

ImmPACT® VIP

Peroxidase Substrate

Cat. No.	SK-4605
Storage	Store reagents in original bottles at 2-8 °C. Avoid storing reagents or working solution in strong direct light.
Description	This kit contains all of the reagents necessary to prepare a working solution. In the presence of peroxidase (HRP) enzyme, the ImmPACT VIP Substrate produces a purple reaction product.

Components

Product Name	Volume
ImmPACT VIP Diluent	120 ml
ImmPACT VIP Reagent 1	3 ml
ImmPACT VIP Reagent 2	2.5 ml
ImmPACT VIP Reagent 3	2.5 ml
ImmPACT VIP Reagent 4	3 ml

Preparation of Substrate Working Solution

To 5 ml of ImmPACT VIP Diluent:

- Add 3 drops (≈ 90 µl*) of ImmPACT VIP Reagent 1
- Add 3 drops (≈ 75 µl*) of ImmPACT VIP Reagent 2
- Add 3 drops (≈ 75 µl*) of ImmPACT VIP Reagent 3
- Add 3 drops (≈ 120 µl*) of ImmPACT VIP Reagent 4

Mix well before use.

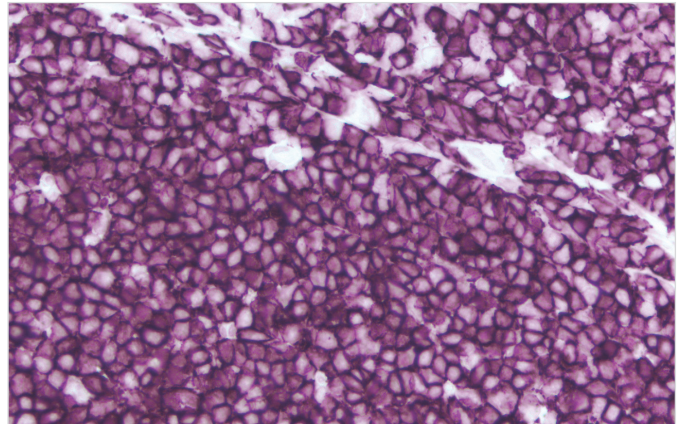
* Drop volumes differ due to solvent compositions.

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.

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.



Tonsil: CD20 (m) detected with ImmPRESS® Polymer Reagent (HRP) and ImmPACT VIP Substrate (purple). No counterstain.

Counterstain, if desired, with Vector® Hematoxylin, Vector Hematoxylin QS or Vector Methyl Green. Note: ImmPACT VIP is not compatible with all methyl green formulations.

For permanent, non-aqueous mounting: Dehydrate, clear, and coverslip using a non-aqueous mounting media, such as VectaMount® Mounting Medium (H-5000) or VectaMount® Express Mounting Medium (H-5700). Do not use inferior alcohol grades with this product.

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 VIP working solution to reduce surface tension (e.g. for use in an automated stainer).
- 3) Prolonged incubation in alcohol or use of alcohol-based differentiating solutions may decrease sensitivity.
- 4) 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