

## Levamisole Solution

**Cat. No.:** SP-5000

**Storage:** Refrigerate 2-8 °C



### Description:

Levamisole will inhibit most forms of alkaline phosphatase other than the intestinal isoenzyme. Since the intestinal form of the enzyme is used as the marker enzyme in VECTASTAIN ABC-AP Kits, levamisole can be used to inhibit endogenous alkaline Phosphatase in most cases. Because the inhibition produced by levamisole is reversible, the inhibitor must be added to the substrate solution.

### COMPONENTS

<u>Product Name</u>	<u>Volume</u>
Levamisole Solution	18 ml

### Guidelines for Use:

When used in conjunction with Vector Laboratories' alkaline phosphatase substrate kits, add one drop of Levamisole Solution to 5 ml of alkaline phosphatase substrate buffer and mix well. Then add substrate reagents and continue as outlined in the substrate kit instructions.

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