

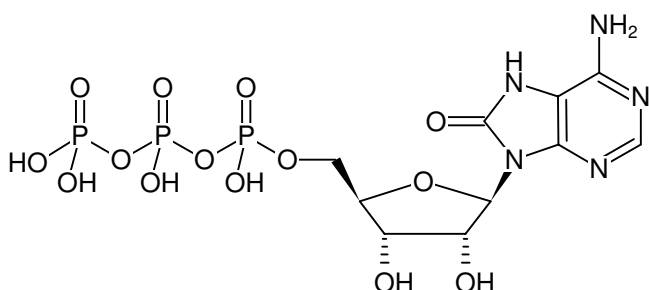
**8-Oxo-ATP**

8-Hydroxy-ATP

8-Oxoadenosine-5'-triphosphate, Sodium salt

8-Hydroxyadenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-953S	10 µl (100 mM)
NU-953L	5 x 10 µl (100 mM)



Structural formula of 8-Oxo-ATP

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:** C₁₀H₁₆N₅O₁₄P₃ (free acid)**Molecular Weight:** 523.18 g/mol (free acid)**Exact Mass:** 522.99 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} 268 nm, ε 12.2 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)