



N4-Methyl-dCTP

N4-Methyl-2'-deoxycytidine-5'-triphosphate, Sodium salt

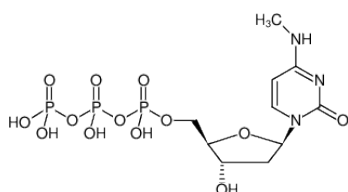
Cat. No.	Amount
NU-219S	50 µl (100 mM)
NU-219L	5 x 50 µl (100 mM)

Related Products:

Selected References:

Flores-Juarez *et al.* (2013) Capacity of N4-methyl-2'-deoxycytidine 5'-triphosphate to sustain the polymerase chain reaction using various thermostable DNA polymerases. *Analytical Biochemistry* **438**(1):73.

Li *et al.* (1993) Elimination of band compression in sequencing gels by the use of N4-methyl-2'-deoxycytidine 5' triphosphate. *Analytical Biochemistry* **21**(11):2709.



Structural formula of N4-Methyl-dCTP

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: C₁₀H₁₈N₃O₁₃P₃ (free acid)

Molecular Weight: 481.18 g/mol (free acid)

Exact Mass: 481.01 g/mol (free 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} 270 nm, ε 11.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)