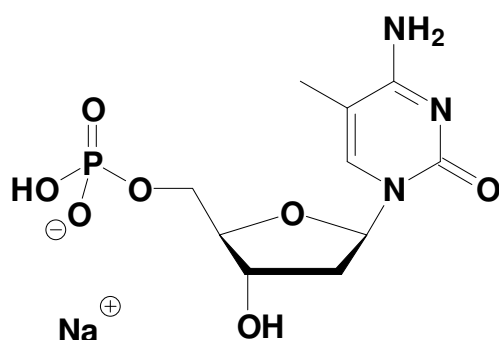


## 5-Methyl-dCMP

5-Methyl-2'-Deoxy-Cytidine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-1137S	100 Units
NU-1137L	500 Units



**Cat. No.:** NU-1137

**Molecular Formula:** C<sub>10</sub>H<sub>15</sub>N<sub>3</sub>O<sub>7</sub>P (Anion)

**Molecular Weight:** 320.21 (Anion)

**Purity:** > 95%, clear aqueous solution, pH 7.5

### Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

\* 1 unit = 1 µl of a 10 mM solution

**For research use only!**

### 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.