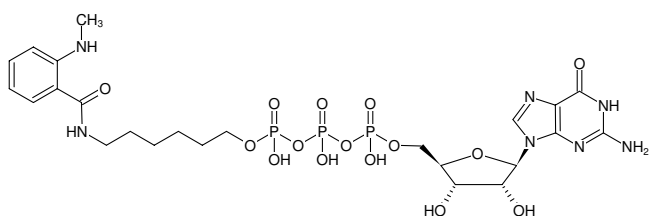




γ -(6-Aminoethyl)-GTP-MANT

γ -(6-Aminoethyl)-guanosine-5'-triphosphate, labeled with MANT, Triethylammonium salt

Cat. No.	Amount
NU-834-MNT	200 μ l (5 mM)



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

For research use only!

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: C₂₄H₃₆N₇O₁₅P₃ (free acid)

Molecular Weight: 755.50 g/mol (free acid)

Exact Mass: 755.15 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_{exc} 335 nm, λ_{em} 440 nm, ϵ 2.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)