

ImmPACT NovaRED®

Peroxidase Substrate

Cat. No.	SK-4805
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 NovaRED Substrate produces a red reaction product.

Components

Product Name	Volume
ImmPACT NovaRED Diluent	120 ml
ImmPACT NovaRED Reagent 1	3 ml
ImmPACT NovaRED Reagent 2	1.75 ml
ImmPACT NovaRED Reagent 3	1.75 ml
ImmPACT NovaRED Reagent 4	2.5 ml

Preparation of Substrate Working Solution

To 5 ml of ImmPACT NovaRED Diluent:

- Add 3 drops (≈ 80 µl*) of ImmPACT NovaRED Reagent 1
- Add 2 drops (≈ 50 µl*) of ImmPACT NovaRED Reagent 2
- Add 2 drops (≈ 50 µl*) of ImmPACT NovaRED Reagent 3
- Add 2 drops (≈ 80 µl*) of ImmPACT NovaRED Reagent 4

Mix well before use.

* Drop volumes differ due to solvent compositions.

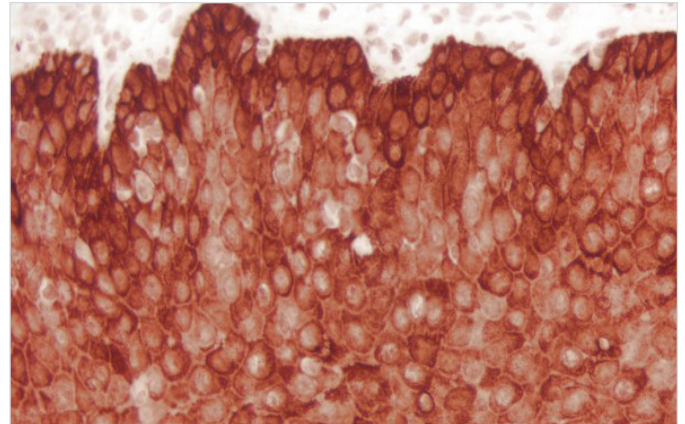
IMPORTANT: Little is known about the toxicity and carcinogenicity of the substrate kit components. Appropriate care should be exercised when using this reagent including gloves, eye protection, lab coats, and good laboratory procedures. Dispose in accordance with local regulations.

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.

Counterstain, if desired, with Vector® Hematoxylin, Vector Hematoxylin QS or Vector Methyl Green.



Tonsil: Cytokeratin (AE1/AE3 (m) detected with ImmPRESS® Reagent (HRP) and ImmPACT NovaRED Substrate (red). No counterstain.

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 NovaRED 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