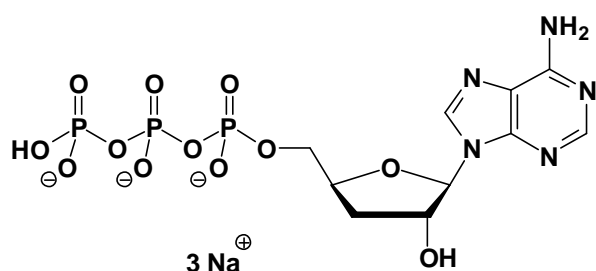


3'-dATP

3'-Deoxy-adenosine-5'-triphosphate, Sodium salt
(Cordycepin triphosphate)

Cat. No.	Amount
NU-1123S	50 Units
NU-1123L	250 Units



Cat. No.: NU-1123

Molecular Formula: C₁₀H₁₃N₅O₁₂P₃ (Anion)

Molecular Weight: 488.16 (Anion)

Purity: > 95%, clear aqueous solution, pH 7.5

Spectroscopic properties: λ_{max} 259 nm; ε 15400

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C.

For research use only!

1 unit = 1 μl of a 10 mM solution

Selected References:

Shuman (1995) Vaccinia virus DNA ligase: specificity, fidelity, and inhibition. *Biochemistry*. **34(49)**:16138.

Labryere *et al.* (1990) Characterization of ATP and calmodulin-binding properties of a truncated form of Bacillus anthracis adenylate cyclase. *Biochemistry*. **29(20)**:4922.

Freist *et al.* (1983) Isoleucyl-tRNA synthetase from Baker's yeast. Catalytic mechanism, 2',3'-specificity and fidelity in aminoacylation of tRNA^{Leu} with isoleucine and valine investigated with initial-rate kinetics using analogs of tRNA, ATP and amino acids. *Eur J Biochem*. **131(1)**:65.