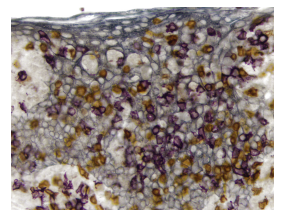
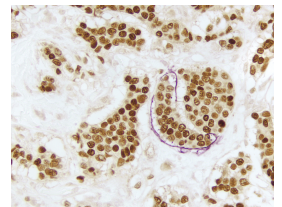
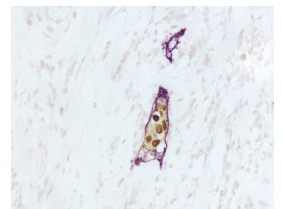


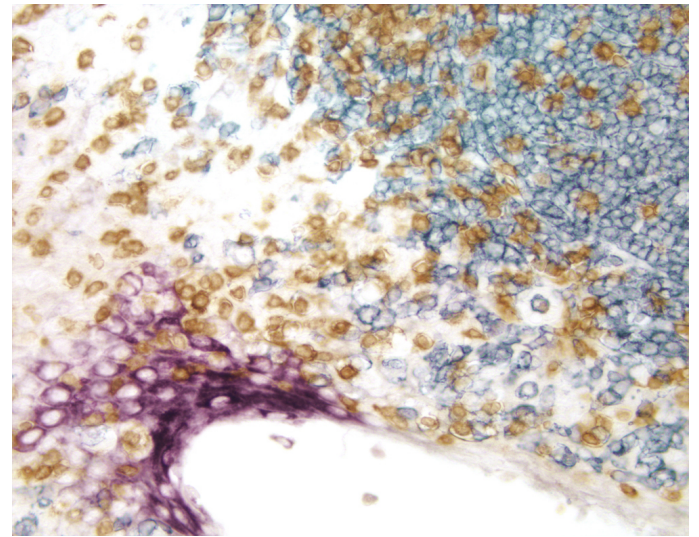
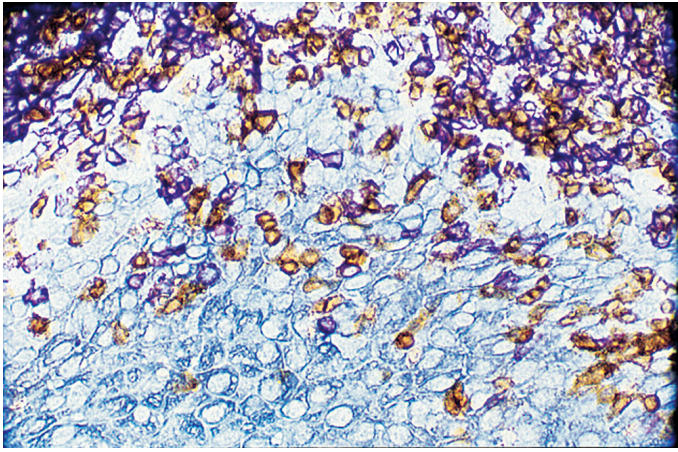


DAB Substrate Kit, Peroxidase (HRP), with Nickel, (3,3'-diaminobenzidine)

SK-4100

[Product Images](#)





Short Description

This kit contains all of the reagents necessary to prepare either a DAB or a DAB/nickel substrate working solution. Vector DAB Substrate (3,3'-diaminobenzidine) produces a brown reaction product in the presence of peroxidase (HRP) enzyme. This kit also includes nickel chloride, which, when added to the substrate working solution, results in a gray-black reaction product. Vector DAB Substrate is suitable for darkfield and electron microscopy (EM).

Features of this kit:

- Ideal for IHC/ICC/ISH and blots
- Heat Stable
- Permanent mounting
- Ideal for single and multiple labeling
- 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	Brown, Gray-Black
Microscopy	Brightfield, Darkfield, Electron, Spectral Imaging
Additional Substrate Properties	Heat Resistant, Multiple Labeling
Mounting	Aqueous (Hardening), Non-Aqueous (Permanent)
Safety Title	Cancer

