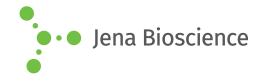
# **DATA SHEET**





# **ATP-acetyl-hex-Biotin**

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

Cat. No.	Amount
NU-277	16 x 0.01 μmol (16 x approximately 11.5 μg)

Structural formula of ATP-acetyl-hex-Biotin

## For general laboratory use.

**Shipping:** shipped on dry ice

Storage Conditions: store at -80 °C

**Shelf Life:** 6 months after date of delivery **Molecular Formula:** C<sub>26</sub>H<sub>41</sub>N<sub>8</sub>O<sub>16</sub>P<sub>3</sub> (free acid) **Molecular Weight:** 846.64 g/mol (free acid)

Exact Mass: 846.16 g/mol (free acid)

**Purity:** ≥ 75 % (HPLC)

Form: solid

Color: colorless to white

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:

- 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.