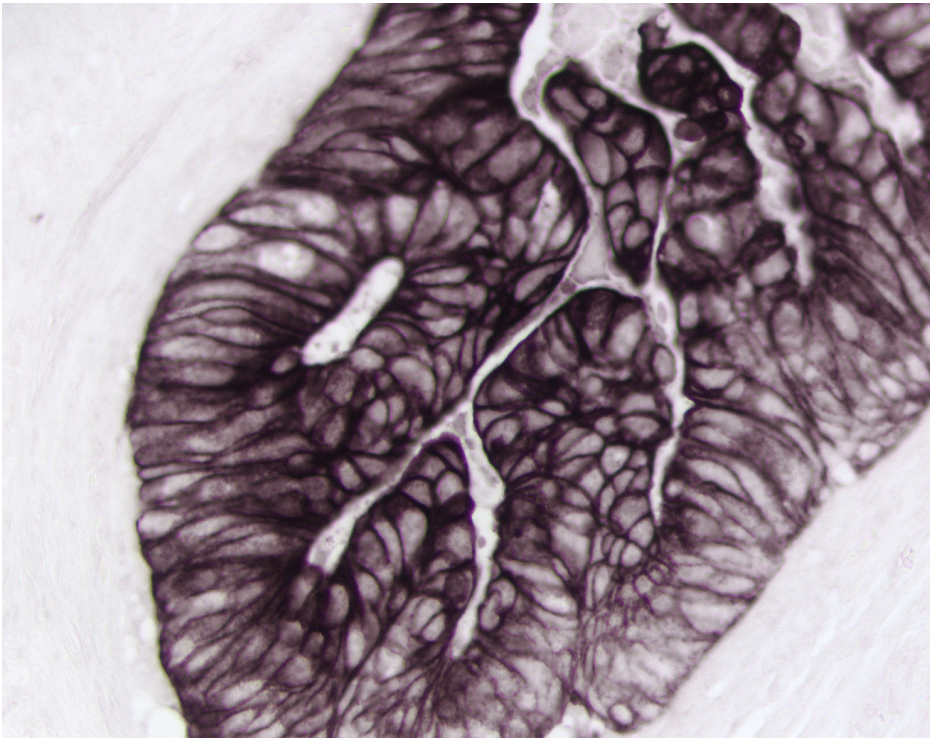




Vector® Black Substrate Kit, Alkaline Phosphatase (AP)

SK-5200

[Product Images](#)



Short Description

Vector Black AP Substrate Kit forms a brown-black precipitate that can be dehydrated, cleared, and permanently mounted. This chromogen can be used in immunohistochemical or blotting applications.

Features:

- Greater sensitivity than conventional substrates
- Consistent and reliable
- 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 200 ml of working solution

Additional Information

Unit Size	1 kit
Applications	Immunohistochemistry / Immunocytochemistry, Blotting Applications, Elispot
Detection Enzyme(s)	Alkaline Phosphatase
Substrate Reaction Product	Precipitating Reaction Product
Chromogen Color	Brown-Black
Microscopy	Brightfield
Mounting	Non-Aqueous (Permanent)
Safety Title	Reproductive Harm

