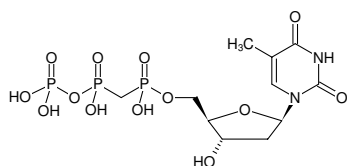


**dTpCpp**

(dTMP CPP)

2'-Deoxythymidine-5'-[( $\alpha,\beta$ )-methylene]triphosphate, Sodium salt

Cat. No.	Amount
NU-895S	50 $\mu$ l (10 mM)
NU-895L	5 x 50 $\mu$ l (10 mM)



Structural formula of dTpCpp

**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:** C<sub>11</sub>H<sub>19</sub>N<sub>2</sub>O<sub>13</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 480.20 g/mol (free acid)**Exact Mass:** 480.01 g/mol (free acid)**Purity:**  $\geq$  95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5  $\pm$  0.5**Spectroscopic Properties:**  $\lambda_{\max}$  267 nm,  $\epsilon$  9.6 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)**Selected References:**

Gowda *et al.* (2012) Low fidelity bypass of O(2)-(3-pyridyl)-4-oxobutylthymine, the most persistent bulky adduct produced by the tobacco specific nitrosamine 4-(methylnitrosamino)-1-(3-pyridyl)-1-butanone by model DNA polymerases. *Chem. Res. Toxicol.* **25** (6):1195.

Beese *et al.* (1993) Crystal structures of the Klenow fragment of DNA polymerase I complexed with deoxynucleoside triphosphate and pyrophosphate. *Biochemistry* **32** (51):14095.