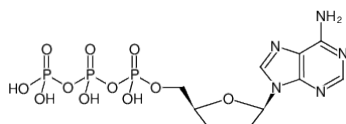




ddATP - high concentration

2',3'-Dideoxyadenosine-5'-triphosphate, Lithium salt

Cat. No.	Amount
NU-291S	10 µl (100 mM)
NU-291L	5 x 10 µl (100 mM)



Structural formula of ddATP - high concentration

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: C₁₀H₁₆N₅O₁₁P₃ (free acid)

Molecular Weight: 475.18 g/mol (free acid)

Exact Mass: 475.01 g/mol (free acid)

CAS#: 24027-80-3

Purity: ≥ 98 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 100 mM - 110 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{max} 259 nm, ε 15.1 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Selected References:

Koji *et al.* (2008) In situ detection of methylated DNA by histo endonuclease-linked detection of methylated DNA sites: a new principle of analysis of DNA methylation. *Histochem Cell Biol* **130**:917.

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