



ImmPACT™ AMEC *Red*
Peroxidase Substrate
Cat. No. SK-4285

Store at 2-8 °C

Burlingame, CA USA

ImmPACT™ AMEC *Red* Substrate

The ImmPACT™ AMEC *Red* Substrate contains all of the reagents necessary to prepare a working solution for staining tissue sections or cells, or for use on nylon and PVDF membranes. 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.

REAGENTS:

120 ml ImmPACT™ AMEC *Red* Diluent
3 ml ImmPACT™ AMEC *Red* Reagent 1
3 ml ImmPACT™ AMEC *Red* Reagent 2

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 (approximately 90 µl*) of ImmPACT™ AMEC *Red* Reagent 1
 - Add 2 drops (approximately 80 µl*) of ImmPACT™ AMEC *Red* Reagent 2
- Mix well before use

INSTRUCTIONS FOR USE:

After incubation with a peroxidase (HRP) detection system, rinse well. 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 (Cat. No. H-3401) or Vector® Hematoxylin QS (Cat. No. H-3404). (See counterstain compatibility chart on reverse side.) Do not dehydrate. Coverslip with an aqueous-based mounting medium such as *VectaMount™ AQ* (Cat. No. H-5501).

NOTES:

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.

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).

This product is not recommended for use on nitrocellulose membranes.

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.

* Drop volumes differ due to solvent compositions.