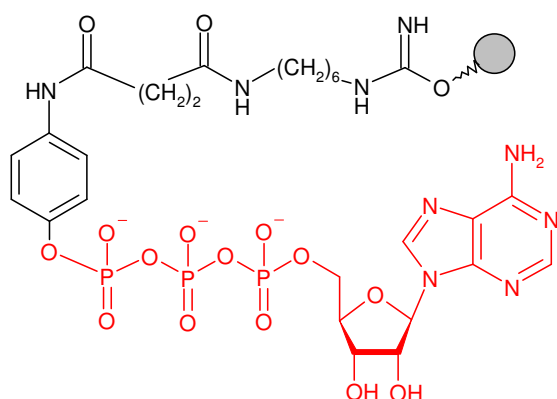


γ -Aminophenyl-ATP-Sepharose[®] (C₁₀-spacer)

Cat. No.	Amount
AC-101S	1 ml (bulk material)
AC-101L	5 ml (bulk material)
AC-101C	1 ml (syringe column)
AC-101SC	0.2 ml (scr. column)
AC-101MP	1 ml (MPLC column)



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

Degree of substitution: 20 μ mol AP-ATP/ml Gel

Linker: C₁₀

Storage buffer: 20% Ethanol

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:

- Haystead *et al.* (1993) γ -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.
- Drewes *et al.* (1995) Microtubule-associated Protein/Mic-rotubule Affinity-regulating Kinase (p110^{mark}). *J. Biol. Chem.* **270** (13): 7679.