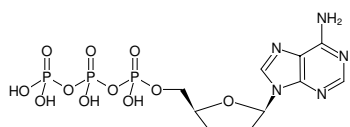


**ddATP**

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

Cat. No.	Amount
NU-1015S	100 µl (1 µmol)
NU-1015L	5 x 100 µl (5 µmol)



Structural formula of ddATP

For in vitro use only!**Shipping:** shipped on blue ice**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:** clear aqueous solution**Concentration:** 10 mM - 11 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.