

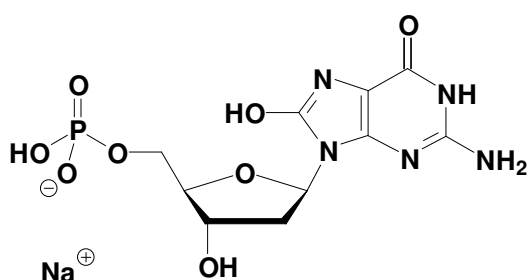
8-Oxo-dGMP

8-Oxo-2'-Deoxy-Guanosine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-1153S	50 Units
NU-1153L	250 Units

Selected References:

Kaminsky *et al.* (2007) The c-Myc Target Gene Rcl (C6orf108) Encodes a Novel Enzyme, Deoxynucleoside 5'-monophosphate N-Glycosidase. *J. Biol. Chem.* **282** (11): 8150.



Cat. No.: NU-1153

Molecular Formula: C₁₀H₁₃N₅O₈P (Anion)

Molecular Weight: 362.21 (Anion)

Purity: > 95%, clear aqueous solution, pH 7.5

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

* 1 unit = 1 µl of a 10 mM solution

For research use only!