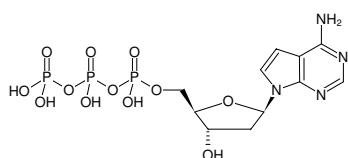


**7-Deaza-dATP**

7-Deaza-2'-deoxyadenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1175S	100 µl (10 mM)
NU-1175L	5 x 100 µl (10 mM)



Structural formula of 7-Deaza-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:** C₁₁H₁₇N₄O₁₂P₃ (free acid)**Molecular Weight:** 490.19 g/mol (free acid)**Exact Mass:** 490.01 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution**Concentration:** 10 mM - 11 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ_{max} 270 nm, ε 11.3 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Selected References:**Pivonková *et al.* (2010) Direct voltammetric analysis of DNA modified with enzymatically incorporated 7-deazapurines. *Anal Chem.* **82** (16):6807.Fletcher *et al.* (1996) Human Telomerase Inhibition by 7'-Deaza-2'-deoxypurine Nucleoside Triphosphates. *Biochemistry* **35** (49):15611.Yamakawa *et al.* (1997) A DNA cycle sequencing reaction that minimizes compression on automated fluorescent sequencers. *Nucleic Acids Research* **25** (6):1311.