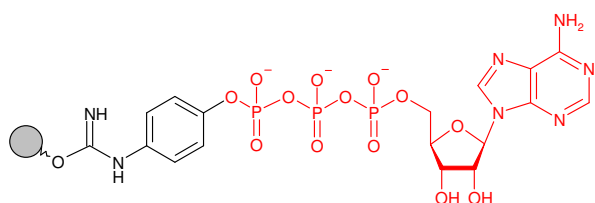


Immobilized γ -Aminophenyl-ATP (no spacer)

Cat. No.	Amount
AC-102S	1 ml (bulk material)
AC-102L	5 ml (bulk material)
AC-102C	1 ml (syringe column)
AC-102SC	0.2 ml (screening col.)
AC-102MP	1 ml (MPLC column)



	Agarose characteristics	Toyopearl AF- 650M characteristics
Bead/Particle size	45-165 μ m	40-90 μ m (mean pore size: 1000 Å)
Recommended linear flow rate	11.5 cm/h	60-600 cm/h
Maximum pressure	0.25 bar (3.6 psi)	3 bar (43.5 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!	

Aminophenyl-Adenosine triphosphate (AP-ATP) immobilized on Agarose, suitable for purification of ATP-binding proteins.

Bulk material, syringe column and screening column based on Agarose. MPLC column based on Toyopearl® AF-650M.

Toyopearl® is a trademark of TOSOH Bioscience.

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

Storage buffer: 20 % Ethanol

Storage: Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at 4°C.

Shelf life: 12 months

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.