

VECTASTAIN® Elite® ABC Universal PLUS Kit Peroxidase

(Horse Anti-Mouse/Rabbit IgG)

Cat. No.: PK-8200

Storage: 2-8 °C

Instructions for Immunohistochemical staining using mouse or rabbit primary antibodies.

DESCRIPTION

The VECTASTAIN Elite ABC Universal PLUS Kit is an enzymatic, avidin/biotin based amplification system that produces crisp, highly sensitive, specific staining with low background.

The reagents in the VECTASTAIN Elite ABC Universal PLUS Kit are ready-to-use — no mixing or titering is necessary to obtain optimal staining. Dilution of the reagents or changes in suggested incubation time may affect performance. The reagents are supplied in convenient dropper bottles. (To remove the drop dispenser tip, press laterally with thumb until tip snaps off).

A number of different wash buffers can be used with the VECTASTAIN Elite ABC System. One of the most common is 10 mM sodium phosphate, pH 7.5, 0.9% saline (PBS).

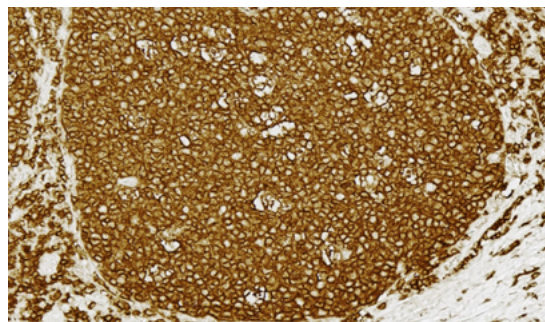
KIT COMPONENTS

Product Name	Volume
BLOXALL® Blocking Solution	50 ml
Normal Horse Serum, 2.5%	50 ml
Horse Anti-Mouse/Rabbit IgG, Biotinylated	50 ml
VECTASTAIN Elite ABC Reagent	50 ml
ImmPACT® DAB EqV Reagent 1 (Chromogen)	25 ml
ImmPACT DAB EqV Reagent 2 (Diluent)	25 ml

The VECTASTAIN Elite ABC Universal PLUS Kit will stain approximately 500 sections.

STORAGE:

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



Tonsil (FFPE): Stained for CD20 antigen using a mouse monoclonal antibody and VECTASTAIN Elite ABC Universal PLUS Kit (PK-8200). No counterstain.

STAINING PROCEDURE

1. 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.

2. 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).
3. If quenching of endogenous peroxidase activity is required, incubate the slides in BLOXALL Blocking Solution for 10 minutes.
4. Wash in buffer for 5 minutes.
5. Incubate for 20 minutes with Normal Horse Serum, 2.5%.
6. Tip off excess serum from sections.
7. Incubate with mouse or rabbit primary antibody diluted in appropriate antibody diluent solution, such as diluted normal horse serum or BSA.
8. Wash for 5 minutes in buffer.
9. Incubate for 30 minutes with prediluted biotinylated horse anti-mouse/rabbit IgG secondary antibody.
10. Wash for 5 minutes in buffer.
11. Incubate for 30 minutes with VECTASTAIN Elite ABC Reagent.
12. Wash for 5 minutes in buffer.
13. Mix ImmPACT DAB EqV solutions in 1:1 ratio for required volume and incubate on section until appropriate stain intensity develops.
14. Rinse in tap water.
15. Counterstain (optional), clear and mount.

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