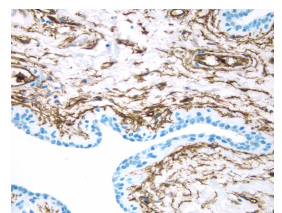
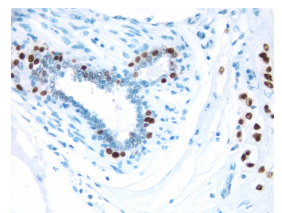
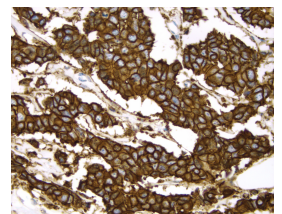
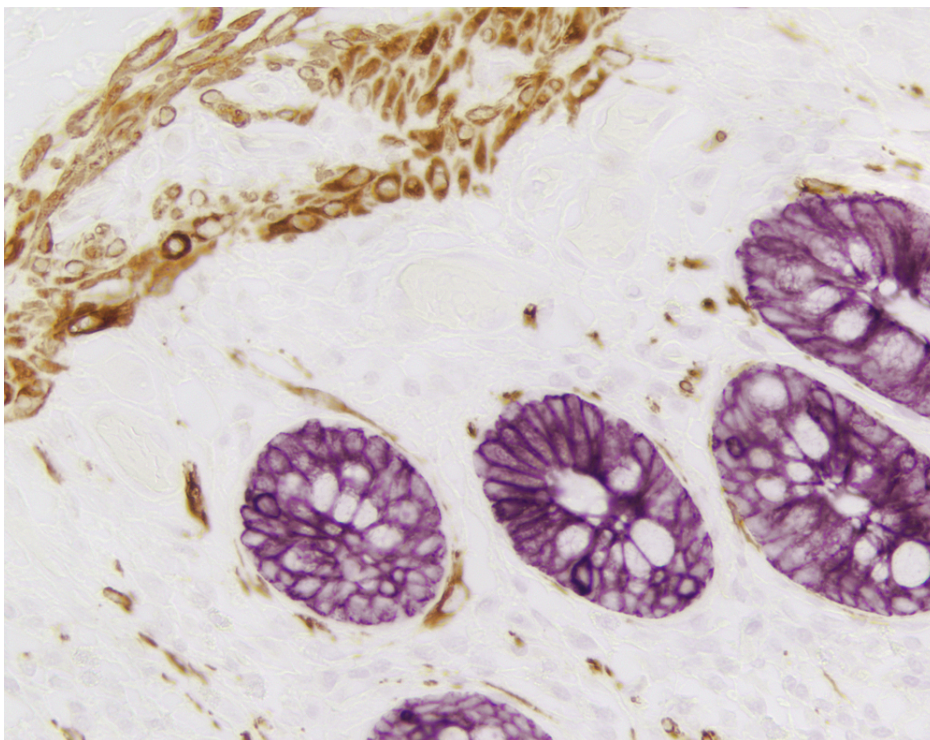




VECTASTAIN® Elite® ABC-HRP Kit, Peroxidase (Mouse IgG)

PK-6102

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. It is available with a biotinylated anti-mouse IgG secondary antibody.

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

Kit contains sufficient reagents to stain approximately 500-1000 sections.

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	Mouse
Format	Concentrate

