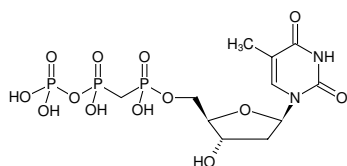


**dTpCcp**

(dTMPCCPP)

2'-Deoxythymidine-5'-[(α,β)-methylene]triphosphate, Sodium salt

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



Structural formula of dTpCcp

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:** C₁₁H₁₉N₂O₁₃P₃ (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:** λ_{\max} 267 nm, ϵ 9.6 L mmol⁻¹ cm⁻¹ (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.