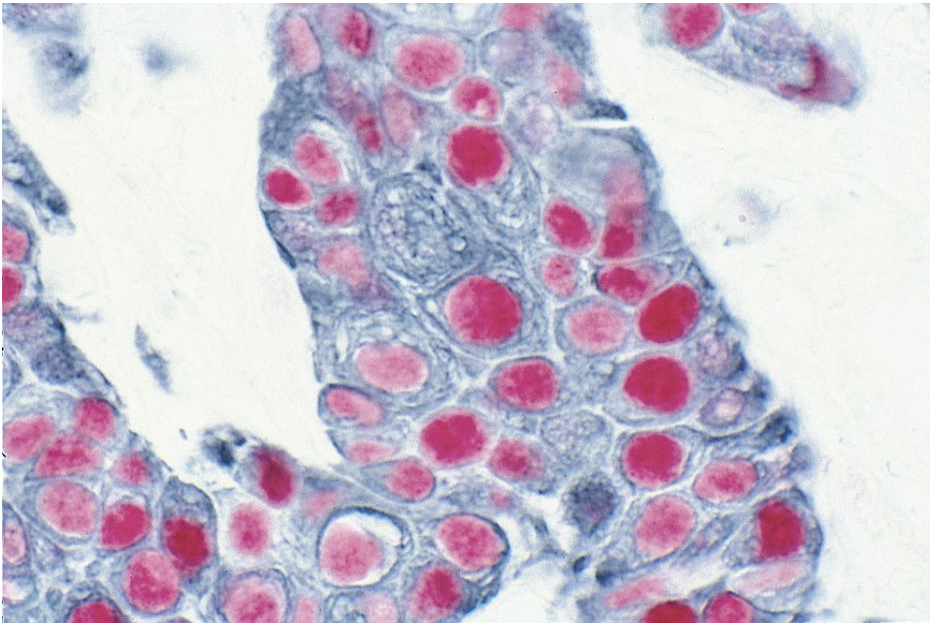




# VECTASTAIN® ABC-AP Kit, Alkaline Phosphatase (Rabbit IgG)

## AK-5001

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 with a biotinylated anti-rabbit IgG 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

## 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	Rabbit
Format	Concentrate

