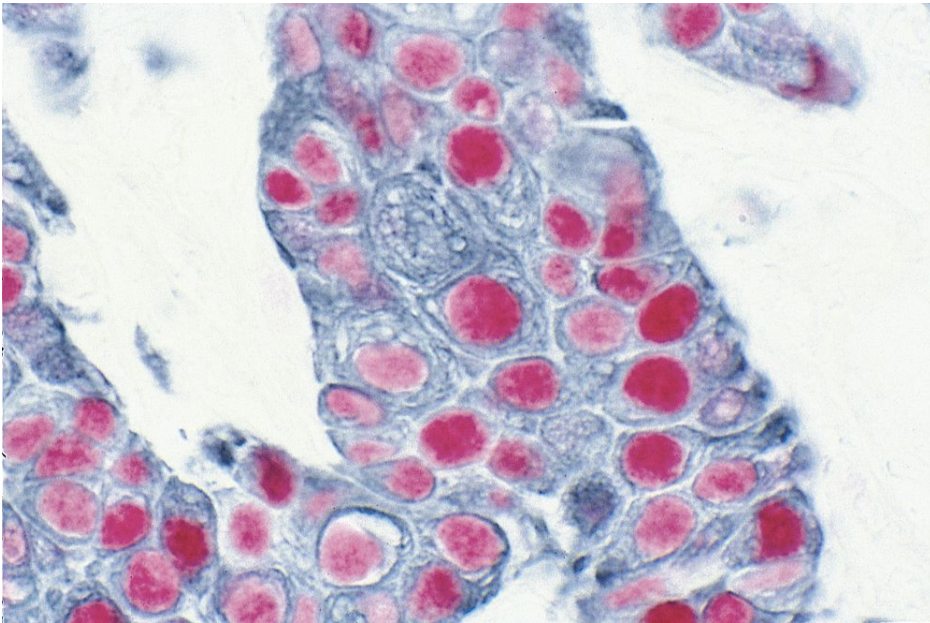




VECTASTAIN® ABC-AP Kit, Alkaline Phosphatase (Standard)

AK-5000

[Product Images](#)



Short Description

The VECTASTAIN ABC Reagent can be used to detect any biotinylated molecule. This is the original avidin/biotin ABC complex formation without a biotinylated secondary antibody.

Features:

- Alkaline phosphatase enzyme detection system
- Avidin/biotin technology
- High sensitivity/low background
- Preferred for tissues that have high endogenous peroxidase activity
- Offers additional substrate choices
- Used as a second label in multiple antigen labeling applications

Product review on [Biocompare.com](https://www.biocompare.com)

Additional Information

Unit Size	1 kit
Applications	Immunohistochemistry / Immunocytochemistry, In situ hybridization, Blotting Applications, Elispot, ELISAs
Technology	ABC (Avidin/Biotin) Systems
Detection Enzyme(s)	Alkaline Phosphatase
Target Species	No antibody included
Format	Concentrate

