



# VECTASTAIN® ABC-HRP Kit, Peroxidase (Standard)

## PK-4000

[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:

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

## Additional Information

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

