

ImmPACT® SG
Peroxidase Substrate

Cat. No.: SK-4705

Storage: 2-8 °C

DESCRIPTION

This kit contains all of the reagents necessary to prepare a working solution. In the presence of peroxidase (HRP) enzyme, the ImmPACT SG Substrate produces a blue/gray reaction product.

COMPONENTS

<u>Product Name</u>	<u>Volume</u>
ImmPACT SG Diluent	120 ml
ImmPACT SG Reagent 1	2.5 ml
ImmPACT SG Reagent 2	3 ml

STORAGE:

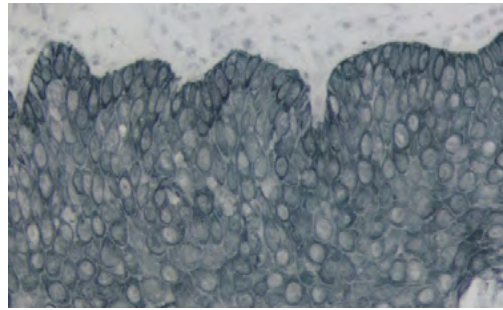
- Store reagents in original bottles at 2-8 °C
- Avoid storing reagents or working solution in strong direct light

PREPARATION OF SUBSTRATE WORKING SOLUTION:

- To 5 ml of ImmPACT SG Diluent:
 - Add 3 drops (≈ 90 µl*) of ImmPACT SG Reagent 1
 - Add 3 drops (≈ 120 µl*) of ImmPACT SG Reagent 2
- Mix well before use.

* Drop volumes differ due to solvent compositions.

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.



Tonsil: Cytokeratin AE1/AE3 (m) detected with ImmPRESS® Polymer Reagent (HRP) and ImmPACT SG Substrate (blue-gray). 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 2-15 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 for best contrast. (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. 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 µm filter.
- 2) 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).
- 3) This product is not recommended for use on nitrocellulose membranes.

Detailed product listings, specifications, protocols and additional information is available on our website: vectorlabs.com