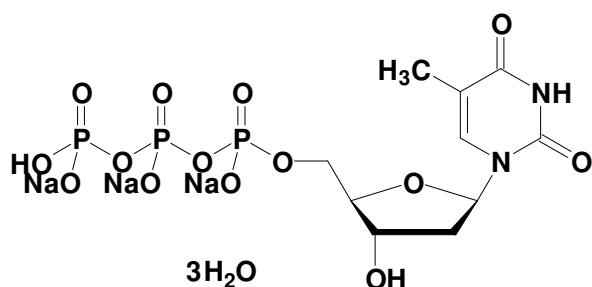


dTTP, lyophilized

2'-Deoxythymidine 5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1004-10	10 mg, lyophilized



Molecular Formula: $C_{10}H_{14}N_2Na_3O_{14}P_3 \times 3H_2O$

Molecular Weight: 602.16

Spectroscopic Properties: $\lambda_{max} = 267 \text{ nm}$;
 $\epsilon = 9.6 \text{ mmol}^{-1} \text{ cm}^{-1}$ (pH 7.5)

Purity: $\geq 95\%$ (HPLC)

Storage conditions:

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

For research use only!

Selected References:

Erlich *et al.* (1988) Primer-directed enzymatic amplification of DNA with a thermostable DNA polymerase. *Science* **29 (239)**:487.

Gelfand *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.