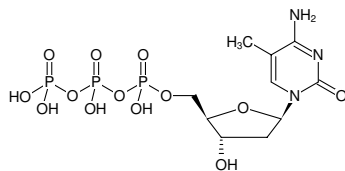


**5-Methyl-dCTP**

mdCTP, 5-mdCTP

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

Cat. No.	Amount
NU-1125S	10 µl (100 mM)
NU-1125L	5 x 10 µl (100 mM)



Structural formula of 5-Methyl-dCTP

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:** C₁₀H₁₈N₃O₁₃P₃ (free acid)**Molecular Weight:** 481.18 g/mol (free acid)**Exact Mass:** 481.01 g/mol (free acid)**CAS#:** 22003-12-9**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 100 mM - 110 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)**Applications:**Incorporation into DNA by
- PCR with *Taq* polymerase ^{in-house data, [1-2]}**Description:**5-methylated DNA probes can be used as methylation reference fragment ^[1-2] or for pull-down of 5-hmC binding proteins from cellular lysate^[3].**Selected References:**

- [1] Szwagierczak *et al.* (2011) Characterization of PvuRts1I endonuclease as a tool to investigate genomic 5-hydroxymethylcytosine. *Nucl. Acids Res.* **39**(12):5149.
- [2] Szwagierczak *et al.* (2010) Sensitive enzymatic quantification of 5-hydroxymethylcytosine in genomic DNA. *Nucl. Acids Res.* **38**(19):e181.
- [3] Lafaye *et al.* (2014) DNA binding of the p21 repressor ZBTB2 is inhibited by cytosine hydroxymethylation. *Biochem. Biophys. Res. Commun.* **446**:341.
- 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.
- Wu *et al.* (2007) Crystal structure of RS21-C6, involved in nucleoside triphosphate pyrophosphohydrolysis. *J. Mol. Biol.* **367**:1405.
- Dezhurov *et al.* (2005) A new highly efficient photoreactive analogue of dCTP. Synthesis, characterization, and application in photoaffinity modification of DNA binding proteins. *Bioconjug. Chem.* **16** (1):215.
- Song *et al.* (2015) Human dCTP pyrophosphatase 1 promotes breast cancer cell growth and stemness through the modulation on 5-methyl-dCTP metabolism and global hypomethylation. *Oncogenesis* **4**: e159.
- Frauer *et al.* (2009) A versatile non-radioactive assay for DNA methyltransferase activity and DNA binding. *Nucleic Acids Research* **37** (3):e22.
- Zeschnigk *et al.* (2009) Massive parallel bisulfite sequencing of CG-rich DNA fragments reveals that methylation of many X-chromosomal CpG islands in female blood DNA is incomplete. *Human Molecular Genetics* **18** (8):1439.
- Holliday *et al.* (2002) DNA methylation and epigenetic inheritance. *Methods.* **27** (2):179.
- Asamizu *et al.* (1999) A large scale structural analysis of cDNAs in a unicellular green alga, *Chlamydomonas reinhardtii*. I. Generation of 3433 non-redundant expressed sequence tags. *DNA Res.* **6** (6):369.
- Ito *et al.* (1998) Systematic method to obtain novel genes that are regulated by mi transcription factor: impaired expression of granzyme B and tryptophan hydroxylase in mi/mi cultured mast cells. *Blood* **91** (9):3210.
- Lefaucheur *et al.* (1998) Evidence for three adult fast myosin heavy chain isoforms in type II skeletal muscle fibers in pigs. *J. Anim. Sci.* **76** (6):1584.
- Wong *et al.* (1997) A novel method for producing partial restriction digestion of

**5-Methyl-dCTP**

mdCTP, 5-mdCTP

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

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.

[3] Lafaye *et al.* (2014) DNA binding of the p21 repressor ZBTB2 is inhibited by cytosine hydroxymethylation. *Biochem. Biophys. Res. Commun.* **446**:341.

Nelson *et al.* (1993) Restriction endonuclease cleavage of 5-methyl-deoxycytosine hemimethylated DNA at high enzyme-to-substrate ratios. *Nucl. Acid. Res.* **21 (3)**:681.