

**dATP - Solution**

100 mM Sodium salt solution

Cat. No.	Amount
NU-1001L	1 ml (100 mM)
NU-1001-10ML	10 ml (100 mM)
NU-1001-100ML	100 ml (100 mM)
NU-1001-200ML	200 ml (100 mM)

For general laboratory use.**Shipping:** shipped on gel packs**Storage Conditions:** store at -20 °C**Shelf Life:** 12 months**Molecular Formula:** C₁₀H₁₆N₅O₁₂P₃ (free acid)**Molecular Weight:** 491.18 g/mol (free acid)**CAS#:** 1927-31-7**Purity:** ≥ 99 % (HPLC)**Form:** clear aqueous solution**Concentration:** 100 mM - 110 mM**pH:** 8.5 ± 0.2 (22 °C)**Spectroscopic Properties:** λ_{max} 259 nm, ε 15.1 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.0)**Description:**

dATP, 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.