

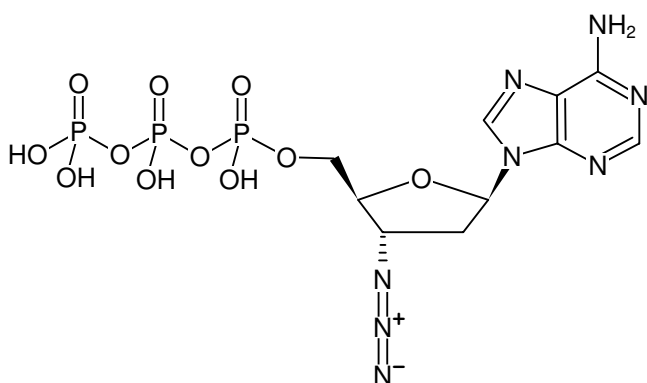
**3'-Azido-2',3'-ddATP**

3'-Azido-2',3'-dideoxyadenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-882S	10 µl (100 mM)
NU-882L	5 x 10 µl (100 mM)

Applications:Enzymatic 3'-click-labeling of RNA^[1]**Selected References:**

[1] Winz *et al.* (2012) Site-specific terminal and internal labeling of RNA by poly (A) polymerase tailing and copper-catalyzed or copper-free strain-promoted click chemistry. *Nucleic Acids Res.* **40** (10):e78.



Structural formula of 3'-Azido-2',3'-ddATP

For research 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:** C₁₀H₁₅N₈O₁₁P₃ (free acid)**Molecular Weight:** 516.19 g/mol (free acid)**Exact Mass:** 516.01 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 100 mM - 110 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ_{max} 258 nm, ε 15.4 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)