



## Sodium Orthovanadate (1000x)

ATPase Inhibitor  
Sodium orthovanadate

Cat. No.	Amount
AK-102V-S	100 µl
AK-102V-L	500 µl

**For research use only!** Not intended for human or animal diagnostic or therapeutic uses.

**Shipping:** shipped on blue ice

**Storage Conditions:** store at -20 °C

**Shelf Life:** 12 months after date of delivery

**Molecular Formula:** Na<sub>3</sub>VO<sub>4</sub>

**Molecular Weight:** 183.91 g/mol

**CAS#:** 13721-39-6

**pH:** 10.0

### Description:

Sodium Orthovanadate is an inhibitor of ATPase, alkaline phosphatase and tyrosine phosphatase that can be used in the purification procedure of ATP binding proteins (for further information see the manual of the ATP Affinity Test Kit #AK-102).

### Content:

200 mM activated Sodium Orthovanadate dissolved in water, pH adjusted to 10.0

### Usage:

Dilute 1:1000 in purification buffers.

### Activity:

Inhibitor of ATPase, alkaline phosphatase and tyrosine phosphatase

### Related Products:

Aminophenyl-ATP-Agarose, C10-spacer pre-swollen in 20% ethanol #AC-101

8-[(6-Amino)hexyl]-amino-ATP-Agarose pre-swollen in 20% ethanol #AC-127

N6-(6-Amino)hexyl-ATP-Agarose pre-swollen in 20% ethanol #AC-129

2'/3'-EDA-ATP-Agarose pre-swollen in 20% ethanol #AC-131

PBS Tablets #AK-102P

100x Protease Inhibitor Mix #AK-102I

5x Binding Buffer #AK-102B

5x Wash Buffer #AK-102W

5x Elution Buffer #AK-102E

### Selected References:

Haystead *et al.* (1993) Gamma-phosphate-linked ATP-Sepharose for the affinity purification of protein-kinases - rapid purification to homogeneity of skeletal-muscle mitogen-activated protein-kinase. *Eur. J. Biochem.* **214** (2):459.

Jenö *et al.* (1989) Purification and Characterization of a 40 S Ribosomal Protein S6 Kinase from Vanadate-stimulated Swiss 3T3 Cells. *J. Biol. Chem.* **264**:1293.

McNutt *et al.* (1981) Comparison of cell peripheries in the human colon carcinoma cell lines SW480 and SW620 grown in floating chamber culture, cover slip culture, athymic (nude) mice, and BALB/c mice. *Lab. Invest.* **44**:309.

Trayer *et al.* (1974) Affinity Chromatography of Nicotinamide Nucleotide-Dependent Dehydrogenases on Immobilized Nucleotide Derivates. *Biochem. J.* **139**:609.

Scherer *et al.* (1954) Studies on the propagation in vitro of poliomyelitis viruses. IV. Viral multiplication in a stable strain of human malignant epithelial cells (strain HeLa) derived from an epidermoid carcinoma of the cervix. *J. Exp. Med.* **97**:695.