

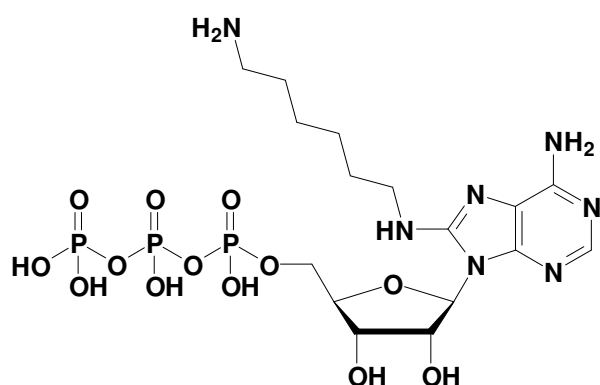
8-[(6-Amino)hexyl]-amino-ATP

8-[(6-Amino)hexyl]-amino-adenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-807S	50 Units
NU-807L	250 Units

Selected References:

Komori *et al.* (2008) Measurement system for simultaneous observation of myosin V chemical and mechanical events. *BioSystems* **93**:48.



Cat. No.: NU-807

Molecular Formula: C₁₆H₃₀N₇O₁₃P₃ (free acid)

Molecular Weight: 621.37 (free acid)

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

Spectroscopic Properties: λ_{max} 273nm;
 ε 17000

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 μl of a 10 mM solution