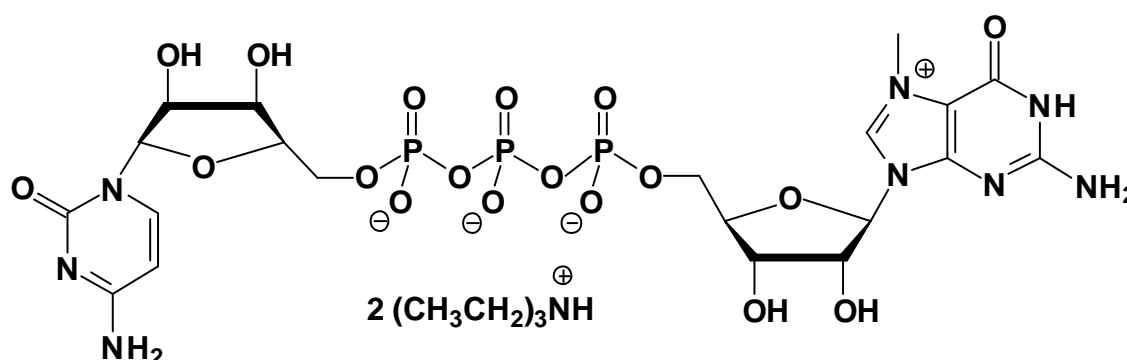


m⁷GP₃C

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

Cat. No.	Amount
NU-512S	50 Units
NU-512L	250 Units



Catalogue-No.: NU-512

Molecular Formula: C₂₀H₂₇N₈O₁₈P₃ (Anion)

Molecular Weight: 760.39 (Anion)

Purity: > 95%, clear aqueous solution, pH 7.5

Spectroscopic properties: λ_{max} 260 nm
ε 17300

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

For research use only!

1 unit = 1 μl of a 10 mM solution

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.