

Vector® NovaRED™ Peroxidase Substrate

Cat. No.: SK-4800

Storage: 2-8 °C

DESCRIPTION

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

COMPONENTS

<u>Product Name</u>	<u>Volume</u>
Vector NovaRED Reagent 1	6 ml
Vector NovaRED Reagent 2	3.5 ml
Vector NovaRED Reagent 3	3.5 ml
Vector NovaRED Reagent 4	5 ml

STORAGE:

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

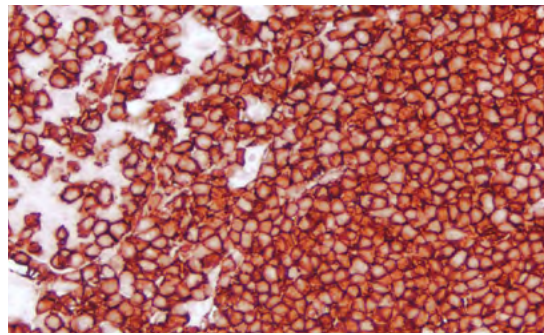
PREPARATION OF SUBSTRATE WORKING SOLUTION:

To 5 ml of distilled water
 Add 3 drops (≈ 81 μl[†]) of Vector NovaRED Reagent 1
 Add 2 drops (≈ 50 μl[†]) of Vector NovaRED Reagent 2
 Add 2 drops (≈ 50 μl[†]) of Vector NovaRED Reagent 3
 Add 2 drops (≈ 80 μl[†]) of Vector NovaRED 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: CD20 (m) detected with ImmPRESS® HRP Anti-Mouse IgG and Vector NovaRED Substrate (red). No counterstain.

INSTRUCTIONS FOR USE:

After incubation with a peroxidase (HRP) detection system, rinse sections in buffer. Incubate with the substrate working solution for 5-15 minutes. Optimal development times should be determined by the investigator.

Wash slides with 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). Do not mount with aqueous mounting media.

NOTES

- 1) We recommend using glass-distilled water in the preparation of the substrate. Deionized water may contain inhibitors of the peroxidase reaction.
- 2) The color of the reagent solutions may darken with time. This will have no effect on the quality or intensity of the staining.
- 3) Prolonged incubation in alcohol or use of alcohol-based differentiating solutions may decrease sensitivity.

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