DATA SHEET





5-Methyl-CTP

m⁵CTP 5-Me-CTP 5-Methylcytidine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1138S	10 μl (100 mM)
NU-1138L	5 x 10 μl (100 mM)



Structural formula of 5-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)

CAS#: 327174-86-7 (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} 277 nm, ϵ 9.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Applications:

Functional influence of modified bases^[1-5]

Incorporation by PCR^[3]

Related Products:

HighYield T7 RNA Synthesis Kit, #RNT-101

Selected References:

[1] Adams et al. (2014) 185 CELLULAR VIABILITY AND EXPRESSION OF GREEN FLUORESCENT PROTEIN IN BOVINE FETAL FIBROBLASTS FOLLOWING TRANSFECTION OF SYNTHETIC mRNA INCLUDING MODIFIED BASES. *Reprod. Fertil. Dev.* **26**:207.

[2] Tchigvintsev *et al.* (2013) Biochemical and Structural Studies of Conserved Maf Proteins Revealed Nucleotide Pyrophosphatases with a Preference for Modified Nucleotides. *Chem. Biol. Oxford U.K.* **20**:1386.

[3] Wong *et al.* (1997) A novel method for producing partial restriction digestion of DNA fragments by PCR with 5-methyl-CTP. *Nucleic Acids Res.* **25**:4169.

[4] Kormann *et al.* (2011) Expression of therapeutic proteins after delivery of chemically modified mRNA in mice. *Nature Biotechnology* **29**:154.

[5] Warren *et al.* (2010) Highly efficient reprogramming to pluripotency and directed differentiation of human cells with synthetic modified mRNA. *Cell Stem Cell* **7** (5):618.

