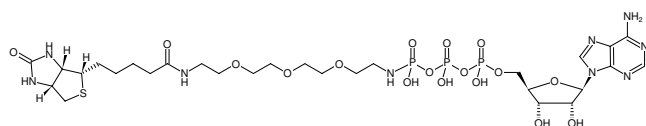




γ -[(PEG₃-Amino)-imido]-ATP - Biotin

Adenosine-5'-triphosphate [γ]-biotinyl-3,6,9-trioxaundecanediamine; ATP- γ -Biotin-LC-PEO-amine
 γ -[(PEG₃-Amino)-imido]-adenosine-5'-triphosphate - Biotin, Sodium salt

Cat. No.	Amount
NU-970-BIO	50 μ l (1 mM)



Structural formula of γ -[(PEG₃-Amino)-imido]-ATP - Biotin

For research use only!

Shipping: shipped on blue ice

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: C₂₈H₄₈N₉O₁₇P₃S (free acid)

Molecular Weight: 907.72 g/mol (free acid)

Purity: \geq 95 % (HPLC)

Form: clear aqueous solution

Concentration: 1.0 mM - 1.1 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)

Please note: This compound contains a phosphoramidate linkage which is hydrolyzed at pH <7.

Selected References:

Arora *et al.* (2017) Unexpected biotinylation using ATP-c-Biotin-LC-PEO-amine as a kinase substrate. *Biochem. Biophys. Res. Commun.* **432**:287.

Embogama *et al.* (2017) K-BILDS: A Kinase Substrate Discovery Tool. *Chembiochem.* **18** (1):136.

Ma *et al.* (2010) Highly sensitive detection of DNA phosphorylation by counting single nanoparticles. *Anal. Bioanal. Chem.* **397** (6):2279.

Seneviranthne *et al.* (2012) Kinase-Catalyzed Biotinylation. *Curr. Protoc. Chem. Biol.* **4** (1):83.

Seneviranthne *et al.* (2016) The generality of kinase-catalyzed biotinylation. *Biorg. Med. Chem.* **24** (1):12.

Green *et al.* (2007) Kinase-catalyzed biotinylation for phosphoprotein detection. *J. Am. Chem. Soc.* **129** (1):10.