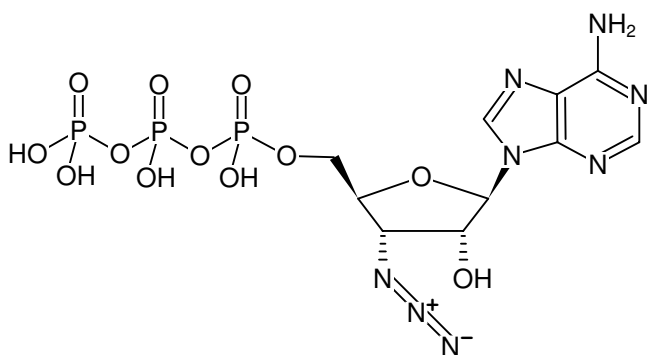


**3'-Azido-3'-dATP**

3'-Azido-3'-deoxyadenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-990	100 µl (10 mM)



Structural formula of 3'-Azido-3'-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:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 259 nm, ε 15.4 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)