



ImmPACT™ DAB EqV **Peroxidase Substrate** Cat. No. SK-4103

Store at 2-8 °C

Burlingame, CA USA

ImmPACT™ DAB EqV Substrate

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

REAGENTS:

- 200 ml **ImmPACT™ DAB EqV** Reagent 1
- 200 ml **ImmPACT™ DAB EqV** Reagent 2

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:

- Mix an equal volume of **ImmPACT™ DAB EqV** Reagent 1 with **ImmPACT™ DAB EqV** Reagent 2
- Mix well before use

INSTRUCTIONS FOR USE:

After incubation with a peroxidase (HRP) detection system, rinse well. 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. (See counterstain compatibility chart on reverse side.) Coverslip with either a non-aqueous mounting medium such as **VectaMount™** (Cat. No. H-5000) or an aqueous mounting medium such as **VectaMount™ AQ** (Cat. No. H-5501).

NOTES:

Working solution is stable for several days if stored at 2-8 °C. If a precipitate forms filter through a 0.45 µm filter.

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).

The DAB reaction product can be intensified using a DAB Enhancing Solution, Cat. No. H-2200, after development.

This product is not recommended for use on nitrocellulose membranes. Instead use Vector® DAB (Cat. No. SK-4100) or **ImmPACT™ DAB** (Cat. No. SK-4105).

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.