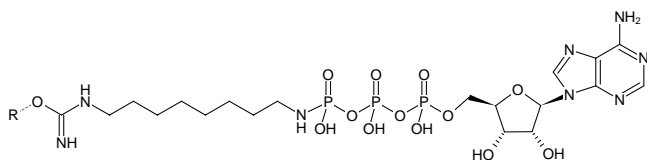




## Immobilized $\gamma$ -Amino-octyl-ATP

Adenosine triphosphate (ATP) immobilized on Agarose  
 $\gamma$ -Amino-octyl-ATP-Agarose

Cat. No.	Amount
AC-105S	1 ml
AC-105L	5 ml



Structural formula of Immobilized  $\gamma$ -Amino-octyl-ATP

	Agarose characteristics
Bead/Particle size	45-165 $\mu$ 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 <b>Not stable in organic solvents!</b>
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:** 5  $\mu$ mol ATP/ml gel

**Storage buffer:** 100 mM Tris, pH 7.5, 50 % Glycerol, 0.02 % Thimerosal

**Please note:** This compound contains a phosphoramidate linkage which is hydrolyzed at pH <7.

**Selected References:**

Suto *et al.* (1998) Synthesis of  $\gamma$ -phosphate linked nucleoside affinity chromatography resins for protein purification, including ribonucleoside triphosphate reductase. *Nucleosides and nucleotides* **17 (8)**:1453.

Tsushima *et al.* (2010) A comprehensive screening system for damaged nucleotide-binding proteins. *Mutat Res.* **703 (1)**:37.

Iyama *et al.* (2010) NUDT16 is a (deoxy)inosine diphosphatase, and its deficiency induces accumulation of single-strand breaks in nuclear DNA and growth arrest. *Nucleic Acids Res.* **38 (14)**:4834.