

ImmPACT[®] DAB EqV

Peroxidase Substrate

Cat. No.	SK-4103
Storage	Store reagents in original bottles at 2-8 °C. Avoid storing reagents or working solution in strong direct light.
Description	The ImmPACT DAB EqV Substrate contains two reagents necessary to prepare a working solution of diaminobenzidine (DAB). In the presence of peroxidase (HRP) enzyme, ImmPACT DAB EqV Substrate produces a brown reaction product.

Components

Product Name	Volume
ImmPACT DAB EqV Reagent 1 (Chromogen)	50 ml or 200 ml
ImmPACT DAB EqV Reagent 2 (Diluent)	50 ml or 200 ml

Preparation of Substrate Working Solution

- Mix an equal volume of ImmPACT DAB EqV Reagent 1 with ImmPACT DAB EqV Reagent 2.
- Mix well before use. Use immediately.

IMPORTANT: DAB is a suspected carcinogen. 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

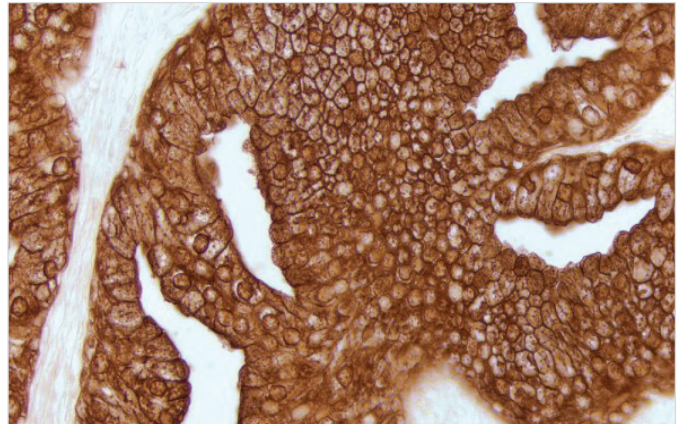
After incubation with a peroxidase (HRP) detection system, rinse sections in buffer. Incubate tissue sections or cells with the substrate working solution at room temperature for 2-10 minutes. Optimal development times should be determined by the investigator.

Wash for 5 minutes in water.

Counterstain if desired.

For permanent, non-aqueous mounting: dehydrate, clear, and coverslip using a non-aqueous mounting media, such as VectaMount[®] Mounting Medium (H-5000) or VectaMount[®] Express Mounting Medium (H-5700).

For aqueous mounting: Coverslip using an aqueous mounting media such as VectaMount[®] AQ Mounting Medium (H-5501).



Prostate: Anti-PSA (m) detected with ImmPRESS[®] Horse Anti-Mouse IgG Kit (HRP) and ImmPACT DAB EqV Substrate (brown). No Counterstain.

Notes

- 1) Working solution is stable for several days if stored at 2-8 °C. If a precipitate forms filter through a 0.45µm filter.
- 2) It is not necessary or recommended to add detergent to ImmPACT DAB EqV working solution to reduce surface tension (e.g. for use in an automated stainer).
- 3) The DAB reaction product can be intensified using a DAB Enhancing Solution (H-2200) after development.
- 4) This product is not recommended for use on nitrocellulose membranes. Instead use Vector DAB or ImmPACT DAB.

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