

## Protease Inhibitor Mix (100x)

### For the purification of proteins

Cat.-No.	Amount
AK-102I-S	0.5 ml
AK-102I-L	2.5 ml

**Description:**

The 100x Protease Inhibitor Mix can be used in the purification procedure of ATP-binding proteins (for further information see data sheet of ATP Affinity Test Kit, cat.# AK-102) or for general protein purification procedures. Contains PMSF, Pefabloc, Aprotinin and Pepstatin A in methanol. Dilute 1:100 in purification buffers.

**Storage conditions:**

Store at -20°C.  
Stable for 1 year.

[Associated products available from Jena Bioscience](#)

For detailed information please view the sections on [www.jenabioscience.com](http://www.jenabioscience.com)

**FOR RESEARCH USE ONLY.  
NOT INTENDED FOR HUMAN OR ANIMAL DIAGNOSTIC OR  
THERAPEUTIC USES.**

Aminophenyl-ATP-Sepharose®, C <sub>10</sub> -linked, pre-swollen in 20% ethanol	Cat.# AC-101
8-[(6-Amino)hexyl]-amino-ATP-Sepharose®, pre-swollen in 20% ethanol	
N <sup>6</sup> -(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
200 mM Sodium Orthovanadate, activated, pH 10.0	Cat.# AK-102V
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.