

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 data sheet of the ATP Affinity Test Kit, cat.# AK-102)

Activated, pH is adjusted to 10.0.

Dilute 1:1000 in purification buffers.

Storage conditions:

Store at -20°C.

Stable for 1 year.

FOR RESEARCH USE ONLY.

NOT INTENDED FOR HUMAN OR ANIMAL DIAGNOSTIC OR THERAPEUTIC USES.

[Associated products available from Jena Bioscience](#)

For detailed information please view the sections on www.jenabioscience.com

Aminophenyl-ATP-Sepharose®, C ₁₀ -linked, pre-swollen in 20% ethanol	Cat.# AC-101
8-[(6-Amino)hexyl]-amino-ATP-Sepharose®, pre-swollen in 20% ethanol	
N ⁶ -(6-Amino)hexyl-ATP -Sepharose®, pre-swollen in 20% ethanol	Cat.# AC-129
2'/3'-EDA-ATP -Sepharose®, 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

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

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.

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.