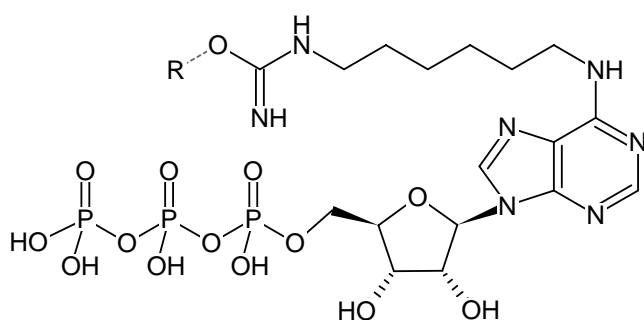




Immobilized N⁶-(6-Amino)hexyl-ATP

Adenosine triphosphate (ATP) immobilized on Agarose
N⁶-(6-Amino)hexyl-ATP-Agarose

Cat. No.	Amount
AC-129S	1 ml
AC-129L	5 ml



Structural formula of Immobilized N⁶-(6-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:

Hewitt *et al.* (2001) The human cytomegalovirus gene product US6 inhibits ATP binding by TAP. *EMBO J.* **20** (3):387.

Koyama *et al.* (1989) Phosphofructokinase from porcine heart, liver and erythrocytes. *Proc Natl Acad Sci U S A.* **93** (3):517.

Rouvière *et al.* (1985) Component A2 of the methylcoenzyme M methylreductase system from *Methanobacterium thermoautotrophicum*. *J Bacteriol.* **162** (1):61.

Robinson *et al.* (1981) Prediction of neutral salt elution profiles for affinity chromatography. *Proc Natl Acad Sci U S A.* **78** (4):2287.

Homcy *et al.* (1978) Affinity purification of cardiac adenylate cyclase: dependence on prior hydrophobic resolution. *Proc Natl Acad Sci U S A.* **75** (1):59.