

ImmPACT[®] AEC

Peroxidase Substrate

Cat. No.	SK-4205
Storage	Store reagents in original bottles at 2-8 °C. Avoid storing reagents or working solution in strong direct light.
Description	This kit contains all of the reagents necessary to prepare a working solution of ImmPACT AEC (3-amino-9-ethylcarbazole). ImmPACT AEC Substrate produces a red reaction product in the presence of peroxidase (HRP) enzyme.

Components

Product Name	Volume
ImmPACT AEC Diluent	120 ml
ImmPACT AEC - Reagent 1	3 ml
ImmPACT AEC - Reagent 2	3 ml
ImmPACT AEC - Reagent 3	3 ml

Preparation of Substrate Working Solution

To 5 ml of ImmPACT AEC Diluent:

- Add 2 drops ($\approx 64 \mu\text{l}^*$) of ImmPACT AEC Reagent 1
- Add 3 drops ($\approx 90 \mu\text{l}^*$) of ImmPACT AEC Reagent 2
- Add 2 drops ($\approx 80 \mu\text{l}^*$) of ImmPACT AEC Reagent 3

Mix well before use.

* Drop volumes differ due to solvent compositions.

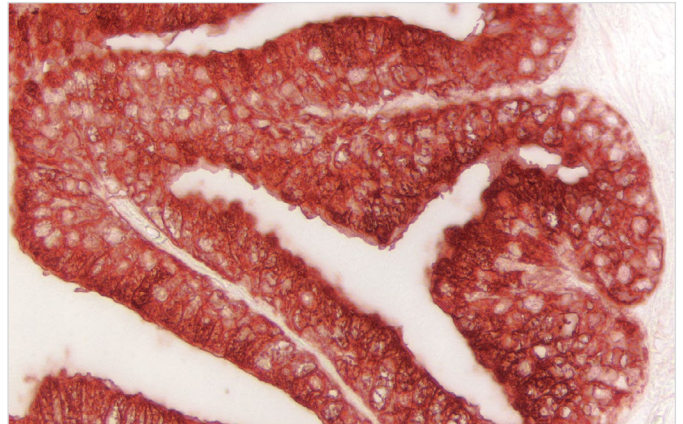
IMPORTANT: Little is known about the toxicity and carcinogenicity of the substrate kit components. Appropriate care should be exercised when using this reagent including gloves, eye protection, lab coats, and good laboratory procedures. Dispose in accordance with local regulations.

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.



Prostate: Prostate Specific Antigen (m) detected with ImmPRESS[®] Polymer Reagent (HRP) and ImmPACT AEC Substrate (red). No Counterstain.

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

ImmPACT AEC reaction product is soluble in alcohols and other solvents.

Notes

- 1) Unused Working Solution is stable up to 14 days if stored at 2-8 °C. Working solution may change color during storage but this will have no effect on the quality or intensity of the staining. If a precipitate forms, filter through a 0.45 μm filter.
- 2) It is not necessary or recommended to add detergent to ImmPACT AEC working solution to reduce surface tension (e.g. for use in an automated stainer).
- 3) This product is not recommended for use on nitrocellulose membranes.

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