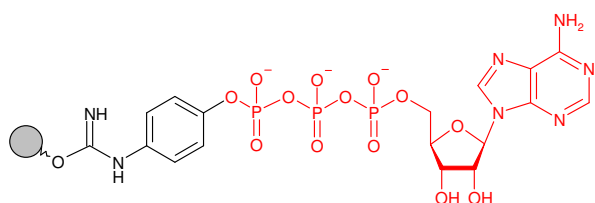


γ -Aminophenyl-ATP - Sepharose[®] (no spacer)

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 (scr. column)
AC-102MP	1 ml (MPLC column)



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

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

Linker: none

Storage buffer: 20% Ethanol

Available as prepacked column (AC-102C)

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

Long term storage at 4°C.

Expiry: 12 month

[®] Sepharose is a trademark of Amersham-Biosciences.

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:

Legisa *et al.* (2006) Citrate Inhibition-Resistant Form of δ -Phosphofructo-1-Kinase from *Aspergillus niger*. *AEM* **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.