



ImmPACT™ SG Peroxidase Substrate

Cat. No. SK-4705

Store at 2-8 °C

Burlingame, CA USA

ImmPACT™ SG Substrate

The **ImmPACT™ SG** Substrate contains all of the reagents necessary to prepare a working solution for staining tissue sections or cells, or for use on nylon and PVDF membranes. In the presence of peroxidase (HRP) enzyme, the **ImmPACT™ SG** Substrate produces a blue/gray reaction product.

REAGENTS:

- 120 ml **ImmPACT™ SG** Diluent
- 2.5 ml **ImmPACT™ SG** Reagent 1
- 3 ml **ImmPACT™ SG** 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:

- To 5 ml of **ImmPACT™ SG** Diluent:
 - Add 3 drops (approximately 90 μ l*) of **ImmPACT™ SG** Reagent 1
 - Add 3 drops (approximately 120 μ l*) of **ImmPACT™ SG** Reagent 2
- 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 2-15 minutes. Optimal development times should be determined by the investigator.

Wash slides for 5 minutes in water.

Counterstain, if desired, with Vector® Nuclear Fast Red (Cat. No. H-3403) for best contrast. (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:

Unused working solution is stable up to 14 days if stored at 2-8 °C. Working solution may change color during storage but this will have no effect on the quality or intensity of the staining. If a precipitate forms filter through a 0.45 μ m filter.

It is not necessary or recommended to add detergent to **ImmPACT™ SG** working solution to reduce surface tension (e.g. for use in an automated stainer).

This product is not recommended for use on nitrocellulose membranes.

IMPORTANT: Little is known about the toxicity and carcinogenicity of the substrate kit components. 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.

** Drop volumes differ due to solvent compositions.*