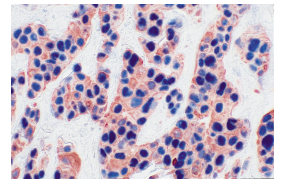
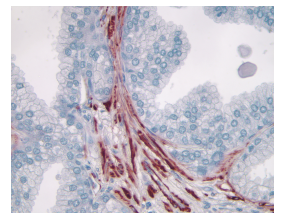
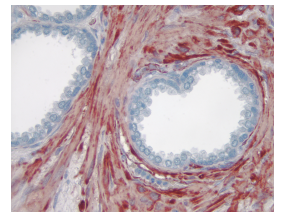
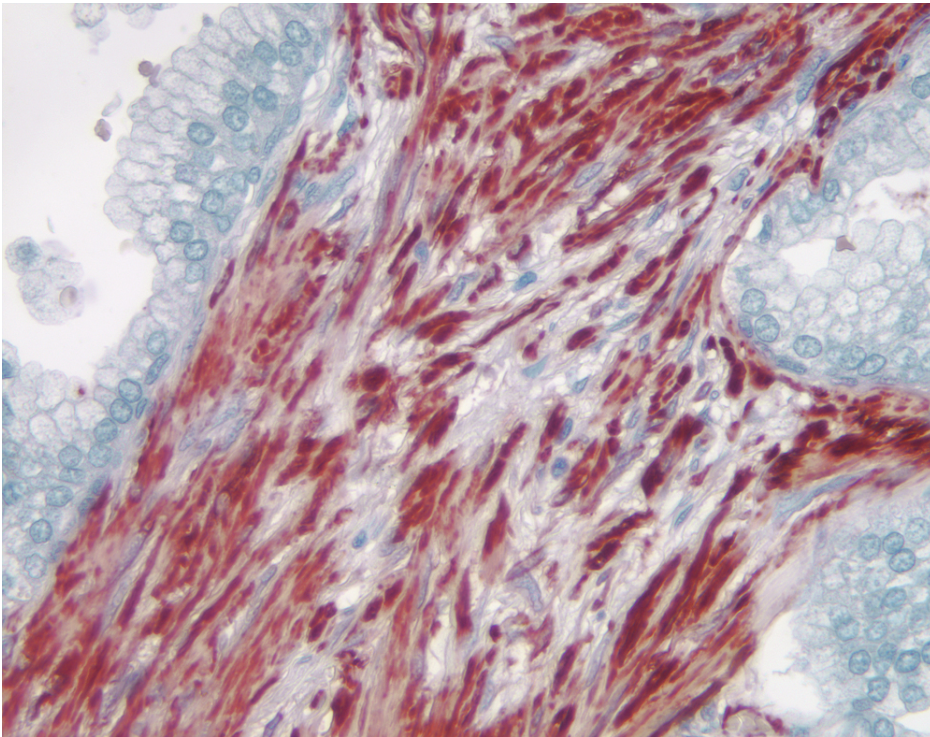




AEC Substrate Kit, Peroxidase (HRP), (3-amino-9-ethylcarbazole)

SK-4200

Product Images



Short Description

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.

Features:

- Greater sensitivity than conventional substrates
- Consistent and reliable
- Suitable for IHC/ICC and blots
- Stock solutions supplied in convenient dropper bottles promoting ease of handling
- No wait times for mixing and dissolving powders or tablets
- Sufficient reagents to produce 300 ml of working solution

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

