

Vector® AEC Peroxidase Substrate

Cat. No.: SK-4200

Storage: 2-8 °C

DESCRIPTION

This kit contains all of the reagents necessary to prepare the substrate working solution. Vector AEC Substrate (3-amino-9-ethylcarbazole) produces a red reaction product in the presence of peroxidase (HRP) enzyme. Vector AEC Substrate must be aqueously mounted.

COMPONENTS

Product Name	Volume
Vector AEC Reagent 1	6 ml
Vector AEC Reagent 2	6 ml
Vector AEC Reagent 3	6 ml

STORAGE:

- Store reagents in original bottles at 2-8 °C
- Avoid storing reagents or working solution in strong direct light

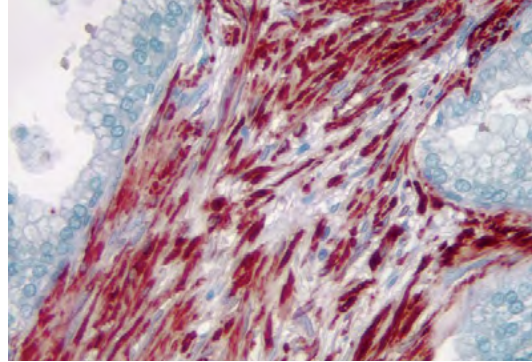
PREPARATION OF SUBSTRATE WORKING SOLUTION:

- To 5 ml of distilled water:
- Add 2 drops ($\approx 72 \mu\text{l}^*$) of Vector AEC Reagent 1
- Add 3 drops ($\approx 90 \mu\text{l}^*$) of Vector AEC Reagent 2
- Add 2 drops ($\approx 80 \mu\text{l}^*$) of Vector AEC Reagent 3

- Mix well before use. Use immediately.

* Drop volumes differ due to solvent compositions.

IMPORTANT: Little is known about the toxicity and carcinogenicity of the substrate kit components. Appropriate care should be exercised when using this reagent including gloves, eye protection, lab coats, and good laboratory procedures. Dispose in accordance with local regulations.



Prostate: Smooth muscle actin (m) detected with ImmPRESS® Anti-Mouse IgG Kit and Vector AEC Substrate (red). Counterstained with Vector Hematoxylin QS (blue).

INSTRUCTIONS FOR USE:

After incubation with a peroxidase (HRP) detection system, rinse sections in buffer. Incubate with the substrate working solution at room temperature for 10-30 minutes. Optimal development times should be determined by the investigator.

Wash for 5 minutes in water.

Counterstain, if desired, with Vector Hematoxylin or Vector Hematoxylin QS. (See counterstain compatibility chart on reverse side.)

For aqueous mounting: Coverslip using an aqueous mounting media such as VectaMount AQ Mounting Medium (H-5501). Do not dehydrate in ethanol.

AEC reaction product is soluble in alcohols and other solvents.

NOTES

We recommend using glass-distilled water in the preparation of the substrate buffer. Deionized water may contain inhibitors of the peroxidase reaction.

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