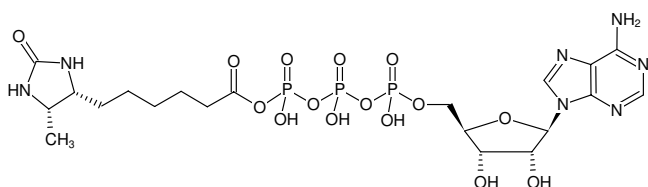




ATP-acetyl-Desthiobiotin

Desthiobiotin-ATP probe
 Desthiobiotin-acyl-ATP (DBAcATP)
 Tri(triethylammonium) salt

| Cat. No. | Amount |
|----------|---|
| NU-276 | 16 x 0.01 μ mol (16 x approximately 10.1 μ g) |



Structural formula of ATP-acetyl-Desthiobiotin

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: C₂₀H₃₂N₇O₁₅P₃ (free acid)

Molecular Weight: 703.43 g/mol (free acid)

Exact Mass: 703.12 g/mol (free acid)

Purity: \geq 75 % (HPLC)

Form: lyophilised

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:

Okerberg *et al.* (2019) Chemoproteomics Using Nucleotide Acyl Phosphates Reveals an ATP Binding Site at the Dimer Interface of Procaspase-6. *Biochemistry*. doi:10.1021/acs.biochem.9b00290.

Nordin *et al.* (2015) ATP Acyl Phosphate Reactivity Reveals Native Conformations of Hsp90 Paralogs and Inhibitor Target Engagement. *Biochemistry*. **54** (19):3024.

Adachi *et al.* (2014) Proteome-wide discovery of unknown ATP-binding proteins and kinase inhibitor target proteins using an ATP probe. *J. Proteome Res.* **13** (12):5461.

Edgar *et al.* (2014) A combination of SILAC and nucleotide acyl phosphate labelling reveals unexpected targets of the Rsk inhibitor BI-D1870. *Biosci. Rep.* **34** (1):e00091.