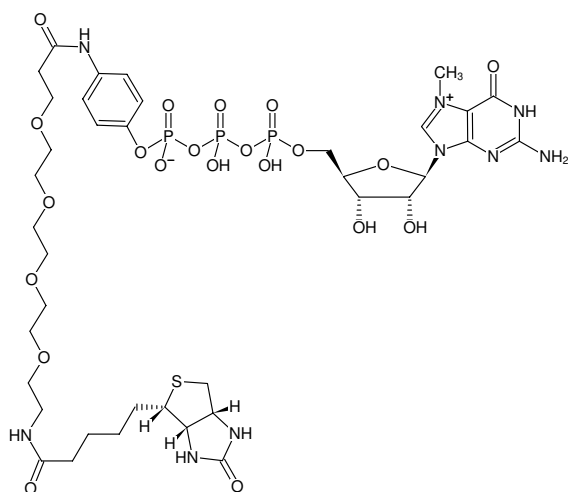



m⁷GTP-γ-aminophenyl-PEG₄-Biotin

 7-Methylguanosine-5'-[γ-(4-aminophenyl)-PEG₄]-5'-triphosphate-Biotin, Sodium salt

Cat. No.	Amount
NU-869-BIO	50 µl (5 mM)


 Structural formula of m⁷GTP-γ-aminophenyl-PEG₄-Biotin

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: C₃₈H₅₈N₉O₂₁P₃S (free acid)

Molecular Weight: 1101.91 g/mol (free acid)

Exact Mass: 1101.27 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 5.0 mM - 5.5 mM

pH: 7.0 ±0.5

Spectroscopic Properties: λ_{max} 260 nm, ε 10.4 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)