

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

Aminophenyl-ATP-Agarose, C₁₀-spacer pre-swollen in 20% ethanol	Cat.# AC-101
8-[(6-Amino)hexyl]-amino-ATP-Agarose pre-swollen in 20% ethanol	Cat.# AC-127
N⁶-(6-Amino)hexyl-ATP-Agarose pre-swollen in 20% ethanol	Cat.# AC-129
2'/3'-EDA-ATP-Agarose pre-swollen in 20% ethanol	Cat.# AC-131
PBS Tablets	Cat.# AK-102P
100x Protease Inhibitor Mix	Cat.# AK-102I
5x Binding Buffer	Cat.# AK-102B
5x Wash Buffer	Cat.# AK-102W
5x Elution Buffer	Cat.# AK-102E