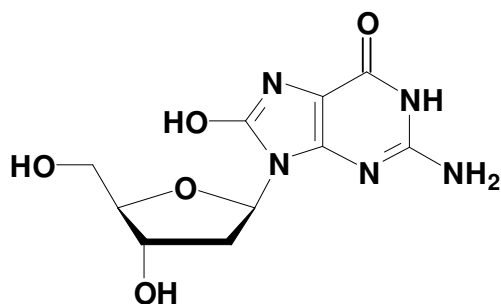


8-Oxo-2'-deoxyguanosine

Cat. No.	Amount
N-1045	25 mg

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



Cat. No.: N-1045

Molecular Formula: C₁₀H₁₃N₅O₅

Molecular Weight: 283.24

Purity: > 98%

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at +4°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

For research use only!

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.