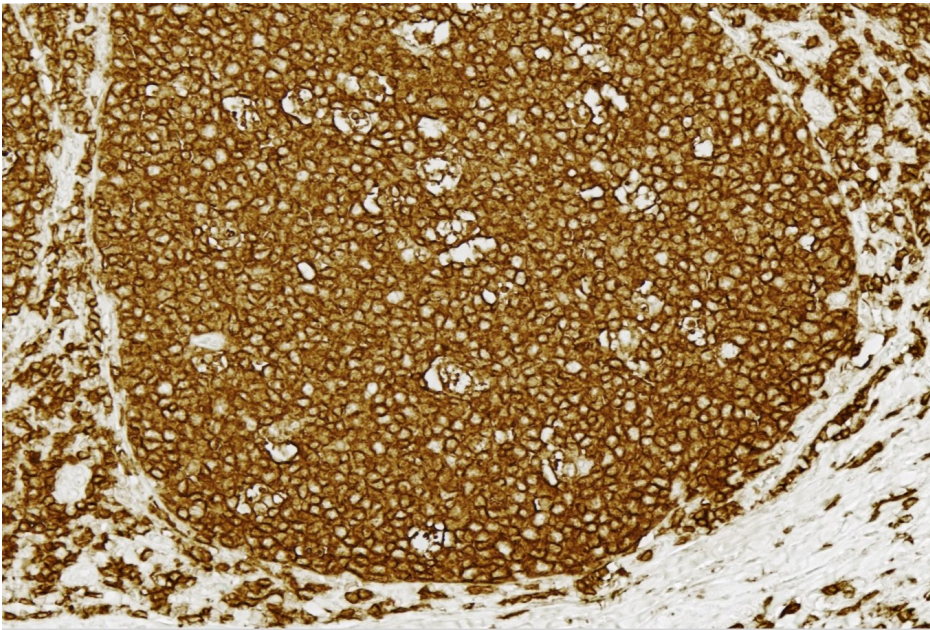




# VECTASTAIN® Elite® ABC Universal PLUS Kit, Peroxidase (Horse Anti-Mouse/Rabbit IgG)

## PK-8200

[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 with the same low background. 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. **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.**

### 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
- **More complete detection systems:** Inclusion of BLOXALL<sup>®</sup> endogenous enzyme quenching solution and ImmPACT<sup>®</sup> DAB EqV substrate in the PLUS kit provides added convenience with optimized, matched reagents in the one product.

## 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

