



N⁶-(6-Aminohexyl)-ATP-Biotin

N⁶-(6-Aminohexyl)-adenosine-5'-triphosphate-Biotin, Triethylammonium salt

Cat. No.	Amount
NU-805-BIO	50 µl (5 mM)

Applications:

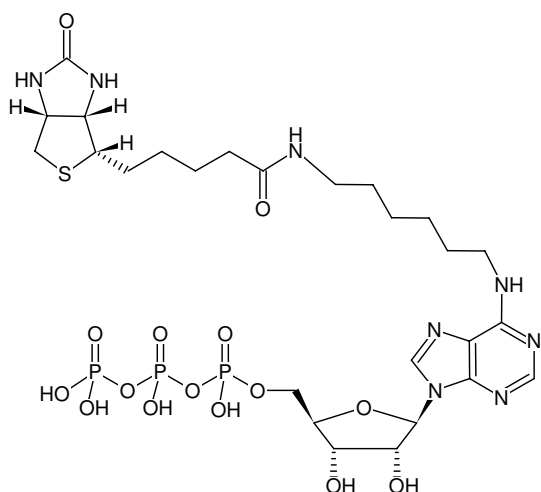
Incorporation into RNA by

- 3'-end labeling with yeast poly(A) polymerase^[1], *S. pombe* Cid1^[2] poly(A) polymerase and *E. coli* poly(A) polymerase^[2]

Selected References:

[1] Martin *et al.* (1998) Tailing and 3'-end labeling of RNA with yeast poly (A) polymerase and various nucleotides. *RNA*. **4** (2):226.

[2] Moritz *et al.* (2014) Simple methods for the 3' biotinylation of RNA. *RNA*. **20** (3):421.



Structural formula of N⁶-(6-Aminohexyl)-ATP-Biotin

For research use only!

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: C₂₆H₄₃N₈O₁₅P₃S (free acid)

Molecular Weight: 832.65 g/mol (free acid)

Exact Mass: 832.18 g/mol (free acid)

CAS#: 126320-25-0

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 5.0 mM - 5.5 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{max} 266 nm, ε 16.2 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)