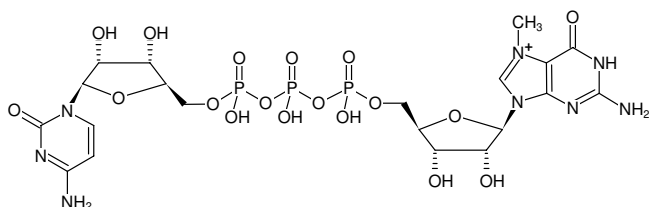


**m⁷GP₃C**P¹-(5'-7-methyl-guanosyl) P³-(5'-(cytidyl))triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-512S	50 µl (10 mM)
NU-512L	5 x 50 µl (10 mM)

Selected References:

Guranowski *et al.* (1996) Dinucleoside 5', 5''-P₁, P₃-triphosphate hydrolase from yellow lupin (*Lupinus luteus*) seeds: purification to homogeneity and hydrolysis of mRNA 5'-cap analogs. *Protein Expr Purif.* 1996 **8** (4):416.

Structural formula of m⁷GP₃C**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:** C₂₀H₂₉N₈O₁₈P₃ (free acid)**Molecular Weight:** 762.41 g/mol (free acid)**Exact Mass:** 762.08 g/mol (free acid)**CAS#:** 70441-86-0**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 260 nm, ε 17.3 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)