

PBS Tablets

Phosphate Buffered Saline, tablets

Cat. No.	Amount
AK-102P-S	10 tablets
AK-102P-L	50 tablets

Description

Phosphate Buffered Saline – for basic applications in Molecular and Cell Biology.

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).

Usage

Dissolve a 5 g tablet in 500 ml ddH₂O to get 1x PBS (140 mM NaCl, 10 mM Phosphate Buffer, 3 mM KCl, pH 7.45).

Storage and Stability

Store at ambient temperature.
Quality guaranteed for 24 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
200 mM Sodium Orthovanadate, activated, pH 10.0	Cat.# AK-102V
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