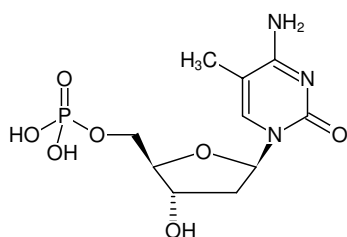




5-Methyl-dCMP

5-Methyl-2'-deoxycytidine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-1137S	100 µl (10 mM)
NU-1137L	5 x 100 µl (10 mM)



Structural formula of 5-Methyl-dCMP

For research use only!

Shipping: shipped on blue ice

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: C₁₀H₁₆N₃O₇P (free acid)

Molecular Weight: 321.23 g/mol (free acid)

Exact Mass: 321.07 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: clear aqueous solution

Concentration: 10 mM - 11 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)

Selected References:

Kamiya *et al.* (2004) Important amino acids in the phosphohydrolase module of *Escherichia coli* Orf135. *Biochem Biophys Res Commun.* **323** (3):1063.

Holliday *et al.* (2002) DNA methylation and epigenetic inheritance. *Methods.* **27** (2):179.

Kaito *et al.* (2001) Activation of the maternally preset program of apoptosis by microinjection of 5-aza-2'-deoxycytidine and 5-methyl-2'-deoxycytidine-5'-triphosphate in *Xenopus laevis* embryos. *Dev Growth Differ.* **43** (4):383.

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** (20):4169.

Chen *et al.* (1993) Direct induction of DNA hypermethylation in sea urchin embryos by microinjection of 5-methyl dCTP stimulates early histone gene expression and leads to developmental arrest. *Dev Biol.* **155** (1):75.

Gehrke *et al.* (1984) Quantitative reversed-phase high-performance liquid chromatography of major and modified nucleosides in DNA. *J. Chromatogr.* **301**:199.

Eick *et al.* (1983) Quantitative determination of 5-methylcytosine in DNA by reverse-phase high-performance liquid chromatography. *Anal. Biochem.* **135**:165.