



# VECTASTAIN® ABC-HRP Kit, Peroxidase (Guinea Pig IgG )

## PK-4007

[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 biotinylated anti-guinea pig IgG secondary antibody.

### Features:

- Good sensitivity
- Low background
- Low cost
- Kit contains sufficient reagents to stain approximately 1000-2000 sections

## Additional Information

---

Unit Size	1 kit
Applications	Immunohistochemistry / Immunocytochemistry, Blotting Applications, Elispot, ELISAs
Technology	ABC (Avidin/Biotin) Systems
Detection Enzyme(s)	Peroxidase
Target Species	Guinea Pig
Format	Concentrate

