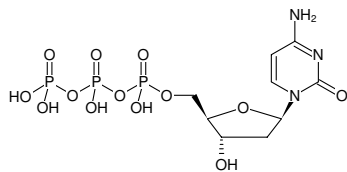




dCTP - Solution

100 mM Sodium salt solution

Cat. No.	Amount
NU-1002L	1 ml (100 mM)
NU-1002-10ML	10 ml (100 mM)
NU-1002-100ML	100 ml (100 mM)



Structural formula of dCTP - Solution

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Shelf Life: 12 months

Molecular Formula: C₉H₁₆N₃O₁₃P₃ (free acid)

Molecular Weight: 467.15 g/mol (free acid)

CAS#: 102783-51-7

Purity: ≥ 99 % (HPLC)

Form: clear aqueous solution

Concentration: 100 mM - 110 mM

pH: 8.5 ± 0.2 (22 °C)

Spectroscopic Properties: λ_{max} 271 nm, ε 8.9 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Description:

dCTP, PCR-grade is supplied as ultrapure aqueous solution (pH 8.5) and suitable for all molecular biology applications including PCR/qPCR, reverse transcription, DNA labeling and DNA sequencing.

Selected References:

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

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.

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