

VECTASTAIN® Universal Quick Kit

Peroxidase, R.T.U.

Cat. No.	PK-7800
Storage	Store reagents in original bottles at 2-8 °C. Do not freeze.
Description	Instructions for Immunohistochemical staining.

The VECTASTAIN Universal Quick Kit was developed using the most advanced biotin/streptavidin complex technology. The components in this kit have been designed to provide the high sensitivity of other biotin/streptavidin systems, but with an accelerated staining protocol utilizing a biotinylated secondary antibody that recognizes mouse, rat, rabbit, goat, sheep, and bovine primary antibodies. The VECTASTAIN Universal Quick Kit allows a primary antibody from different species to be used at high dilution with a rapid staining protocol.

The R.T.U. VECTASTAIN Universal Quick Kit contains three reagents in ready-to-use form. No dilution is necessary.

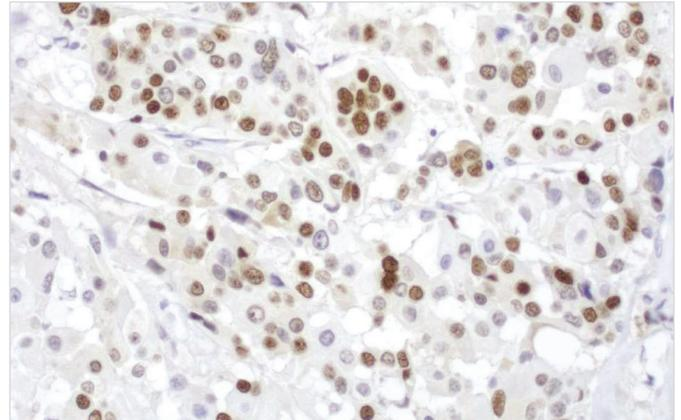
Kit Components

Product Name	Volume
Normal Horse Serum, 2.5%	50 ml
Biotinylated Pan-Specific Universal Antibody (Horse Anti-Mouse/Rabbit/Goat IgG)	50 ml
Streptavidin/Peroxidase Complex	50 ml

The VECTASTAIN Universal Quick Kit will stain approximately 500 sections.

Staining Procedure

- 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.
- If antigen unmasking is required, perform this procedure using a Vector® Antigen Unmasking Solution, Citrate-based (H-3300) or Tris-based (H-3301).
- If quenching of endogenous peroxidase activity is required, incubate the slides in BLOXALL® Blocking Solution (SP-6000) for 10 minutes.



Breast (FFPE): SP4 Cyclin D1 (r) detected with VECTASTAIN Universal Quick Kit, R.T.U. and DAB (brown). Counterstained with Hematoxylin QS (blue).

- Wash in buffer for 5 minutes.
- Incubate sections for 10 minutes with Normal Horse Serum, 2.5%.
- Tip off excess serum from sections.
- Incubate sections in primary antibody diluted buffer containing 1.5% blocking serum.
- Wash in buffer for 5 minutes.
- Incubate sections in Biotinylated Pan-Specific Universal Antibody for 10 minutes.
- Wash in buffer for 5 minutes.
- Incubate sections in Streptavidin/Peroxidase Complex for 5 minutes.
- Wash in buffer for 5 minutes.
- Incubate in a peroxidase substrate solution (not included in kit) until desired stain intensity develops.
- Rinse in tap water.
- Counterstain (optional), clear and mount.

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