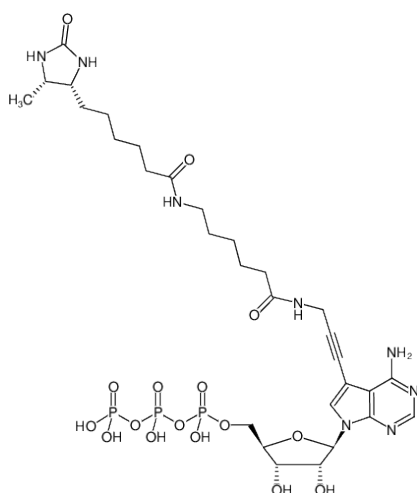




Desthiobiotin-11-ATP

7-Deaza-7-propargylamino-adenosine-5'-triphosphate, labeled with Desthiobiotin, Triethylammonium salt

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



Structural formula of Desthiobiotin-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: C₃₀H₄₇N₈O₁₆P₃ (free acid)

Molecular Weight: 868.67 g/mol (free acid)

Exact Mass: 868.23 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)