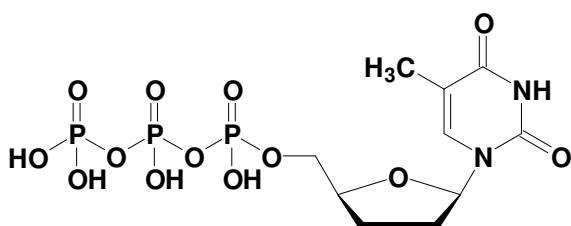


ddTTP 10 mM

Cat. No.	Amount
NU-1018S	200 μ l (2 μ mol)
NU-1018L	1 ml (10 μ mol)



For *in vitro* use only

Quality guaranteed for 12 months

Store at -20°C, short term (up to one week) exposure to ambient temperature possible

Quality control

The dideoxynucleotide solution has a purity of >99% and is functionally tested by chain termination sequencing reactions.

Description

Clear aqueous solution of ultrapure ddTTP at a concentration of 10-11 mM and pH 8.2 +/- 0.2.

ddTTP

2',3'-Dideoxy-thymidine-5'-triphosphate, sodium salt

Molecular formula: C₁₀H₁₇N₂O₁₃P₃ (free acid)

Molecular weight: 466.17 (free acid)

Selected References:

Koji *et al.* (2008) In situ detection of methylated DNA by histone endonuclease-linked detection of methylated DNA sites: a new principle of analysis of DNA methylation. *Histochem Cell Biol* **130**:917.

Sanger *et al.* (1977) DNA sequencing with chain-terminating inhibitors. *Proc. Natl. Acad. Sci. USA* **74**:5463.