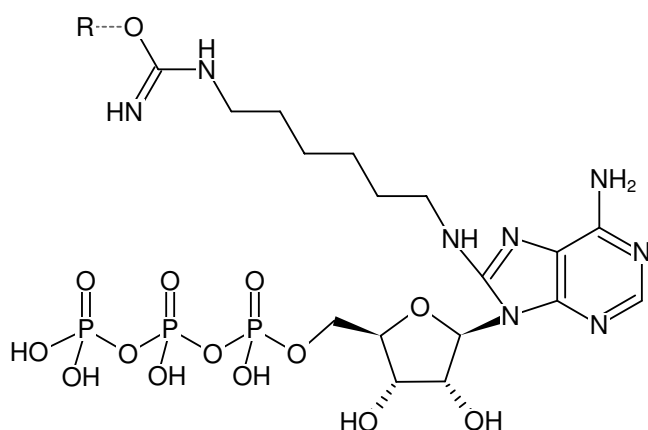




Immobilized 8-Amino-hexyl-ATP

Adenosine triphosphate (ATP) immobilized on Agarose
8-Amino-hexyl-ATP-Agarose

Cat. No.	Amount
AC-127S	1 ml
AC-127L	5 ml



Structural formula of Immobilized 8-Amino-hexyl-ATP

	Agarose characteristics
Bead/Particle size	45-165 µm
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: 5 µmol - 7 µmol ATP/ml gel

Storage buffer: 50% glycerol (cont. 0.02 % thimerosal)

Selected References:

Huang *et al.* (2003) Evolution of aptamers with a new specificity and new secondary structures from an ATP aptamer. *RNA*. **9 (12)**:1456.

Motohashi *et al.* (1994) Isolation of the stable hexameric DnaK.DnaJ complex from *Thermus thermophilus*. *J. Biol. Chem.* **269 (43)**:27074.

Welch *et al.* (1985) Rapid purification of mammalian 70,000-dalton stress proteins: affinity of the proteins for nucleotides. *Mol Cell Biol.* **5 (6)**:1229.