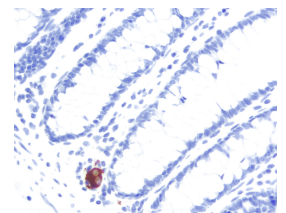
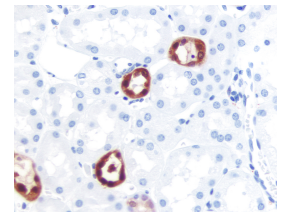
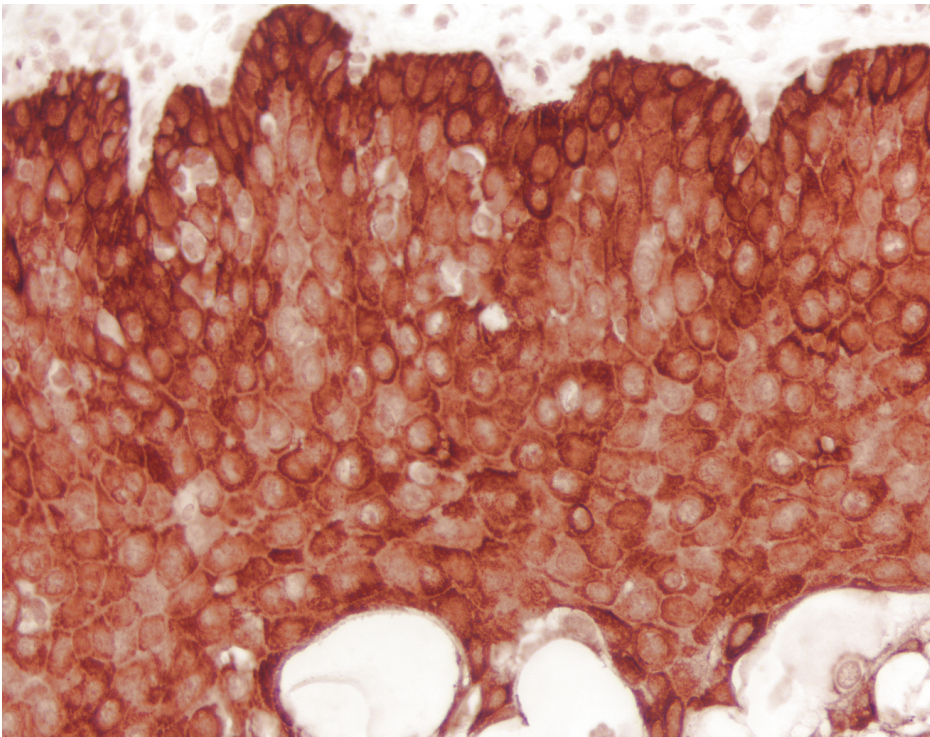




# ImmPACT NovaRED® Substrate Kit, Peroxidase (HRP)

## SK-4805

[Product Images](#)



## Short Description

---

The ImmPACT® NovaRED HRP substrate produces a red reaction products. It produces 2-4 fold greater sensitivity than the original substrate kit (SK-4800).

### Features:

- Optimized for IHC/ICC/ISH and blots
- 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 stock solutions in dropper bottles

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

