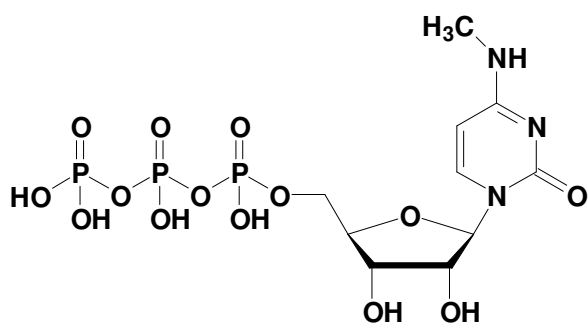


4-N-Methyl-CTP 4-N-Methyl-cytidine-5'-triphosphate

Cat. No.	Amount
NU-1192S	20 Units
NU-1192L	100 Units



Cat. No.: NU-1192

Molecular Formula: C₁₀H₁₈N₃O₁₄P₃ (free acid)

Molecular Weight: 497.18 (free acid)

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

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.

* 1 unit = 1 µl of a 10 mM solution

For research use only!