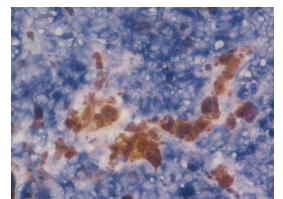
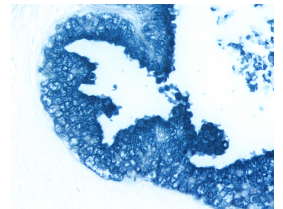
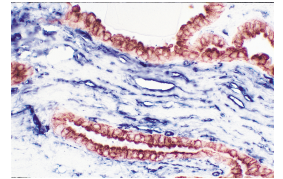
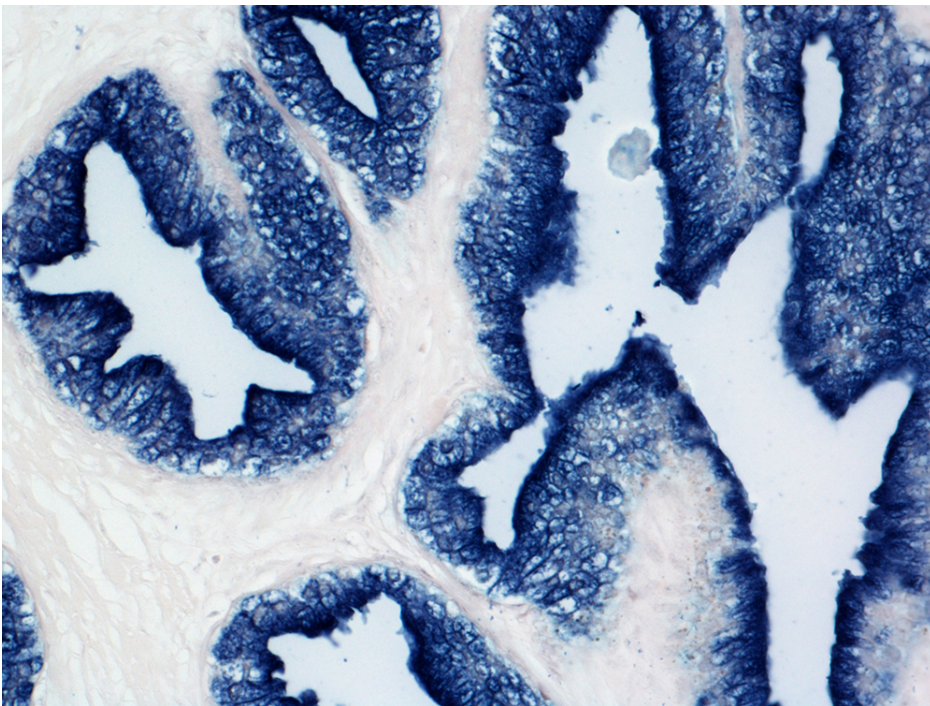
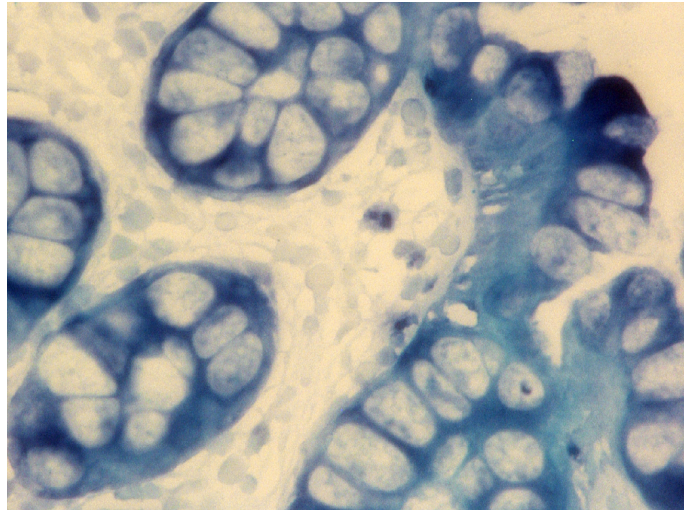
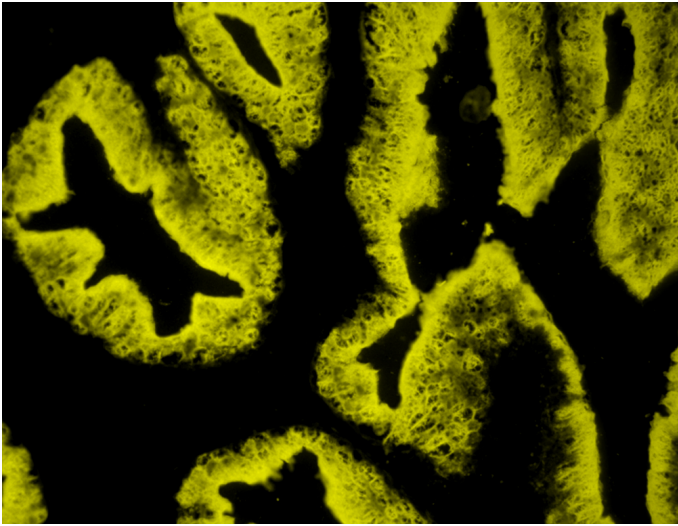


# BCIP/NBT Substrate Kit, Alkaline Phosphatase (AP), (5-bromo-4-chloro-3-indolyl phosphate/nitroblue tetrazolium)

## SK-5400

[Product Images](#)





## Short Description

---

BCIP/NBT AP Substrate Kit forms an indigo, heat stable reaction product. This chromogen can be used singly or in combination with other alkaline phosphatase or peroxidase substrates for multiple label applications.

### Features:

- Greater sensitivity than conventional substrates
- Consistent and reliable
- Ideal for IHC/ICC/ISH and blots
- Fluorescent
- Heat Stable
- Permanent mounting
- No wait times for mixing and dissolving powders or tablets
- Sufficient reagents to produce 200 ml of working solution
- Excitation: 645-685 nm\*
- Emission: peaks at 823 and 855 nm\*

\*Biotechniques (2007) 42:756-759 and BioTechniques (2014) 57:254-256

## Additional Information

---

Unit Size	1 kit
Applications	Immunohistochemistry / Immunocytochemistry, Immunofluorescence, In situ hybridization, Blotting Applications, Elispot
Detection Enzyme(s)	Alkaline Phosphatase
Substrate Reaction Product	Precipitating Reaction Product
Chromogen Color	Indigo
Color of Fluorescence	Far Red
Microscopy	Brightfield, Fluorescence, Spectral Imaging
Additional Substrate Properties	Contrast with Tissue Pigments, Fluorescent, Heat Resistant, Multiple Labeling
Mounting	Aqueous (Hardening), Aqueous (Non-Hardening), Non-Aqueous (Permanent)
Safety Title	Reproductive Harm

