

Sodium Orthovanadate (PHIVA-1mL)

Stable and Potent Phosphatase Inhibitor

DESCRIPTION

Sodium Orthovanadate (CAS Number 13721-39-6, Na_3VO_4) is a general inhibitor for protein phosphotyrosyl phosphatases (PTPs). Inhibition of vanadate is competitive and can be reversed by addition of EDTA or by dilution. Vanadate is commonly used to preserve the protein tyrosyl phosphorylation state in cells and lysates. BioAssay Systems' Vanadate reagent is activated for maximum inhibition.

Suggested Working Concentration: 1-10 mM

Contents: Catalog#: PHIVA-1mL

Sodium Orthovanadate: 1 mL 100 mM aqueous solution (pH 10.0)

Storage conditions: this product is shipped at ambient temperature. Store at -20°C . Shelf life of 3 years after receipt.

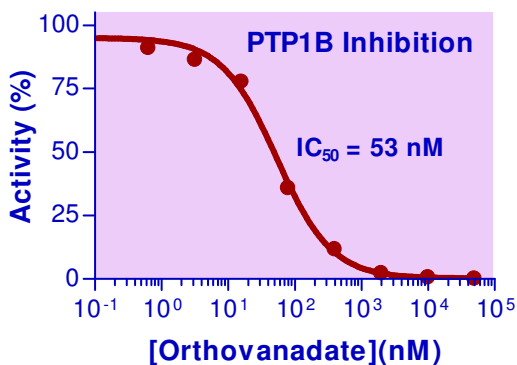
Bulk order of any size (>10 mL) is available. Please contact info@bioassaysys.com.

GENERAL CONSIDERATIONS

EDTA and reducing agents such as DTT can interact with Vanadate and reduce its potency.

LITERATURE

- Gordan, JA (1991) *Methods in Enzymology* 201: 477-82.
- Huyer, G et al (1997) *J. Biol. Chem.* 272: 843-51.



Inhibition of PTP1B Activity

PRECAUTIONS

Reagents are for research use only. Normal precautions for laboratory reagents should be exercised while using the reagents.

Avoid contact and inhalation. Standard laboratory safety procedures should be followed when handling this product. Safety procedures include wearing OSHA approved safety glasses, gloves and protective clothing.

If eye or skin contact occurs, immediately wash affected area with soap and copious amounts of water for at least 15 minutes. Assure adequate flushing by separating the eyelids with fingers. Call a physician to seek medical advice.

If swallowed, wash out mouth with water provided person is conscious and call a physician to seek medical advice.

If inhaled, move individual to fresh air and call a physician to seek medical advice.

TRANSPORT INFORMATION:

DOT (US):	Not regulated
UN Number:	Not regulated
Class:	Not regulated
Packaging Group:	Not regulated

