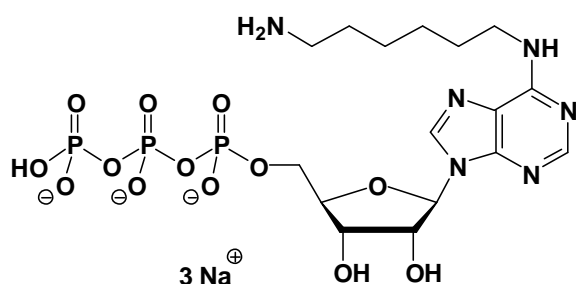


N⁶-(6-Amino)hexyl-ATP

N⁶-(6-Amino)hexyl-adenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-805S	50 Units
NU-805L	250 Units



Cat. No.: NU-805

Molecular Formula: C₁₆H₂₆N₆O₁₃P₃ (Anion)

Molecular Weight: 603.33 (Anion)

Purity: > 95%, clear aqueous solution, pH 7.5

Spectroscopic properties:

λ_{\max} 266 nm; ϵ 16200

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

For research use only!

1 unit = 1 μ l of a 10 mM solution

Selected References:

Ito *et al.* (1998) Modified nucleic acid for systematic evolution of RNA ligands by exponential enrichment. *J. Bioact. Compat. Pol.* **13** (2):114.

Trayer (1974) Affinity chromatography of some adenosine phosphate-requiring systems. *Biochem. Soc. T.* **2** (6):1302.

Trayer *et al.* (1974) Preparation of adenosine nucleotide derivatives suitable for affinity chromatography. *Biochem. J.* **139** (3):609.