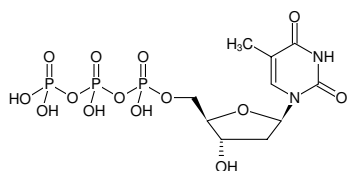


**dTTP - Solid**

2'-Deoxythymidine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1046-20	20 µmol
NU-1046-100	5 x 20 µmol



Structural formula of dTTP - Solid

For research use only!**Shipping:** shipped on blue ice**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:** C₁₀H₁₇N₂O₁₄P₃ (free acid)**Molecular Weight:** 482.17 g/mol (free acid)**CAS#:** 18423-43-3**Purity:** ≥ 95 % (HPLC)**Form:** lyophilised**Spectroscopic Properties:** λ_{max} 267 nm, ε 9.6 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Selected References:**Erlich *et al.* (1988) Primer-directed enzymatic amplification of DNA with a thermostable DNA polymerase. *Science* **29** (239):487.Holland *et al.* (1991) Detection of specific polymerase chain reaction product by utilizing the 5'→3' exonuclease activity of *Thermus aquaticus* DNA polymerase. *Proc. Natl. Acad. Sci. USA* **88** (16):7276.Sanger *et al.* (1977) DNA sequencing with chain-terminating inhibitors. *Proc. Natl. Acad. Sci. USA* **74**:5463.