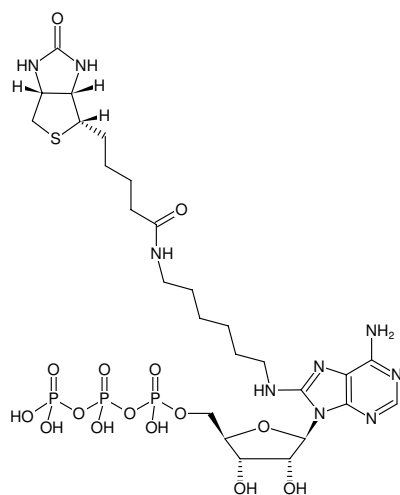




8-(6-Aminoheptyl)-amino-ATP-Biotin

8-(6-Aminoheptyl)-amino-adenosine-5'-triphosphate-Biotin, Triethylammonium salt

Cat. No.	Amount
NU-807-BIO	50 µl (5 mM)



Structural formula of 8-(6-Aminoheptyl)-amino-ATP-Biotin

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: C₂₆H₄₄N₉O₁₅P₃S (free acid)

Molecular Weight: 847.66 g/mol (free acid)

Exact Mass: 847.19 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 5.0 mM - 5.5 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{max} 273 nm, ε 17.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)