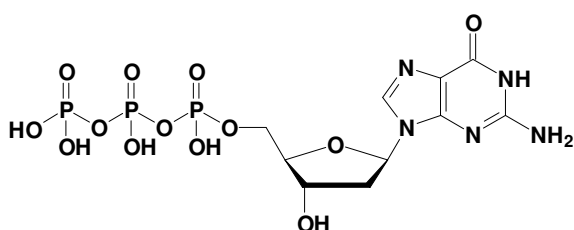


dGTP, solution 100 mM

Cat. No.	Amount
NU-1003	1 ml (100 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

Concentration

100-110 mM

Form

clear aqueous solution, pH 8.5 +/-0.1 (4°C)

Purity

≥99%

Molecular Formula

C₁₀H₁₆N₅O₁₃P₃ (free acid)

Molecular Weight

507.18 (free acid)

Absorbance

absorbance max: 252 nm (pH 7)
ε at absorbance max: 13.7 mmol⁻¹ cm⁻¹

Quality Control Specifications

18 kb long range PCR (template dilution series):
suitable
contamination with bacterial and human DNA:
not detectable
activity of DNase, Protease or Phosphatase:
not detectable

Description

Ultrapure dGTP supplied as clear aqueous solution (pH 8.5).

Selected References:

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.

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

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