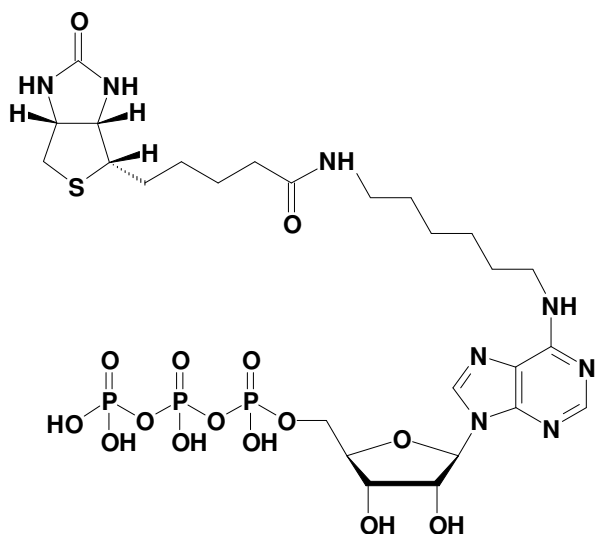


N6-(6-Amino)hexyl-ATP - Biotin

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

Cat. No.	Amount
NU-805-BIO	50 μ l / 5 mM



Applications:

RNA-labeling^[1]

Isolation and characterization of ligase ribozyme^[2]

Selected References:

[1] Folsom *et al.* (1989) Detection of DNA targets with biotinylated and fluoresceinated RNA probes. Effect of the extent of derivatization on detection sensitivity. *Analytic. Biochem.* **182**:309.

[2] Teramoto *et al.* (2000) In vitro selection of a ligase ribozyme carrying alkylamino groups in the side chains. *Bioconjugate Chem.* **11**:744.

Cat. No.: NU-805-BIO

Molecular Formula: C₂₆H₄₃N₈O₁₅P₃S (free acid)

Molecular Weight: 832.65 (free acid)

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

Spectroscopic Properties: λ_{\max} 266nm;
 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!