

## ImmPACT® AMEC Red Peroxidase Substrate

**Cat. No.:** SK-4285

**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 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

<u>Product Name</u>	<u>Volume</u>
ImmPACT AMEC Red Diluent	120 ml
ImmPACT AMEC Red Reagent 1	3 ml
ImmPACT AMEC Red 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:

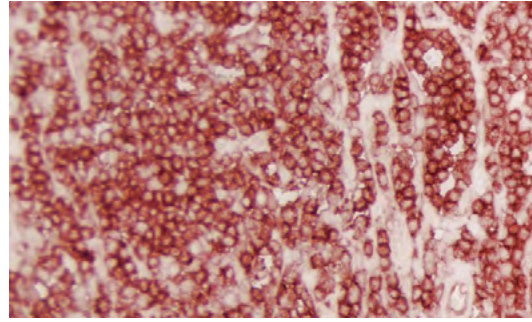
To 5 ml of ImmPACT AMEC Red Diluent:

- Add 3 drops ( $\approx 90 \mu\text{l}^*$ ) of ImmPACT AMEC Red Reagent 1
- Add 2 drops ( $\approx 80 \mu\text{l}^*$ ) 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.



*Tonsil: Leukocyte Common Antigen (LCA) (m) detected with ImmPRESS® Anti-mouse IgG Polymer Kit (HRP) and ImmPACT AMEC Red (red). 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 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. (See counterstain compatibility chart on reverse side.)

**For aqueous mounting:** Coverslip using an aqueous mounting media such as VectaMount AQ Mounting Medium (H-5501). 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  $\mu\text{m}$  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)