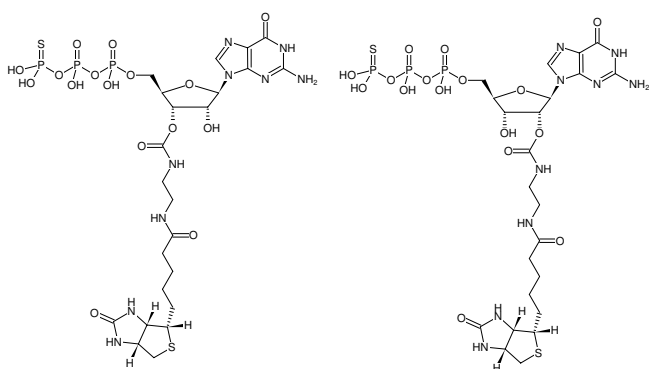




EDA-GTPyS-Biotin

2'/3'-O-(2-Aminoethyl-carbamoyl)-guanosine-5'-(γ-thio)-triphosphate-Biotin, Triethylammonium salt

Cat. No.	Amount
NU-1610-BIO	100 µl (1 mM)



Structural formula of EDA-GTPyS-Biotin

For general laboratory use.

Shipping: shipped on dry ice

Storage Conditions: store at -20 °C

Shelf Life: 6 months after date of delivery

Molecular Formula: C₂₃H₃₆N₉O₁₆P₃S₂ (free acid)

Molecular Weight: 851.63 g/mol (free acid)

Exact Mass: 851.09 g/mol (free acid)

Purity: ≥ 90 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 1.0 mM - 1.1 mM

pH: > 7.5

Spectroscopic Properties: λ_{max} 252 nm, ε 13.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)