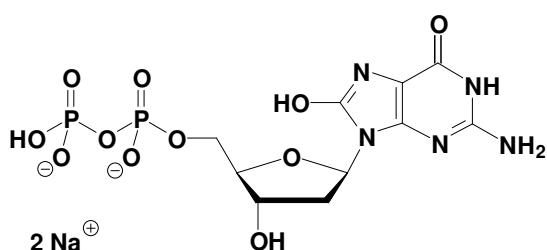


## 8-Oxo-dGDP

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

Cat. No.	Amount
NU-1158S	20 Units
NU-1158L	100 Units



**Cat. No.:** NU-1158

**Molecular Formula:** C<sub>10</sub>H<sub>13</sub>N<sub>5</sub>O<sub>11</sub>P<sub>2</sub> (Anion)

**Molecular Weight:** 441.18 (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!**

### 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.