

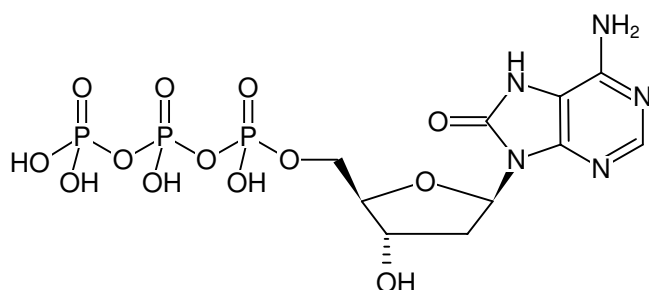
**8-Oxo-dATP**

8-Hydroxy-dATP

8-Oxo-2'-deoxyadenosine-5'-triphosphate, Sodium salt

8-Hydroxy-2'-deoxyadenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-889S	10 µl (100 mM)
NU-889L	5 x 10 µl (100 mM)



Structural formula of 8-Oxo-dATP

For research use only!**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:** C₁₀H₁₆N₅O₁₃P₃ (free acid)**Molecular Weight:** 507.18 g/mol (free acid)**Exact Mass:** 507.00 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} 267 nm, ε 13.6 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Selected References:**

Malvezzi *et al.* (2015) Quantification of pyrophosphate as a universal approach to determine polymerase activity and assay polymerase inhibitors. *Anal. Biochem.* **478**:1.

Kamiya (2010) Mutagenicity of oxidized DNA precursors in living cells: Roles of nucleotide pool sanitization and DNA repair enzymes, and translesion synthesis DNA polymerases. *Mutat. Res.* **703 (1)**:32.

McLennan (2006) The Nudix hydrolase superfamily. *Cell. Mol. Life Sci.* **63 (2)**:123.

Hori *et al.* (2005) In vivo mutagenicities of damaged nucleotides produced by nitric oxide and ionizing radiation. *Biol. Pharm. Bull.* **28 (3)**:520.

Fujikawa *et al.* (1999) The oxidized forms of dATP are substrates for the human MutT homologue, the hMTH1 protein. *J. Biol. Chem.* **274 (26)**:18201.