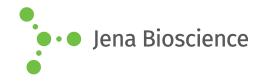
DATA SHEET





■ 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 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:** $C_{10}H_{16}N_3O_7P$ (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.