

## ImmPACT® Vector® Red Alkaline Phosphatase Substrate

**Cat. No.:** SK-5105

**Storage:** 2-8 °C

### DESCRIPTION

This kit contains all of the reagents necessary to prepare a working solution. The ImmPACT Vector Red Substrate produces a red reaction product in the presence of alkaline phosphatase (AP) enzyme. This reaction product is also highly fluorescent and non-fading and can be viewed with Texas Red or rhodamine filter systems.

### COMPONENTS

<u>Product Name</u>	<u>Volume</u>
ImmPACT Vector Red Diluent	120 ml
ImmPACT Vector Red Reagent 1	2.5 ml
ImmPACT Vector Red Reagent 2	2.5 ml

### STORAGE:

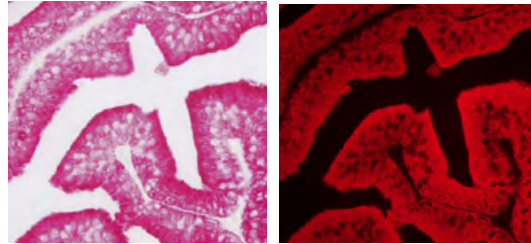
- Store reagents in original bottles at 2-8 °C
- Avoid storing reagents or working solution in strong direct light

### PREPARATION OF SUBSTRATE WORKING SOLUTION:

- To 5 ml ImmPACT Vector Red Diluent:  
Add 2 drops (≈ 80 µl\*) of ImmPACT Vector Red Reagent 1  
Add 2 drops (≈ 60 µl\*) of ImmPACT Vector Red Reagent 2†
- Mix well before use. Use immediately.

\* 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.



Prostate: Anti-PSA (m) detected with ImmPRESS™-AP Anti-Mouse IgG Reagent and ImmPACT Vector Red Alkaline Phosphatase Substrate. Viewed under brightfield and fluorescence microscopy.

### INSTRUCTIONS FOR USE:

After incubation with an AP detection system, rinse sections in buffers. Incubate with the substrate working solution for 20-30 minutes. Optimal development times should be determined by the investigator. Improved staining may be obtained by developing the substrate in the dark.

Wash slides in buffer for 5 minutes. Rinse in water and counterstain, if desired (see chart on reverse).

**For permanent, non-aqueous mounting:** Dehydrate, clear and coverslip using a non-aqueous mounting media, such as VectaMount® Mounting Medium (H-5000).

**For aqueous mounting:** Coverslip using an aqueous mounting media such as VectaMount AQ Mounting Medium (H-5501) or a VECTASHIELD® Antifade Mounting Medium. To increase the intensity of ImmPACT Vector Red fluorescence, rinse slides 2-5 minutes in 100% ethanol before mounting.

Detailed product listings, specifications, protocols and additional information is available on our website: [vectorlabs.com](http://vectorlabs.com)