

**2'-Azido-2'-dATP**

2'-Azido-dATP

2'-Azido-2'-deoxyadenosine-5'-triphosphate, Sodium salt

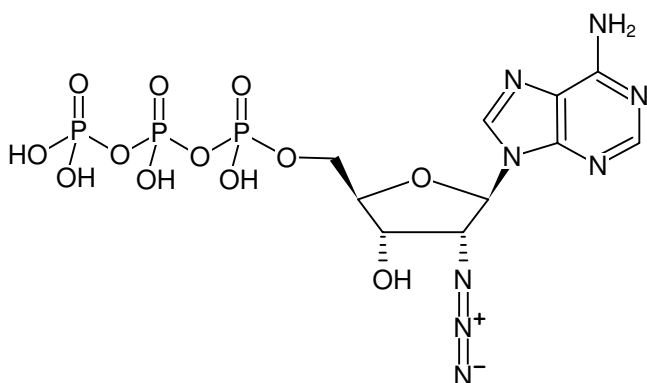
Cat. No.	Amount
NU-976S	10 µl (100 mM)
NU-976L	5 x 10 µl (100 mM)

Selected References:

Randrianjatovo-Gbalou *et al.* (2018) Enzymatic synthesis of random sequences of RNA and RNA analogues by DNA polymerase theta mutants for the generation of aptamer libraries. *Nucleic Acids Res.* doi:10.1093/nar/gky413.

Wang *et al.* (2016) DNA-Catalyzed Introduction of Azide at Tyrosine for Peptide Modification. *Angew. Chem. Int. Ed. Engl.* **55** (34):10052.

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 2'-Azido-2'-dATP

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:** C₁₀H₁₅N₈O₁₂P₃ (free acid)**Molecular Weight:** 532.19 g/mol (free acid)**Exact Mass:** 532.00 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution**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)