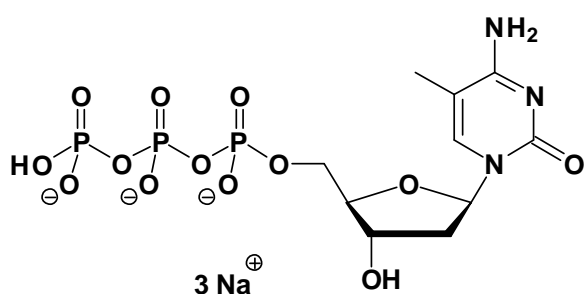


5-Methyl-dCTP

5-Methyl-2'-deoxy-cytidine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1125S	100 Units
NU-1125L	500 Units



Cat. No.: NU-1125

Molecular Formula: C₁₀H₁₅N₃O₁₃P₃ (Anion)

Molecular Weight: 478.16 (Anion)

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

Spectroscopic properties: λ_{max} 277 nm;
ε 9000

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C.

For research use only!

1 unit = 1 μl of a 10 mM solution

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.

Ito, A. *et al.* (1998) *Blood* **91-9**:3210

Lefaucheur, L. *et al.* (1998) *J. Anim. Sci.* **76**:1584

Nelson, P. S. *et al.* (1993) *Nucl. Acid. Res.* **21-3**:681

Asamizu, E. *et al.* (1999) *DNA Research* **6**:369

Wong, K. K. *et al.* (1991) *Nucl. Acid. Res.* **19-5**:1081