



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, nitrocellulose, nylon, or other membranes. The ImmPACT™ DAB substrate produces a brown reaction product.

REAGENTS:

120 ml ImmPACT™ DAB Diluent

3.6 ml ImmPACT™ DAB Chromogen concentrate

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

PREPARATION OF SUBSTRATE WORKING SOLUTION:

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

INSTRUCTIONS FOR USE:

After incubation with a peroxidase (HRP) detection system, rinse well. Incubate tissue sections or blots with the substrate working solution at room temperature until suitable staining develops. Development times should be determined by the investigator but generally 2-10 minutes provides good staining intensity.

Wash for 5 minutes in water.

For tissue sections: Counterstain if desired. Coverslip with either a permanent mounting media (such as **VectaMount™**, H-5000) or an aqueous-based mounting media (such as **VectaMount™ AQ**, H-5501).

For blots: Air dry.

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 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, Cat. No. 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. After completion of the staining procedure, the diluted working solution should be discarded into an equal volume of 3% potassium permanganate (KMnO_4), 2% sodium carbonate (Na_2CO_3) in deionized or distilled water. Any containers used in the staining procedure should be washed in the same solution to neutralize the DAB. Dispose in accordance with local regulations.