

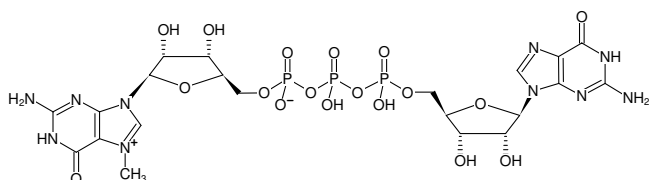

m⁷GP₃G (Monomethylated Cap Analog) - Solution

 m⁷G(5')ppp(5')G

 P¹-(5'-7-methyl-guanosyl) P³-(5'-(guanosyl))triphosphate, Sodium salt

Cat. No.	Amount
NU-852S	10 µl (100 mM)
NU-852L	5 x 10 µl (100 mM)

Selected References:

 Krieg *et al.* (1987) In vitro RNA synthesis with SP6 RNA polymerase. *Methods Enzymol.* **155**:397.

 Structural formula of m⁷GP₃G (Monomethylated Cap Analog) - Solution

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: C₂₁H₂₉N₁₀O₁₈P₃ (free acid)

Molecular Weight: 802.44 g/mol (free acid)

Exact Mass: 802.09 g/mol (free acid)

CAS#: 62828-64-2 (acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 100 mM - 110 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{max} 254 nm, λ 260 nm, ε 21.6 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)