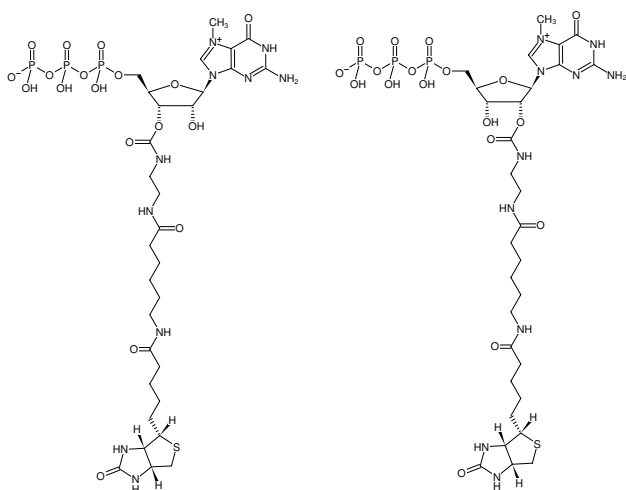



Biotin-11-EDA-m⁷GTP

Biotin-11-2'/3'-O-(2-aminoethyl-carbamoyl)-N7-methyl-guanosine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-824-BIOX	250 µl (1 mM)


 Structural formula of Biotin-11-EDA-m⁷GTP

For research use only!
Shipping: shipped on blue ice

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: C₃₀H₄₉N₁₀O₁₈P₃S (free acid)

Molecular Weight: 962.75 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: clear aqueous solution

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{max} 258/280 nm; ε 9.8/8.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)