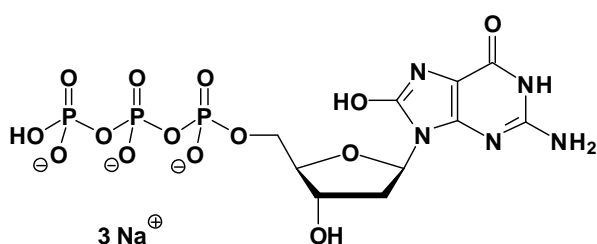


8-Oxo-dGTP

8-Oxo-2'-deoxy-guanosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1117S	30 Units
NU-1117L	150 Units



Cat. No.: NU-1117

Molecular Formula: C₁₀H₁₂N₅O₁₄P₃ (Anion)

Molecular Weight: 519.15 (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.

For research use only!

1 unit = 1 µl of a 10 mM solution

Selected References:

Kai *et al.* (2002) An oxidized nucleotide affects DNA replication through activation of protein kinases in *Xenopus* egg lysates. *Nucl. Acids Res.* **30** (2):569.

Canitrot *et al.* (2000) Nucleotide excision repair DNA synthesis by excess DNA polymerase beta: a potential source of genetic instability in cancer cells. *FASEB J.* **14** (12):1765.

Nampalli *et al.* (2000) Efficient synthesis of 8-oxo-dGTP: A mutagenic nucleotide. *Bioorg. Med. Chem. Lett.* **10** (15):1677.

Zaccolo, M. *et al.* (1996) *J. Mol. Biol.* **255**:589

Lethal mutagenesis:

Loeb, L. *et al.* (1999) *Proc. Natl. Acad. Sci. USA* **96**:1492