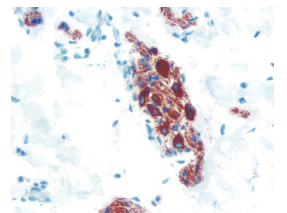
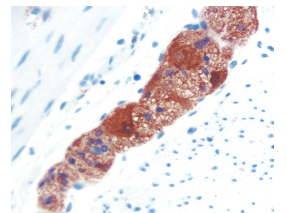
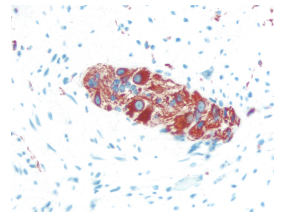
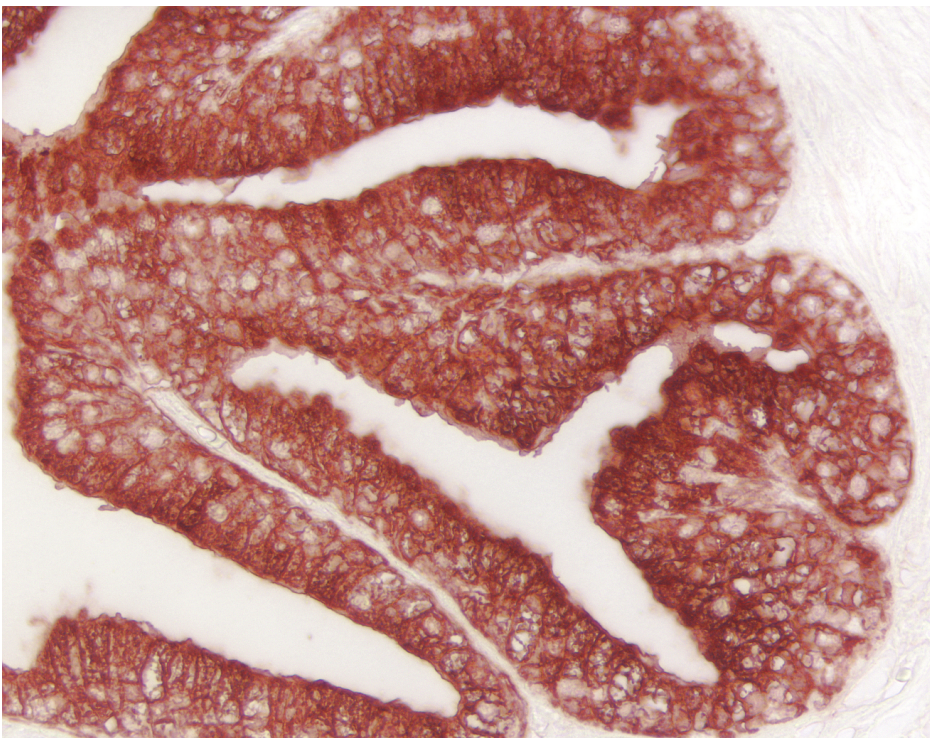




ImmPACT® AEC Substrate Kit, Peroxidase (HRP)

SK-4205

[Product Images](#)



Short Description

ImmPACT AEC (3-amino-9-ethylcarbazole) HRP Substrate produces a red reaction product that can be used as a single label or as a second color for multiple antigen labeling. Stained slides must be aqueously mounted because these reaction products are soluble in organic solvents. It generates a crisper reaction product that is 3-5 times more sensitive than conventional AEC substrates such as AEC Substrate Kit (SK-4200).

Features:

- Optimized for IHC/ICC/ISH and blots
- Working solution 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, Spectral Imaging
Additional Substrate Properties	Contrast with Tissue Pigments, Multiple Labeling
Mounting	Aqueous (Hardening), Aqueous (Non-Hardening)
Safety Title	Cancer

