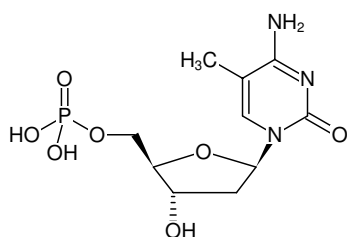


**5-Methyl-dCMP**

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 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:** C₁₀H₁₆N₃O₇P (free acid)**Molecular Weight:** 321.23 g/mol (free acid)**Exact Mass:** 321.07 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} 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.