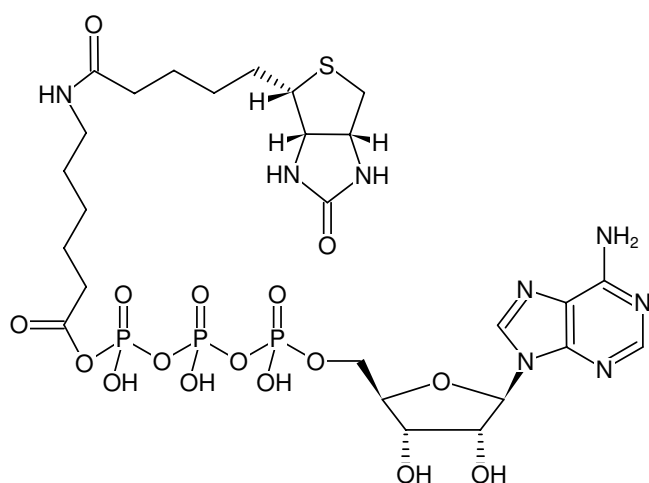




ATP-acetyl-hex-Biotin

Biotin-ATP probe
 Biotin-hex-acyl-ATP (BHAcATP)
 Tri(triethylammonium) salt

Cat. No.	Amount
NU-277	16 x 0.01 μ mol (16 x approximately 11.5 μ g)



Structural formula of ATP-acetyl-hex-Biotin

For research use only!

Shipping: shipped on dry ice

Storage Conditions: store at -80 °C

Shelf Life: 6 months after date of delivery

Molecular Formula: C₂₆H₄₁N₈O₁₆P₃ (free acid)

Molecular Weight: 846.64 g/mol (free acid)

Exact Mass: 846.16 g/mol (free acid)

Purity: \geq 75 % (HPLC)

Form: solid

Color: white to off-white

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

Please note:

- Minimize exposure of the product to ambient temperature and store unused product at -80 °C.
- Avoid repeated freeze/thaw cycles.
- Do not prepare stock solutions for storage.
- Equilibrate vial to room temperature before opening and use it immediately.
- The product is typically not visible in the vial.

Selected References:

Villamor *et al.* (2013) Profiling protein kinases and other ATP binding proteins in Arabidopsis using Acyl-ATP probes. *Mol. Cell. Proteomics* **12** (9):2481.

Xiao *et al.* (2013) Proteome-wide discovery and characterizations of nucleotide-binding proteins with affinity-labeled chemical probes. *Anal. Chem.* **85** (6):3198.

Patricelli *et al.* (2007) Functional interrogation of the kinome using nucleotide acyl phosphates. *Biochemistry* **46** (2):350.