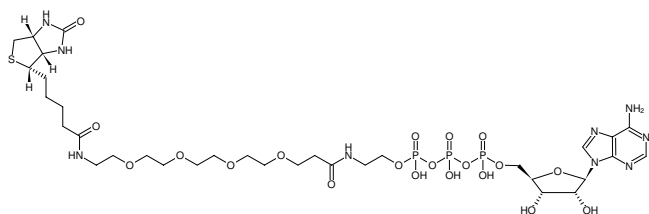




γ -[Ethyl-CONH-(PEG4)]-ATP-Biotin

γ -[Ethyl-CONH-(PEG4)]-Adenosine-5'-triphosphate-Biotin, Triethylammonium salt

Cat. No.	Amount
NU-927-BIO	50 μ l (5 mM)



Structural formula of γ -[Ethyl-CONH-(PEG4)]-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: C₃₃H₅₆N₉O₂₀P₃S (free acid)

Molecular Weight: 1023.83 g/mol (free acid)

Exact Mass: 1023.26 g/mol (free acid)

Purity: \geq 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 5.0 mM - 5.5 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{max} 259 nm, ϵ 15.3 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)