

Phosphatase Inhibitor Cocktail

Catalog No: 37492, 37493

Format: 1 ml, 5 ml

Quality Control:

Active Motif's Phosphatase Inhibitor Cocktail is a broad-spectrum cocktail designed to protect phosphoproteins from dephosphorylation during lysis and extraction procedures from mammalian cells or tissues and for immunoprecipitation and kinase assays. The cocktail solution inhibits both serine/threonine phosphatases and protein tyrosine phosphatases (PTPs). Phosphatase Inhibitor Cocktail is provided as a ready-to-use 100X stock solution. The cocktail composition and inhibitor targets are listed below:

- Sodium Fluoride - Acid Phosphatases
- Sodium Orthovanadate - Protein Tyrosine Phosphatase, Alkaline Phosphatase
- Sodium Pyrophosphate - Serine/Threonine Phosphatases
- β -Glycerophosphate - Serine/Threonine Phosphatases

It is recommended to use 10 μ l of the cocktail solution for every 1 ml of cell lysate. Since phosphatase levels may vary among cell and tissue types, it may be necessary to increase the concentration of inhibitors.

Note: Compounds precipitate out of solution at 4°C. To resuspend, bring the cocktail up to room temperature and shake the solution until any visible crystals completely dissolve into the liquid.

For quality control, each batch of Phosphatase Inhibitor Cocktail is tested for its ability to inhibit alkaline phosphatase activity. (Figure 1) Alkaline phosphatase activity was measured at 0-50 ng/well in the presence or absence of Phosphatase Inhibitor Cocktail. In the presence of Phosphatase Inhibitor Cocktail, no alkaline phosphatase activity is observed, while in the absence of phosphatase inhibitors, increased alkaline phosphatase activity levels are observed with increasing alkaline phosphatase concentrations.

Storage and Guarantee: The Phosphatase Inhibitor Cocktail is provided as a clear solution. The product is shipped at room temperature and should be stored at -20 °C long-term. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.

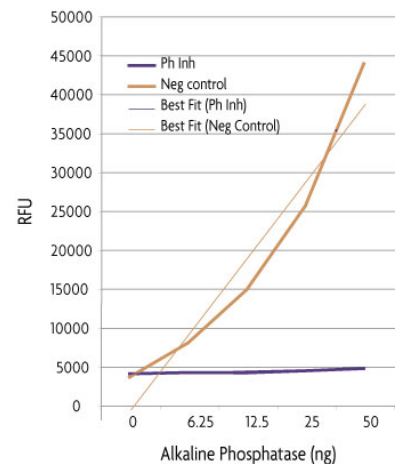


Figure 1: Alkaline Phosphatase Inhibition Assay

Alkaline phosphatase activity was measured at 0-50 ng/well alkaline phosphatase concentrations in the presence (Ph Inh) or absence (Neg control) of Phosphatase Inhibitor Cocktail.