

VECTASTAIN® Universal Quick Kit

Peroxidase

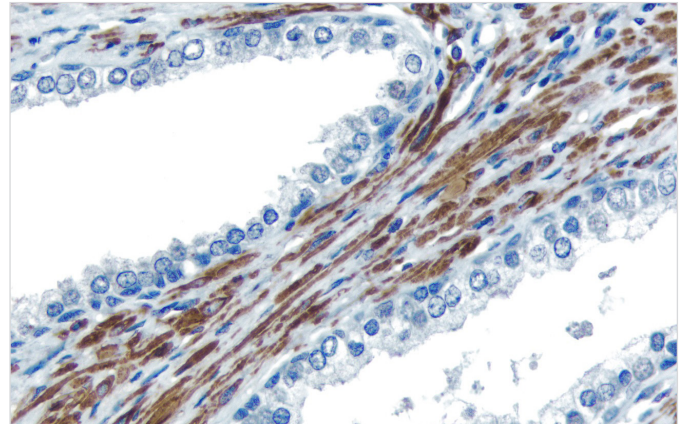
Cat. No.	PK-8800
Storage	Store reagents in original bottles at 2-8 °C. Do not freeze.
Description	Instructions for Immunohistochemical staining.

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. This kit utilizes a biotinylated, universal, pan-specific secondary antibody that recognizes mouse, rat, rabbit, goat, sheep and bovine primary antibodies.

Kit Components

Product Name	Volume
Normal Horse Serum	6 ml
Biotinylated Pan-Specific Universal Antibody (Horse Anti-Mouse/Rabbit/Goat IgG)	2.2 ml
Streptavidin/Peroxidase Complex	1.2 ml

The VECTASTAIN Universal Quick Kit will stain approximately 250-500 sections.



Prostate: Desmin (m) detected with VECTASTAIN Universal Quick Kit, and ImmPACT® DAB (brown). Counterstained with Vector Hematoxylin QS (blue).

Preparation of Working Solutions

Working solutions are prepared in the following manner using PBS (10 mM phosphate, pH 7.5, 0.15M NaCl as the diluent (1 drop ~ 50 µl):

- Blocking solution (Normal Serum): add 1 drop of Normal Horse Serum to 2 ml of PBS.
- Biotinylated Pan-Specific Universal Antibody: add 4 drops of Normal Horse Serum plus 2 drops of Biotinylated Pan-Specific Universal Antibody to 2 ml of PBS.
- Streptavidin/Peroxidase Complex: add 1 drop of Streptavidin/Peroxidase Complex stock solution to 2 ml of PBS.

The working solutions are ready to use immediately after dilution. Unused diluted working solution can be stored under refrigeration for 5 days without appreciable loss in sensitivity.

Staining Procedure

1. For paraffin sections, deparaffinize and hydrate tissue sections through xylenes or other clearing agents and graded alcohol series.

For frozen sections fix with acetone or an appropriate fixative for the antigen under study, if required.

Wash for 5 minutes in tap water.

2. If antigen unmasking is required, perform this procedure using a Vector® Antigen Unmasking Solution, Citrate-based (H-3300) or Tris-based (H-3301).
3. If quenching of endogenous peroxidase activity is required, incubate the slides in BLOXALL® Blocking Solution (SP-6000) for 10 minutes.
4. Wash in buffer for 5 minutes.
5. Incubate sections for 10 minutes with prepared Normal Serum.
6. Tip off excess serum from sections.
7. Incubate sections in primary antibody diluted buffer containing 1.5% blocking serum.
8. Wash in buffer for 5 minutes.
9. Incubate sections with prepared Biotinylated Pan-Specific Universal Antibody for 10 minutes.
10. Wash in buffer for 5 minutes.
11. Incubate sections in prepared Streptavidin/Peroxidase Complex for 5 minutes.
12. Wash in buffer for 5 minutes.
13. Incubate in a peroxidase substrate solution (not included in kit) until desired stain intensity develops.
14. Rinse in tap water.
15. Counterstain (optional), clear and mount.

Detailed product listing, specifications, protocols and additional information is available on our website: vectorlabs.com