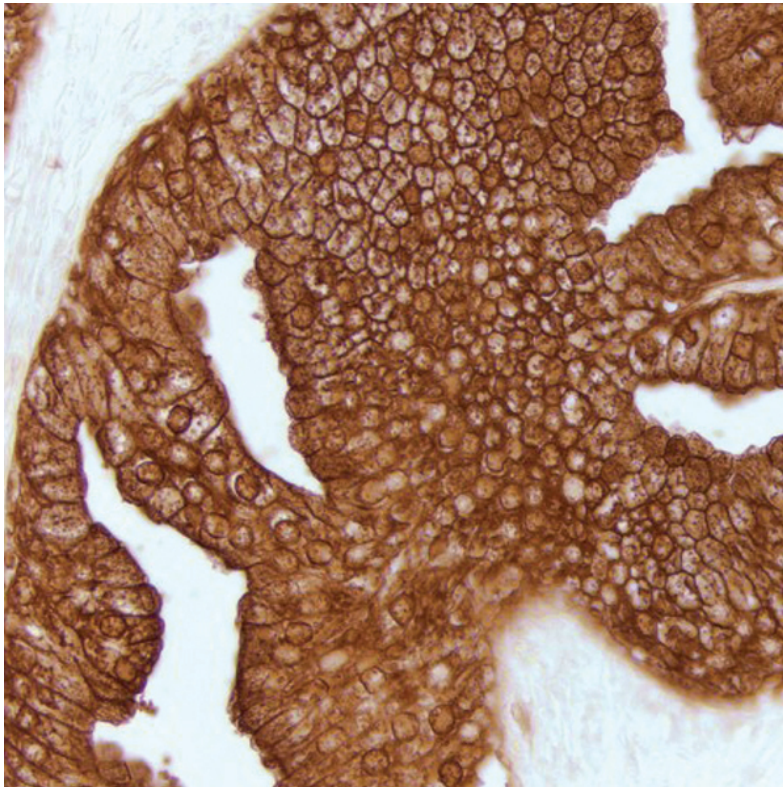




ImmPACT® DAB EqV Substrate Kit, Peroxidase (HRP)

SK-4103

[Product Images](#)



Short Description

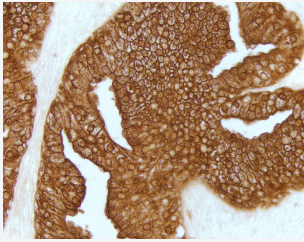
ImmPACT DAB EqV consists of equal volumes of Reagent 1 and Reagent 2, and when mixed together form a working solution. ImmPACT DAB EqV produces a dark brown reaction product and is 3 to 5 times more sensitive than the original DAB substrate.

Features:

- Optimized for IHC/ICC/ISH
- Heat Stable
- Suitable for permanent mounting

Additional Information

Unit Size	100 ml, 400 ml
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)

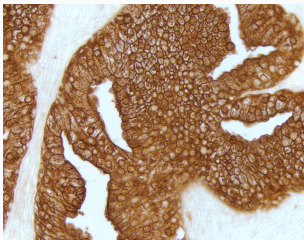


ImmPACT® DAB EqV Substrate Kit, Peroxidase (HRP)

ImmPACT DAB EqV (equal volume) consists of 50 ml each of Reagent 1 and Reagent 2 which when mixed together in equal volumes form the working solution. ImmPACT DAB EqV produces a dark brown reaction product that is 3 to 5 times more sensitive than the original DAB substrate.

Features:

- Optimized for IHC/ICC/ISH
- Heat Stable
- Permanent mounting



ImmPACT® DAB EqV Substrate Kit, Peroxidase (HRP)

ImmPACT DAB EqV (equal volume) consists of 200mL each of Reagent 1 and Reagent 2 which when mixed together in equal volumes form the working solution. ImmPACT DAB EqV produces a dark brown reaction product that is 3 to 5 times more sensitive than the original DAB substrate.

Features:

- Optimized for IHC/ICC/ISH
- Heat Stable
- Suitable for permanent mounting

