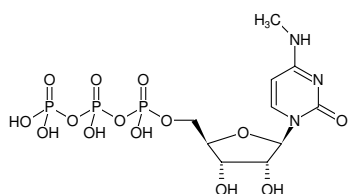


**N⁴-Methyl-CTP**N⁴-Methyl-cytidine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1192S	20 µl (10 mM)
NU-1192L	5 x 20 µl (10 mM)

Selected References:

Miao *et al.* (2015) Base-resolution detection of N⁴-methylcytosine in genomic DNA using 4mC-Tet-assisted-bisulfite-sequencing. *Nucleic Acids Res.* **43** (21):148.

Structural formula of N⁴-Methyl-CTP**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:** C₁₀H₁₈N₃O₁₄P₃ (free acid)**Molecular Weight:** 497.18 g/mol (free acid)**Exact Mass:** 497.00 g/mol (free acid)**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} 274 nm, ε 11.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)