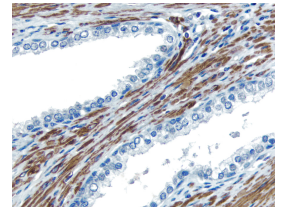
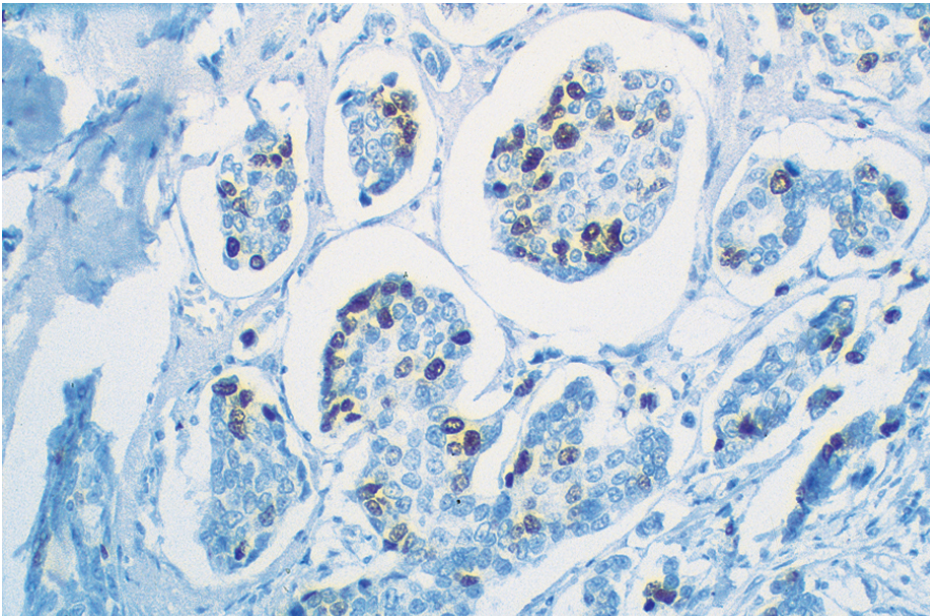




# VECTASTAIN® Universal Quick HRP Kit, Peroxidase, R.T.U.

## PK-7800

[Product Images](#)



## Short Description

---

The VECTASTAIN Universal Quick HRP Kit can detect any biotinylated molecule. Staining can be completed in less than 20 minutes following primary antibody incubation.

### Features:

- **Rapid protocol:** 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.\*
- **Available in Ready-To-Use (R.T.U) format:** Prediluted, stabilized working solutions of VECTASTAIN Universal Quick Kit reagents provide the same high sensitivity and low background as the concentrated VECTASTAIN Universal Kit.

*\*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           | 50 ml                                      |
| Applications        | Immunohistochemistry / Immunocytochemistry |
| Technology          | ABC (Streptavidin/Biotin) Systems          |
| Detection Enzyme(s) | Peroxidase                                 |
| Target Species      | Universal (Mouse/Rabbit/Goat)              |
| Format              | Ready-to-Use                               |

