



VECTASTAIN® Elite® ABC-HRP Kit, Peroxidase, R.T.U. (Universal)

PK-7200

[Product Images](#)



Short Description

The VECTASTAIN Elite ABC system is the most sensitive avidin/biotin-based peroxidase system and is approximately 5 times more sensitive than the original VECTASTAIN ABC Kit. Our Universal Biotinylated Anti-Mouse/Rabbit IgG secondary antibody included in this kit is a cocktail of biotinylated anti-mouse IgG and anti-rabbit IgG, designed for use with both rabbit and mouse primary antibodies.

Features:

- **Advanced avidin/biotin technology:** The Elite ABC complex is smaller, very uniform, and highly active, allowing more accessibility for binding to a biotinylated target
- **Cost effective:** Higher sensitivity leads to lower cost per slide
- **Available in ready-to-use formats:** Prediluted, stabilized working solutions of Elite ABC Kit reagents provide the same high sensitivity and low background as the traditional VECTASTAIN Elite ABC Kit reagents.

Contains sufficient reagent to stain approximately 250-500 sections.

Note: This antibody should not be used to stain tissues from rodents or rabbits due to its reactivity with IgG present in the tissues of these species.

Additional Information

Unit Size	1 kit
Applications	Immunohistochemistry / Immunocytochemistry, In situ hybridization, Blotting Applications, Elispot, ELISAs
Technology	ABC Elite (Avidin/Biotin) Systems
Detection Enzyme(s)	Peroxidase
Target Species	Universal (Mouse/Rabbit)
Format	Ready-to-Use

