

# ImmPACT® AMEC Red

## Peroxidase Substrate

<b>Cat. No.</b>	<b>SK-4285</b>
<b>Storage</b>	Store reagents in original bottles at 2-8 °C. Avoid storing reagents or working solution in strong direct light.
<b>Description</b>	This kit contains all of the reagents necessary to prepare a working solution. In the presence of peroxidase (HRP) enzyme, the ImmPACT AMEC Red Substrate produces a red reaction product that is a superior AEC substitute. Specimens developed with ImmPACT AMEC Red must be aqueously mounted.

### Components

Product Name	Volume
ImmPACT AMEC Red Diluent	120 ml
ImmPACT AMEC Red Reagent 1	3 ml
ImmPACT AMEC Red Reagent 2	3 ml

### Preparation of Substrate Working Solution

To 5 ml of ImmPACT AMEC Red Diluent:

- Add 3 drops (≈ 90 µl\*) of ImmPACT AMEC Red Reagent 1
- Add 2 drops (≈ 80 µl\*) of ImmPACT AMEC Red Reagent 2

Mix well before use.

\* Drop volumes differ due to solvent compositions.

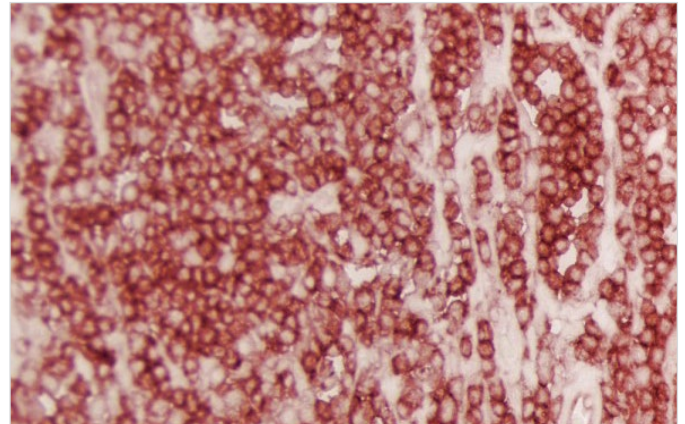
**IMPORTANT:** Little is known about the toxicity and carcinogenicity of the substrate kit components. Appropriate care should be exercised when using this reagent including gloves, eye protection, lab coats, and good laboratory procedures. Dispose in accordance with local regulations.

### 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 10-30 minutes. Optimal development times should be determined by the investigator.

Wash for 5 minutes in water.

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



*Tonsil: Leukocyte Common Antigen (LCA) (m) detected with ImmPRESS® Anti-mouse IgG Polymer Kit (HRP) and ImmPACT AMEC Red (red). No Counterstain.*

For aqueous mounting: Coverslip using an aqueous mounting media such as VectaMount® AQ Mounting Medium. Do not dehydrate in ethanol.

ImmPACT AMEC Red reaction product is soluble in alcohols and other solvents.

### 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 AMEC Red 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](http://vectorlabs.com)