

Vector® VIP
 Peroxidase Substrate

Cat. No.: SK-4600

Storage: 2-8 °C

DESCRIPTION

This kit contains all of the reagents necessary to prepare the substrate working solution. Vector VIP Substrate produces a purple reaction product in the presence of peroxidase (HRP) enzyme. Vector VIP Substrate is also suitable for darkfield and electron microscopy (EM).

COMPONENTS

Product Name	Volume
Vector VIP Reagent 1	7 ml
Vector VIP Reagent 2	5 ml
Vector VIP Reagent 3	5 ml
Vector VIP Reagent 4	7.2 ml

STORAGE:

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

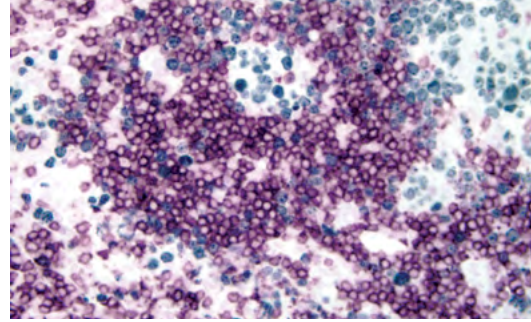
PREPARATION OF SUBSTRATE WORKING SOLUTION:

- To 5 ml of PBS (10 mM sodium phosphate, 0.9% saline, pH 7.5)
- Add 3 drops (≈ 90 µl†) of Vector VIP Reagent 1
- Add 3 drops (≈ 75 µl†) of Vector VIP Reagent 2
- Add 3 drops (≈ 75 µl†) of Vector VIP Reagent 3
- Add 3 drops (≈ 120 µl†) of Vector VIP Reagent 4

-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.



Tonsil (Double Label): CD3 (m) detected with ImmPRESS® Reagent (HRP) Anti-Mouse IgG and Vector VIP Substrate (purple) • Ki67 (m) detected with ImmPRESS Reagent (HR) Anti-Mouse IgG and Vector SG Substrate (blue-gray). No Counterstain.

INSTRUCTIONS FOR USE:

After incubation with a peroxidase (HRP) detection system, rinse sections in buffer. Incubate with the substrate working solution for 2-15 minutes. Optimal development times should be determined by the investigator.

Wash slides for 5 minutes in water.

Counterstain, if desired, with Vector Hematoxylin, Vector Hematoxylin QS, or Vector Methyl Green (See chart on reverse.) Note: Vector VIP is not compatible with all methyl green formulations.

For permanent, non-aqueous mounting: Dehydrate, clear and coverslip using a non-aqueous mounting media, such as VectaMount® Mounting Medium (H-5000). Do not mount with aqueous mounting media.

NOTES

We recommend using glass-distilled water in the preparation of substrate buffer. Deionized water may contain inhibitors of the peroxidase reaction. Solutions containing sodium azide or other inhibitors of peroxidase activity should not be used in diluting the peroxidase substrate.

Detailed product listings, specifications, protocols and additional information are available on our website: vectorlabs.com