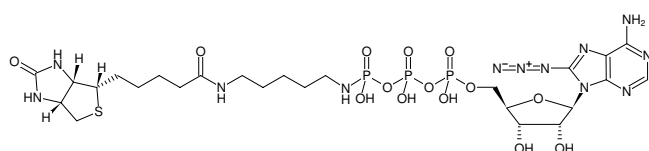




8-Azido-ATP- γ -Biotin

8-N³-ATP[γ]biotinpentylamine, Sodium salt

Cat. No.	Amount
NU-252-BIO	1 mg



Structural formula of 8-Azido-ATP- γ -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: C₂₅H₄₁N₁₂O₁₄P₃S (free acid)

Molecular Weight: 858.65 g/mol (free acid)

Exact Mass: 858.18 g/mol (free acid)

Purity: \geq 95 % (HPLC)

Form: solid

Color: white to off-white

Spectroscopic Properties: λ_{\max} 281 nm, ϵ 13.3 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:

Qui *et al.* (2015) ATP Binding and Hydrolysis Properties of ABCB10 and Their Regulation by Glutathione. *PLoS One*. doi:10.1371.

Mo *et al.* (2013) Assessing ATP binding and hydrolysis by NLR proteins. *Methods Mol. Biol.* **1040**:153.

Yao *et al.* (2009) SVOP is a nucleotide binding protein. *PLoS One*. **4** (4):e5315.

Yao *et al.* (2008) Synaptic Vesicle Protein 2 Binds Adenine Nucleotides. *J. Biol. Chem.* **283** (30):20628.