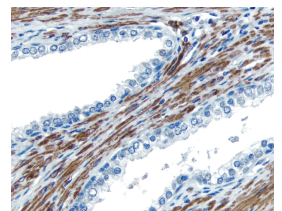
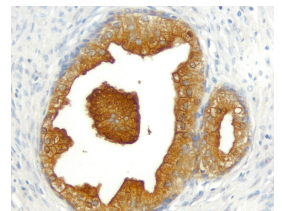
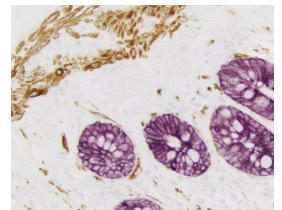
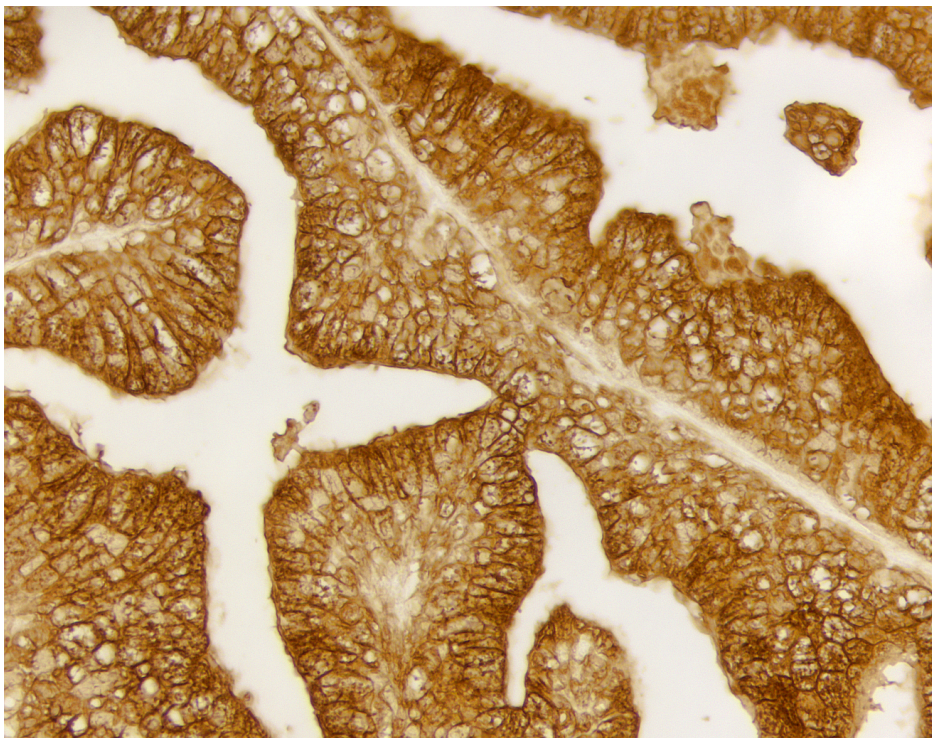


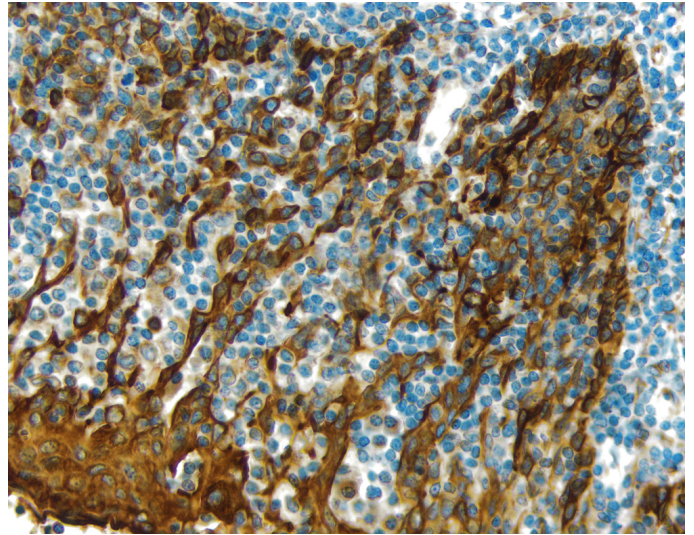
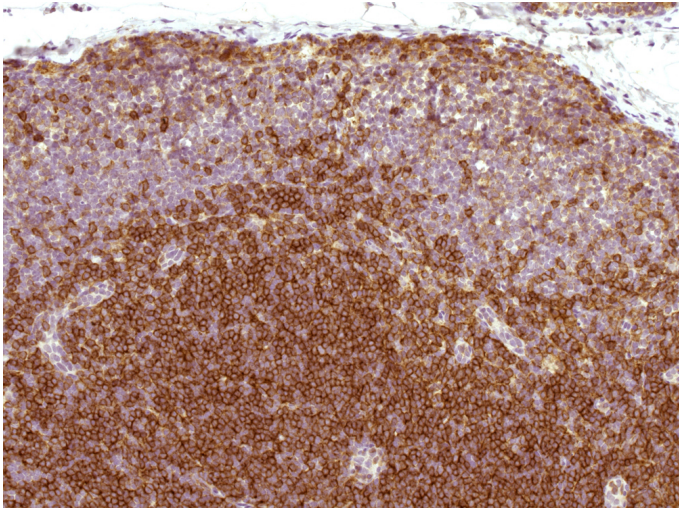


ImmPACT® DAB Substrate Kit, Peroxidase (HRP)

SK-4105

Product Images





Short Description

The ImmPACT DAB Peroxidase (HRP) Substrate reacts with HRP to yield an insoluble brown-colored product. It is 2-4 fold more sensitivity than the original DAB substrate kit.

Features:

- Optimized for IHC/ICC/ISH and blots
- Heat Stable
- Permanent mounting
- Ideal for single and multiple labeling
- Working solutions stable for up to 14 days at 4 °C
- Supplied as 120 ml of diluent and concentrated DAB stock 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	Brown
Microscopy	Brightfield, Darkfield, Electron, Spectral Imaging
Additional Substrate Properties	Heat Resistant, Multiple Labeling
Mounting	Aqueous (Hardening), Aqueous (Non-Hardening), Non-Aqueous (Permanent)

