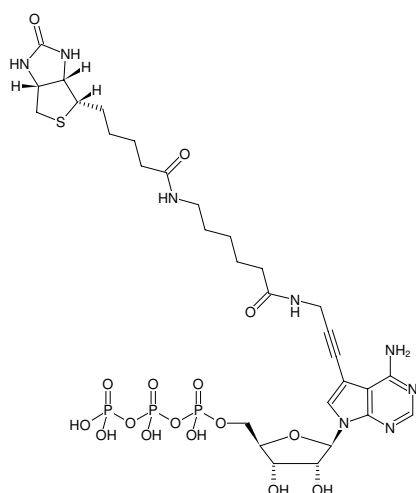




Biotin-11-ATP

Triethylammonium salt

Cat. No.	Amount
NU-957-BIOX-S	10 µl (1 mM)
NU-957-BIOX-L	5 x 10 µl (1 mM)



Structural formula of Biotin-11-ATP

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: C₃₀H₄₅N₈O₁₆P₃S (free acid)

Molecular Weight: 898.71 g/mol (free acid)

Exact Mass: 898.19 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: filtered solution (30 kDa) in 10 mM Tris-HCl

Color: colorless to slightly yellow

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{max} 280 nm, ε 12.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Selected References:

Chu *et al.* (2020) Rocaglates Induce Gain-of-Function Alterations to eIF4A and eIF4F. *Cell Rep.* **30** (8):2481.

Parida *et al.* (2019) Nucleotide Resolution Comparison of Transcription of Human Cytomegalovirus and Host Genomes Reveals Universal Use of RNA Polymerase II Elongation Control Driven by Dissimilar Core Promoter Elements. *mBio* **10** (1):e02047-18.

Mahat *et al.* (2016) Base-pair-resolution genome-wide mapping of active RNA polymerases using precision nuclear run-on (PRO-seq). *Nat. Protoc.* **11** (8):1455.

Nagarkatti *et al.* (2012) Development of an aptamer-based concentration method for the detection of *Trypanosoma cruzi* in blood. *PLoS One* **7** (8):e43533.

Mitobe *et al.* (2008) Involvement of RNA-binding protein Hfq in the post-transcriptional regulation of *invE* gene expression in *Shigella sonnei*. *Int. J. Parasitol.* **283** (9):5738.

Göringer *et al.* (2003) In vitro selection of high-affinity nucleic acid ligands to parasite target molecules. *Int. J. Parasitol.* **33** (12):1309.