

ImmPACT[®] DAB

Peroxidase Substrate

Cat. No.	SK-4105
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 diaminobenzidine (DAB). In the presence of peroxidase (HRP) enzyme, the ImmPACT DAB Substrate produces a brown reaction product.

Components

Product Name	Volume
ImmPACT DAB Reagent 1	3.6 ml
ImmPACT DAB Diluent	120 ml

Preparation of Substrate Working Solution

- Add 1 drop (≈ 30µl) of ImmPACT DAB Reagent 1 to 1 ml ImmPACT DAB Diluent.
- Mix well before use.

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

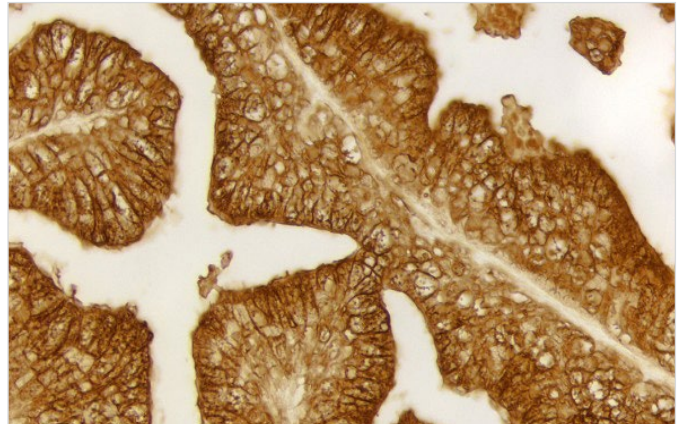
After incubation with a peroxidase (HRP) detection system, rinse sections in buffer. Incubate with the substrate working solution at room temperature for 2-10 minutes. Optimal development times should be determined by the investigator.

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: Prostate Specific Antigen (m) detected with ImmPRESS[®] Reagent (HRP) and ImmPACT DAB (brown). No counterstain.

Notes

- 1) Unused working solution is stable up to 14 days if stored at 2-8 °C or up to 5 days if stored at room temperature (approximately 23 °C). Working solution may change color during storage but this will have no effect on the quality or intensity of the staining. No sample-obscuring precipitate will form.
- 2) It is not necessary or recommended to add detergent to ImmPACT DAB 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.

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