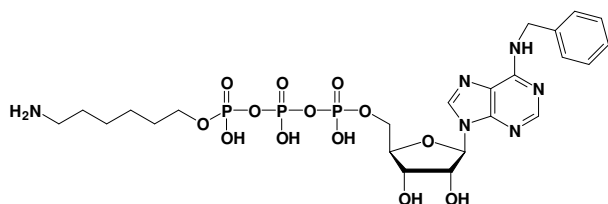


γ -[6-Aminoheptyl]-N⁶-Benzyl-ATP

γ -[6-Aminoheptyl]-N⁶-benzyl-adenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-911-1	1 mg



Cat. No.: NU-911

Molecular Formula: C₂₃H₃₅N₆O₁₃P₃ (free acid)

Molecular Weight: 696.48 (free acid)

Purity: > 95%, lyophilized solid

Spectroscopic Properties: λ_{\max} 269nm;
 ϵ 20,500 M⁻¹cm⁻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!