

## Sodium Orthovanadate, activated (1000x) ATPase Inhibitor

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

### 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, Cat. No. AK-102).

### Composition

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

### Usage

Dilute 1:1000 in purification buffers.

### Storage and Stability

Store at -20°C.

Quality guaranteed for 12 months.

**FOR RESEARCH USE ONLY.  
DO NOT USE FOR HUMAN OR ANIMAL  
DIAGNOSTICS OR THERAPEUTICS.**

### 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 colonic adenocarcinoma 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.

### Associated products available from Jena Bioscience

For detailed information please view the sections on

[www.jenabioscience.com](http://www.jenabioscience.com)

<b>Aminophenyl-ATP-Agarose, C<sub>10</sub>-spacer</b> pre-swollen in 20% ethanol	Cat.# AC-101
<b>8-[(6-Amino)hexyl]-amino-ATP-Agarose</b> pre-swollen in 20% ethanol	Cat.# AC-127
<b>N<sup>6</sup>-(6-Amino)hexyl-ATP-Agarose</b> pre-swollen in 20% ethanol	Cat.# AC-129
<b>2'/3'-EDA-ATP-Agarose</b> pre-swollen in 20% ethanol	Cat.# AC-131
<b>PBS Tablets</b>	Cat.# AK-102P
<b>100x Protease Inhibitor Mix</b>	Cat.# AK-102I
<b>5x Binding Buffer</b>	Cat.# AK-102B
<b>5x Wash Buffer</b>	Cat.# AK-102W
<b>5x Elution Buffer</b>	Cat.# AK-102E