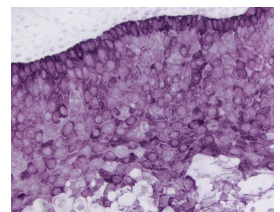
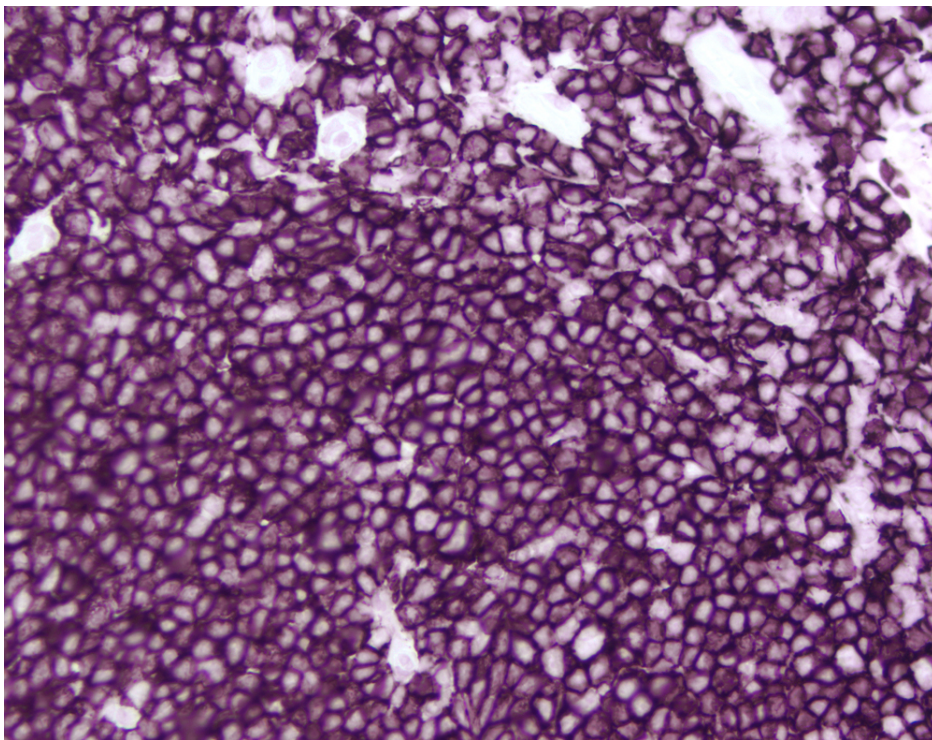




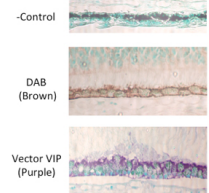
# ImmPACT® VIP Substrate Kit, Peroxidase (HRP)

## SK-4605

[Product Images](#)



EMMPRIN (CD147) Expression in Human Retina



## Short Description

The ImmPACT VIP HRP substrate produces an intense, violet (purple) reaction product, and can be used as an alternative to DAB or as a second color for multiple antigen labeling. This product generates an intense reaction product that is 2-4 times more sensitive than the Vector VIP substrate kit (SK-4600).

### Features:

- 2-4 fold greater sensitivity than original substrate kit
- 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	Purple
Microscopy	Brightfield, Darkfield, Electron, Spectral Imaging
Additional Substrate Properties	Contrast with Tissue Pigments, Multiple Labeling
Mounting	Non-Aqueous (Permanent)

