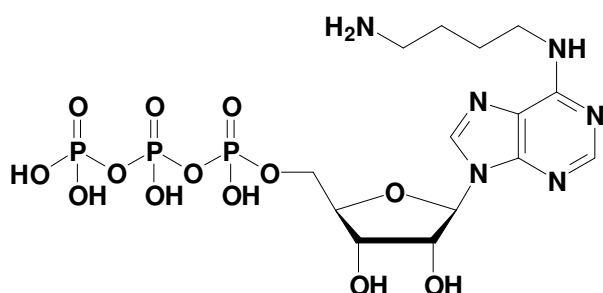


N⁶-(4-Amino)butyl-ATP

N⁶-(4-Amino)butyl-adenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-804S	50 Units
NU-804L	250 Units



Cat. No.: NU-804

Molecular Formula: C₁₄H₂₅N₆O₁₃P₃ (free acid)

Molecular Weight: 578.30 (free acid)

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

Spectroscopic properties:

λ_{\max} 266 nm; ϵ 16,200 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!

1 unit = 1 μ l of a 10 mM solution