

Vector® Red

Alkaline Phosphatase Substrate

Cat. No.: SK-5100

Storage: 2-8 °C

DESCRIPTION

This kit contains all the reagents necessary to prepare the substrate working solution. 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>
Vector Red Reagent 1	4 ml
Vector Red Reagent 2	4 ml
Vector Red Reagent 3	4 ml

STORAGE:

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

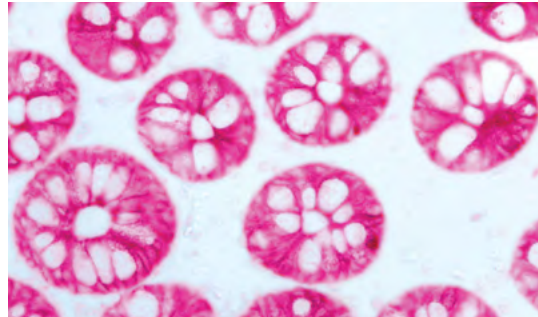
PREPARATION OF SUBSTRATE WORKING SOLUTION:

- To 5 ml of 100 mM - 200 mM Tris-HCl, pH 8.2 - 8.5 buffer*
 - Add 2 drops (≈80 µl) of Vector Red Reagent 1
 - Add 2 drops (≈80 µl) of Vector Red Reagent 2
 - Add 2 drops (≈80 µl) of Vector Red Reagent 3

- Mix well before use. Use immediately.

* 0.1% Tween 20 in the Vector Red working solution is recommended, especially for use with antigen retrieved tissues.

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.



Colon: Cytokeratin AE1/AE3 (m) detected with ImmPRESS®-AP Polymer Detection System and Vector Red Substrate (magenta). No counterstain.

INSTRUCTIONS FOR USE:

After incubation with an AP detection system, rinse sections in buffer. 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 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