



ImmPACT™ DAB Peroxidase Substrate

Cat. No. SK-4105

Store at 2-8 °C

Burlingame, CA USA

ImmPACT™ DAB Substrate

The **ImmPACT™ DAB** Substrate contains all of the reagents necessary to prepare a working solution of diaminobenzidine (DAB) for staining tissue sections or cells, or for use on nitrocellulose, nylon, or other membranes. In the presence of peroxidase (HRP) enzyme, the **ImmPACT™ DAB** Substrate produces a brown reaction product.

REAGENTS:

- 120 ml **ImmPACT™ DAB** Diluent
- 3.6 ml **ImmPACT™ DAB** Chromogen concentrate

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 (approximately 30 μ l) of **ImmPACT™ DAB** Chromogen concentrate to 1 ml **ImmPACT™ DAB** Diluent
- Mix well before use

INSTRUCTIONS FOR USE:

After incubation with a peroxidase (HRP) detection system, rinse well. 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.) Coverslip with either a non-aqueous mounting medium such as **VectaMount™** (H-5000) or an aqueous mounting medium such as **VectaMount™ AQ** (H-5501).

NOTES:

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.

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).

The DAB reaction product can be intensified using a DAB Enhancing Solution (H-2200) after development.

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.

For Laboratory Use

04/14