



γ -(6-Aminoethyl)-ATP-Biotin

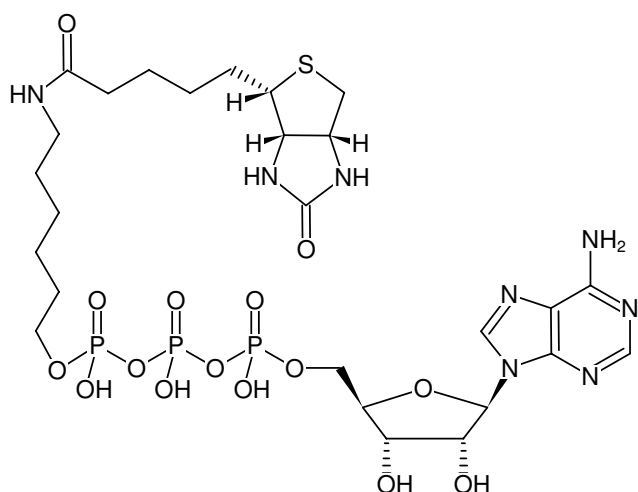
γ -(6-Aminoethyl)-adenosine-5'-triphosphate-Biotin, Triethylammonium salt

Cat. No.	Amount
NU-833-BIO	50 μ l (5 mM)

Selected References:

Tian *et al.* (2019) Ultrasensitive amplification-free detection of protein kinase based on catalyzed assembly and enumeration of gold nanoparticles. *Chem. Commun. (Camb)*. **55 (17)**:2505.

Wang *et al.* (2015) Phosphorylation-directed assembly of a single quantum dot based nanosensor for protein kinase assay. *Anal. Chem.* **87 (9)**:4696.



Structural formula of γ -(6-Aminoethyl)-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: C₂₆H₄₃N₈O₁₅P₃S (free acid)

Molecular Weight: 832.65 g/mol (free acid)

Exact Mass: 832.18 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)