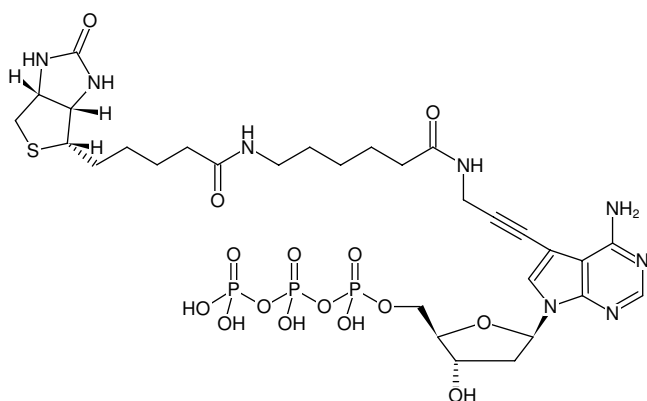



Biotin-11-dATP - high concentration

γ -[N-(Biotin-6-amino-hexanoyl)]-7-propargylamino-2'-deoxy-7-deaza-adenosine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-1175-BIOX-HC	20 μ l (5 mM)



Structural formula of Biotin-11-dATP - high concentration

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: C₃₀H₄₅N₈O₁₅P₃S (free acid)

Molecular Weight: 882.71 g/mol (free acid)

Exact Mass: 882.19 g/mol (free acid)

Purity: \geq 95 % (HPLC)

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

Color: colorless to slightly yellow

Concentration: 5.0 mM - 5.5 mM

pH: 7.5 \pm 0.5

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