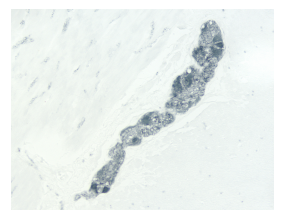
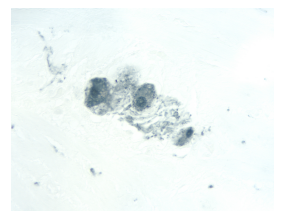
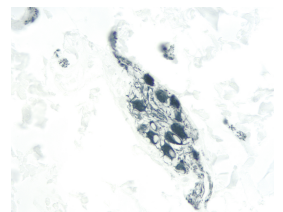
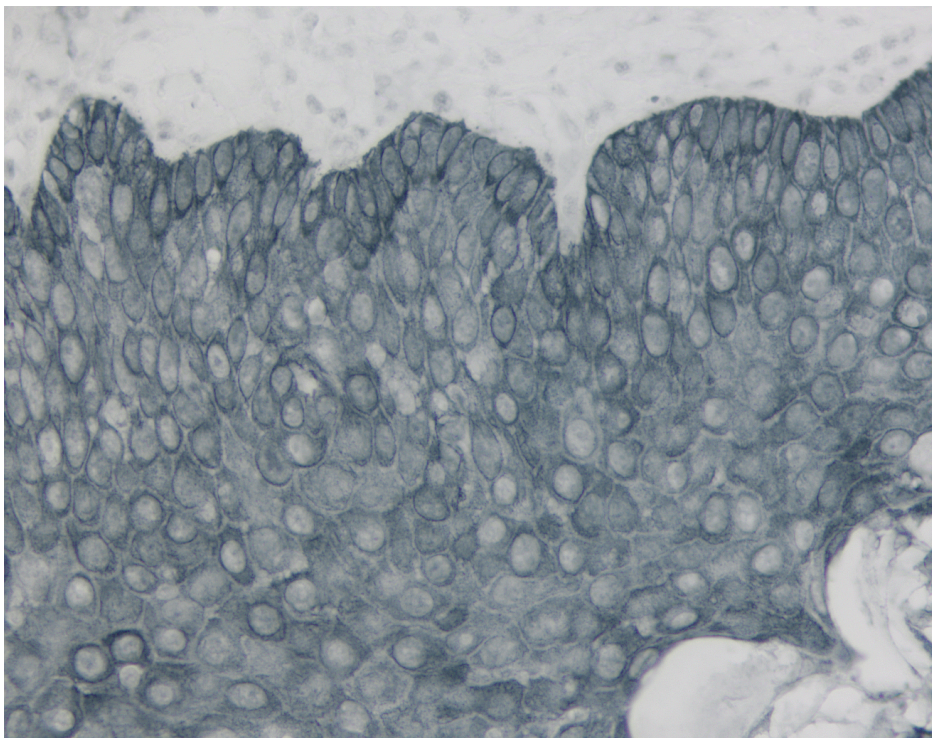




ImmPACT® SG Substrate Kit, Peroxidase (HRP)

SK-4705

[Product Images](#)



Short Description

The ImmPACT SG HRP substrate produces a blue-gray reaction product that can be used as a single label or as a second color for multiple antigen labeling. It is 2-4 fold greater sensitivity than original substrate kit.

Features:

- Optimized for IHC/ICC/ISH and blots
- Permanent mounting
- 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	Blue-Gray
Microscopy	Brightfield, Darkfield, Electron, Spectral Imaging
Additional Substrate Properties	Contrast with Tissue Pigments, Multiple Labeling
Mounting	Aqueous (Hardening), Non-Aqueous (Permanent)

