

## ImmPACT® DAB Peroxidase Substrate

**Cat. No.:** SK-4105

**Storage:** 2-8 °C

### 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

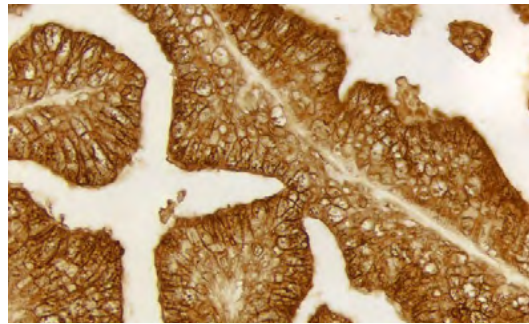
### STORAGE:

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

### PREPARATION OF SUBSTRATE WORKING SOLUTION:

- Add 1 drop (≈ 30 µl) 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.



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

### 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. (See counterstain compatibility chart on reverse side.)

**For permanent, non-aqueous mounting:** Dehydrate, clear and coverslip using a non-aqueous mounting media, such as VectaMount® Mounting Medium (H-5000).

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

### 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](http://vectorlabs.com)