



VECTASTAIN® ABC-HRP Kit, Peroxidase (Rabbit IgG)

PK-4001

[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-rabbit IgG secondary antibody.

Features:

- Good sensitivity
- Low background
- 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	Rabbit
Format	Concentrate

