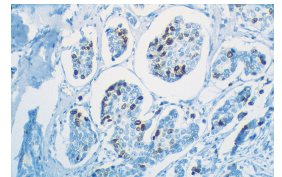
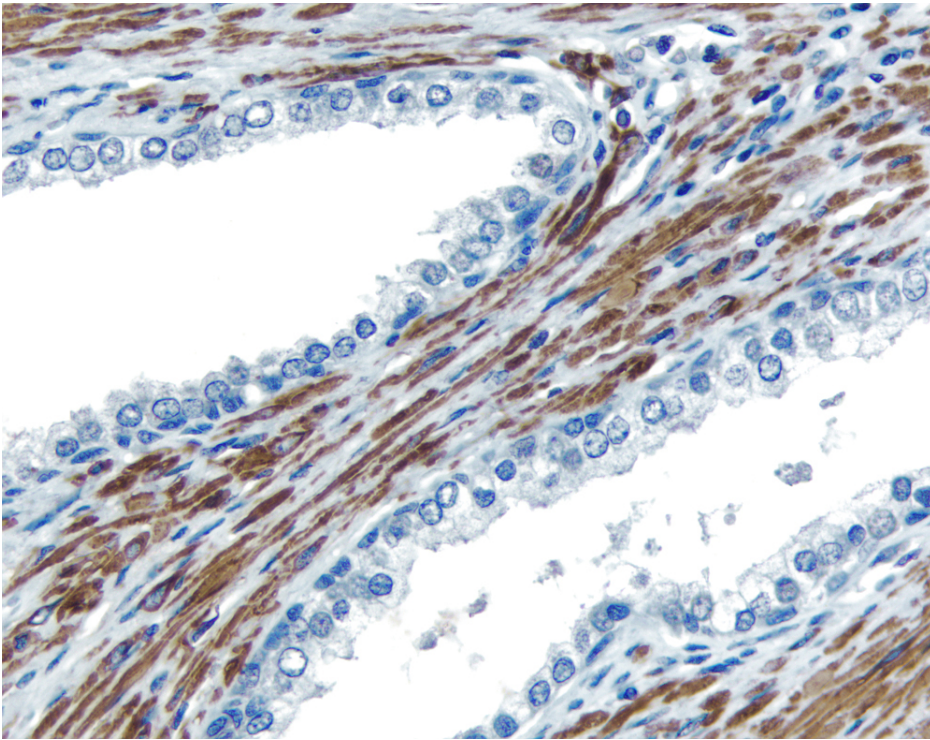




# VECTASTAIN® Universal Quick HRP Kit, Peroxidase (Concentrate)

## PK-8800

[Product Images](#)



## Short Description

The VECTASTAIN Universal Quick Kit is designed for rapid immunohistochemical staining of tissue sections. The kit relies on a proprietary preformed peroxidase-streptavidin complex to achieve outstanding sensitivity with short incubation times.

### Features:

- **Rapid protocol:** Staining in less than 20 minutes following primary antibody incubation. Working solutions can be used immediately after dilution.
- **High sensitivity, low background.**
- **Biotinylated, Universal Pan-Specific secondary** antibody recognizes mouse, rabbit, and goat primary antibodies as well as those from related species such as rat, bovine, and sheep.\*
- **Concentrated format:** Reagents are supplied as liquid concentrates enabling flexibility in dilution and application while maximizing cost-savings per assay.

*\*Do not use the pan-specific secondary antibody to stain rat, mouse or other rodent, rabbit, goat, bovine, or sheep tissue due to potential reactivity with endogenous IgG.*

Sufficient reagents to stain approximately 250-500 sections.

## Additional Information

Unit Size	1 kit
Applications	Immunohistochemistry / Immunocytochemistry
Technology	ABC (Streptavidin/Biotin) Systems
Detection Enzyme(s)	Peroxidase
Target Species	Universal (Mouse/Rabbit/Goat)
Format	Concentrate

