



Levamisole Solution

Cat. No. SP-5000

Concentration 125 mM

Storage Refrigerate

REMARKS:

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 the VECTASTAIN® ABC-AP Kit, 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.

This lot of levamisole was tested on two frozen tissues and was found to completely inhibit endogenous alkaline phosphatase activity.

INSTRUCTIONS 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 instructions for the Alkaline Phosphatase Substrate Kit.