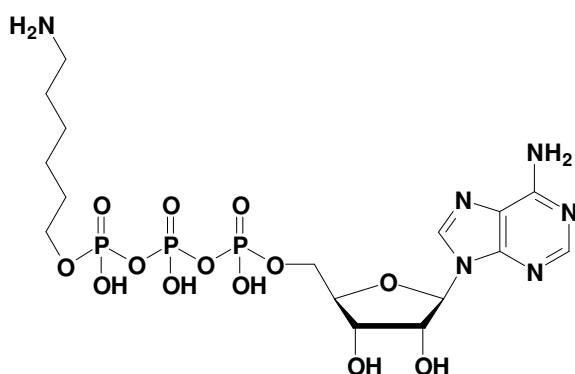


γ -[6-Aminoheptyl]-ATP

γ -[6-Aminoheptyl]-adenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-833S	50 Units
NU-833L	250 Units



Cat. No.: NU-833

Molecular Formula: $C_{16}H_{29}N_6O_{13}P_3$ (free acid)

Molecular Weight: 606.35 (free acid)

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

Spectroscopic Properties: λ_{max} 259nm;
 ϵ 15,300 $M^{-1}cm^{-1}$ Tris HCl pH 7.5

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