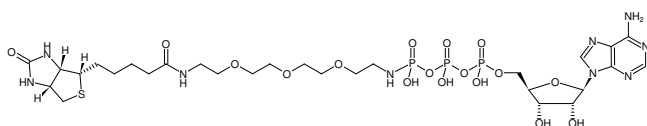




## $\gamma$ -[(PEG<sub>3</sub>-Amino)-imido]-ATP-Biotin

Adenosine-5'-triphosphate [ $\gamma$ ]-biotinyl-3,6,9-trioxaundecanediamine; ATP- $\gamma$ -Biotin-LC-PEO-amine  
 $\gamma$ -[(PEG<sub>3</sub>-Amino)-imido]-adenosine-5'-triphosphate-Biotin, Sodium salt

Cat. No.	Amount
NU-970-BIO	50 $\mu$ l (1 mM)



Structural formula of  $\gamma$ -[(PEG<sub>3</sub>-Amino)-imido]-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:** C<sub>28</sub>H<sub>48</sub>N<sub>9</sub>O<sub>17</sub>P<sub>3</sub>S (free acid)

**Molecular Weight:** 907.72 g/mol (free acid)

**Exact Mass:** 907.21 g/mol (free acid)

**Purity:**  $\geq$  95 % (HPLC)

**Form:** solution in water

**Color:** colorless to slightly yellow

**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5  $\pm$  0.5

**Spectroscopic Properties:**  $\lambda_{\max}$  259 nm,  $\epsilon$  15.3 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

**Please note:** This compound contains a phosphoramidate linkage which is hydrolyzed at pH <7.0.

### 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.