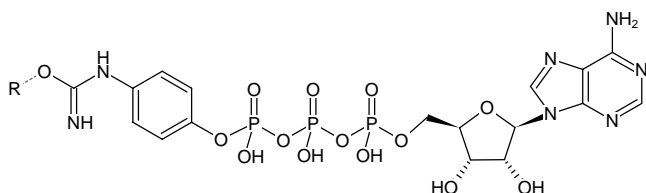




Immobilized γ -Aminophenyl-ATP (no spacer)

Aminophenyl-Adenosine triphosphate (AP-ATP) immobilized on Agarose
 γ -Aminophenyl-ATP-Agarose (no spacer)

Cat. No.	Amount
AC-102S	1 ml
AC-102L	5 ml



Structural formula of Immobilized γ -Aminophenyl-ATP (no spacer)

	Agarose characteristics
Bead/Particle size	45-165 μ m
Recommended linear flow rate	11.5 cm/h
Maximum pressure	0.25 bar (3.6 psi)
pH stability	short term: 4 - 9 / long term: 7.5
Chemical stability	Stable to all solutions commonly used in gel filtration including 8 M urea and 6 M guanidine hydrochloride Not stable in organic solvents!
Sterilization	Not autoclavable!

R= Agarose

For research use only!

Shipping: shipped at 4 °C

Storage Conditions: store at 4 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Applications:

Suitable for purification of ATP-binding proteins.

Degree of substitution: 25 μ mol - 27 μ mol AP-ATP/ml gel

Storage buffer: 20 % Ethanol

Important note: The use of this reagent for the purification of protein kinases is subject to United States Patent No. 5, 536, 822. Such use is prohibited without a license from Serenex, Inc. Information on obtaining a license is available at licensing@serenex.com.

Selected References:

Mlakar *et al.* (2006) Citrate Inhibition-Resistant Form of 6-Phosphofructo-1-Kinase from *Aspergillus niger*. *Applied and Environmental Microbiology* **72** (7):4515.

Drewes *et al.* (1995) Microtubule-associated Protein/Microtubule Affinity-regulating Kinase (p110^{mark}). *J. Biol. Chem.* **270**:7679.

Haystead *et al.* (1993) γ -phosphate linked ATP-Sepharose for the affinity purification of protein-kinases - rapid purification to homogeneity of skeletal-muscle mitogenactivated protein-kinase. *Eur. J. Biochem.* **214**:459.

Usenik *et al.* (2010) Evolution of allosteric citrate binding sites on 6-phosphofructo-1-kinase. *PLoS One.* **5** (11):e15447.