

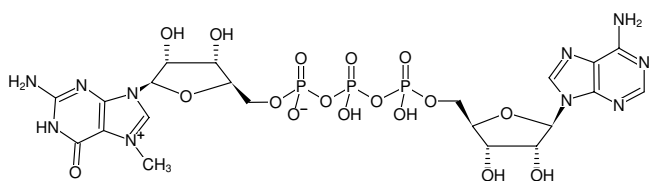
**m⁷GP₃A**m⁷GpppA, m⁷G(5')ppp(5')AP¹-(5'-7-methyl-guanosyl) P³-(5'-(adenosyl))triphosphate, Sodium salt

Cat. No.	Amount
NU-535S	10 µl (100 mM)
NU-535L	5 x 10 µl (100 mM)

Selected References:

Tomoo *et al.* (2003) Structural features of human initiation factor 4E, studied by X-ray crystal analyses and molecular dynamics simulations. *J Mol Biol.* **328(2)**:365-83.

Contreras *et al.* (1982) Simple, efficient in vitro synthesis of capped RNA useful for direct expression of cloned eukaryotic genes. *Nucleic Acids Res.* **10(20)**:6353-62.

Structural formula of m⁷GP₃A**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:** C₂₁H₂₉N₁₀O₁₇P₃ (free acid)**Molecular Weight:** 786.44 g/mol (free acid)**Exact Mass:** 786.09 g/mol (free acid)**CAS#:** 62828-63-1 (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} 260 nm, ε 22.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)