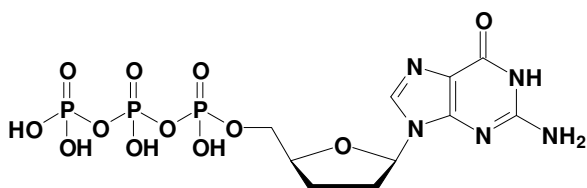


ddGTP 10 mM

Cat. No.	Amount
NU-1017S	200 μ l (2 μ mol)
NU-1017L	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 ddGTP at a concentration of 10-11 mM and pH 8.2 +/- 0.2.

ddGTP

2',3'-Dideoxy-guanosine-5'-triphosphate, sodium salt
Molecular formula: C₁₀H₁₆N₅O₁₂P₃ (free acid)
Molecular weight: 491.18 (free acid)

Selected References:

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