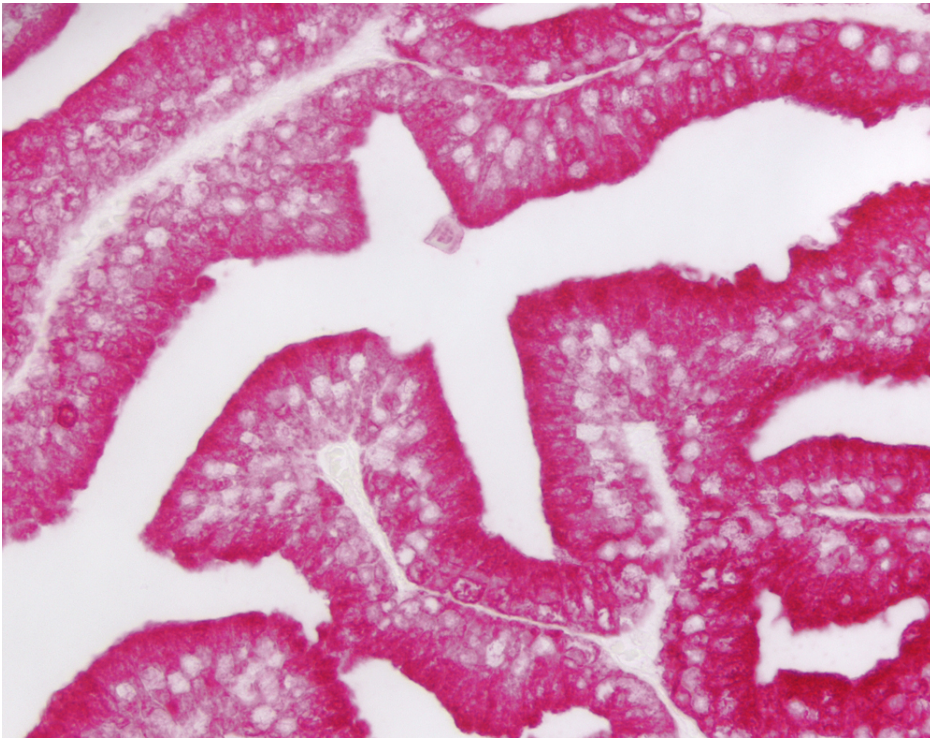


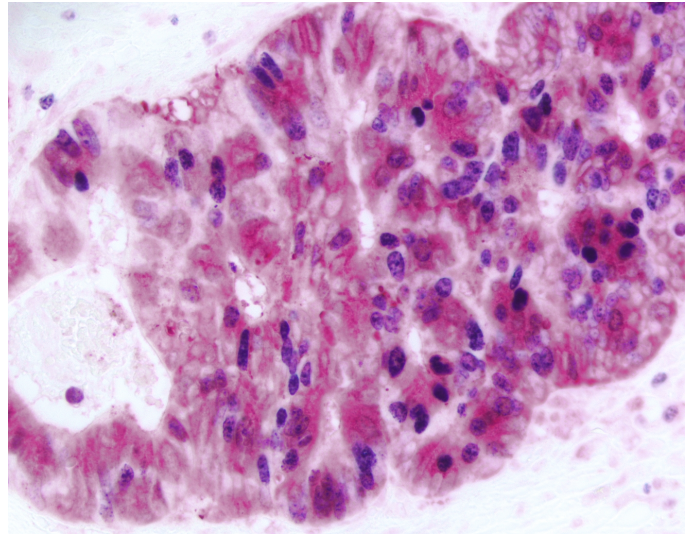
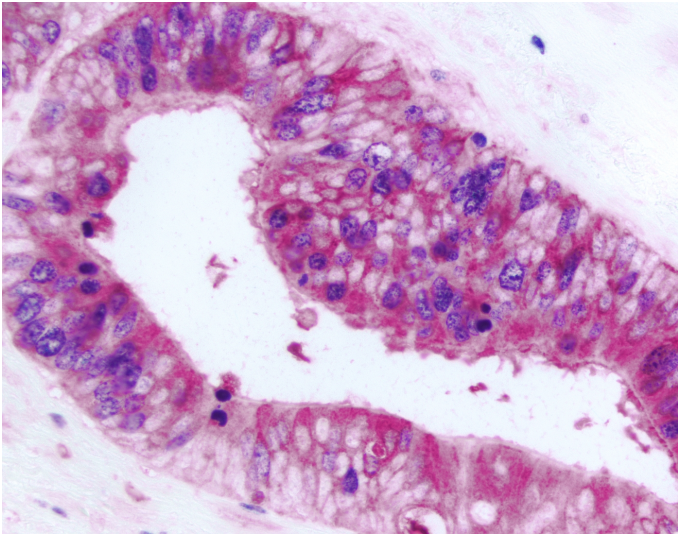


ImmPACT® Vector® Red Substrate Kit, Alkaline Phosphatase (AP)

SK-5105

Product Images





Short Description

ImmPACT Vector Red AP Substrate Kit produces a magenta reaction product that is crisper and is 2-4 times more sensitive than the original Vector® Red substrate kit (SK-5100). It can be used singly or in combination with other alkaline phosphatase or peroxidase substrates for multiple label applications.

Features:

- Ideal for IHC/ICC/ISH and blots
- Fluorescent
- Heat Stable
- Permanent mounting
- Supplied as 120 ml of ready-to-use diluent and stock solutions in dropper bottles
- 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)

