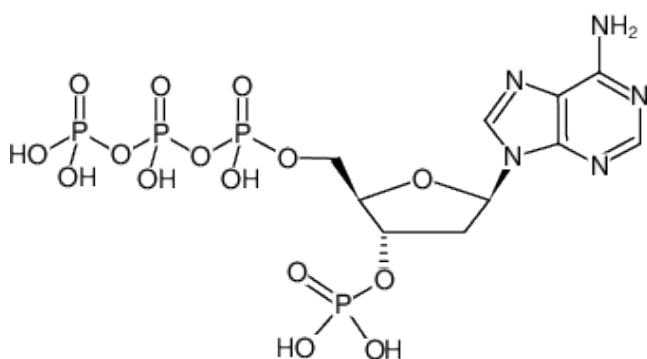


**3'-Phosphate-dATP**

3'-Phosphate-2'-deoxyadenosine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-1221S	200 µl (10 mM)
NU-1221L	5 x 200 µl (10 mM)



Structural formula of 3'-Phosphate-dATP

For general laboratory use.**Shipping:** shipped on gel packs**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:** C₁₀H₁₇N₅O₁₅P₄ (free acid)**Molecular Weight:** 571.16 g/mol (free acid)**Exact Mass:** 570.97 g/mol (free acid)**Purity:** ≥ 98 % (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)