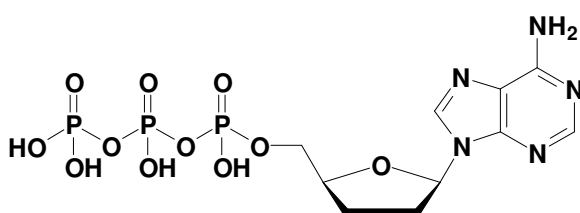


ddATP 10 mM

Cat. No.	Amount
NU-1015S	200 μ l (2 μ mol)
NU-1015L	1 ml (10 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

Quality control

The dideoxynucleotide solution has a purity of >99% and is functionally tested by chain termination sequencing reactions.

Description

Clear aqueous solution of ultrapure ddATP at a concentration of 10-11 mM and pH 8.0 +/-0.2.

ddATP

2',3'-Dideoxy-adenosine-5'-triphosphate, sodium salt
Molecular formula: C₁₀H₁₆N₅O₁₁P₃ (free acid)
Molecular weight: 475.18 (free acid)

Selected References:

- Koji *et al.* (2008) In situ detection of methylated DNA by histone 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.