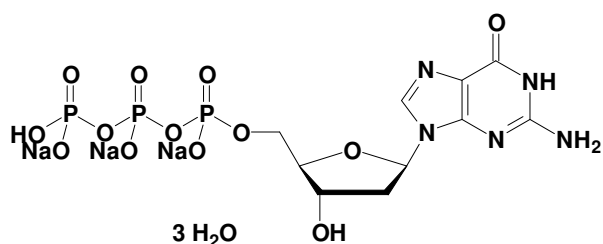


dGTP, lyophilized

2'-Deoxyguanosine 5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1003-10	10 mg, lyophilized



Molecular Formula: C₁₀H₁₃N₅Na₃O₁₃P₃ × 3H₂O

Molecular Weight: 627.16

Spectroscopic Properties: λ_{max} = 252 nm;
 ε = 14.2 mmol⁻¹ cm⁻¹ (pH 7.5)

Purity: at least 95%

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.