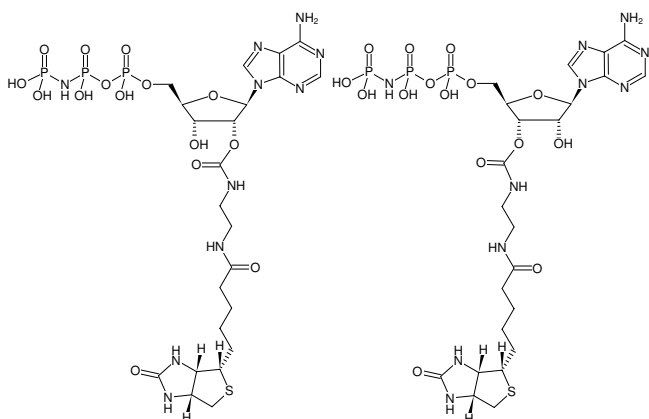




EDA-AppNHp (EDA-AMPPNP)-Biotin

2'/3'-O-(2-Aminoethyl-carbamoyl)-adenosine-5'-[(β,γ)-imido] triphosphate-Biotin, Triethylammonium salt

Cat. No.	Amount
NU-810-BIO	100 µl (5 mM)



Structural formula of EDA-AppNHp (EDA-AMPPNP)-Biotin

For general laboratory use.

Shipping: shipped on dry ice

Storage Conditions: store at -20 °C

Shelf Life: 6 months after date of delivery

Molecular Formula: C₂₃H₃₇N₁₀O₁₅P₃S (free acid)

Molecular Weight: 818.58 g/mol (free acid)

Exact Mass: 818.14 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} 259 nm, ε 15.3 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)