

# VECTASTAIN® Elite® ABC-HRP Reagent

Peroxidase, R.T.U.

**Cat. No.** PK-7100

**Storage** Store reagents in original bottles at 2-8 °C. Do not freeze.

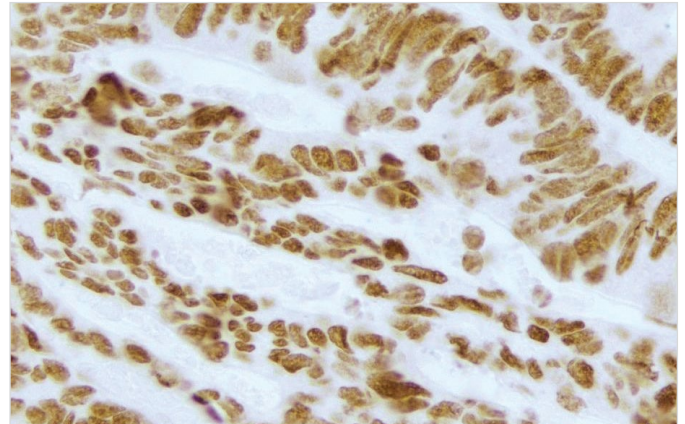
**Description** Instructions for Immunohistochemical staining. The VECTASTAIN Elite ABC System is an enzymatic, avidin/biotin based amplification system that produces crisp, highly sensitive, specific staining with low background. The ready-to-use (R.T.U.) VECTASTAIN Elite ABC Reagent can be used to detect any molecule that is biotinylated.

The R.T.U. reagents are supplied in convenient dropper bottles. No dilution is necessary.

## Kit Components

Product	Volume
VECTASTAIN Elite ABC Reagent, Peroxidase, R.T.U.	50 ml

The VECTASTAIN Elite ABC Reagent will stain approximately 500 sections.



Colon carcinoma (FFPE): p53 (r) detected with R.T.U. VECTASTAIN Elite ABC Reagent and DAB (brown). No counterstain.

## Staining Procedure

- For paraffin sections, deparaffinize and hydrate through xylenes or other clearing agents and graded alcohol series.  
For frozen sections or cell preparations fix with acetone or an appropriate fixative for the antigen under study, if necessary.  
Wash for 5 minutes in tap water.
- If antigen unmasking is required, perform this procedure using a Vector® Antigen Unmasking Solution, Citrate-based, pH 6.0 (H-3300) or Tris-based pH 9.0 (H-3301).
- If quenching of endogenous peroxidase activity is required, incubate the slides in BLOXALL® Blocking Solution (SP-6000) for 10 minutes.
- Wash in buffer for 5 minutes.
- Incubate for 20 minutes with 2.5% normal blocking serum.
- Tip off excess serum from sections.
- Incubate for 30 minutes with primary antibody diluted in buffer with 2.5 % normal serum.
- Wash for 5 minutes in buffer.
- Incubate for 30 minutes with a diluted biotinylated secondary antibody.
- Wash for 5 minutes in buffer.
- Incubate for 30 minutes with the R.T.U. VECTASTAIN Elite ABC Reagent.
- Wash for 5 minutes in buffer.
- 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 listings, specifications, protocols and additional information is available on our website: [vectorlabs.com](http://vectorlabs.com)