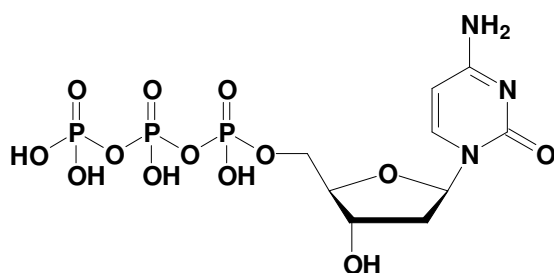


dCTP, solution 100 mM

Cat. No.	Amount
NU-1002	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

C₉H₁₆N₃O₁₃P₃ (free acid)

Molecular Weight

467.15 (free acid)

Spectroscopic Properties: λ_{\max} 271 nm;

ϵ 8900

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 dCTP supplied as clear aqueous solution (pH 8.5).

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.