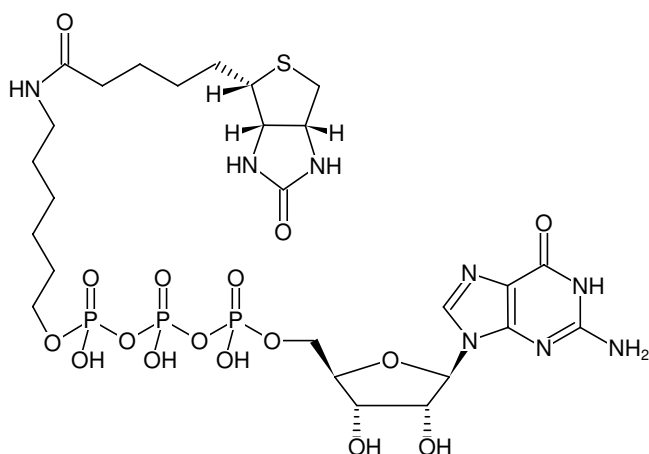




γ -(6-Aminoethyl)-GTP-Biotin

γ -(6-Aminoethyl)-guanosine-5'-triphosphate-Biotin, Triethylammonium salt

Cat. No.	Amount
NU-834-BIO	50 μ l (5 mM)



Structural formula of γ -(6-Aminoethyl)-GTP-Biotin

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: C₂₆H₄₃N₈O₁₆P₃S (free acid)

Molecular Weight: 848.65 g/mol (free acid)

Exact Mass: 848.17 g/mol (free acid)

Purity: \geq 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 5.0 mM - 5.5 mM

pH: 7.5 \pm 0.5

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