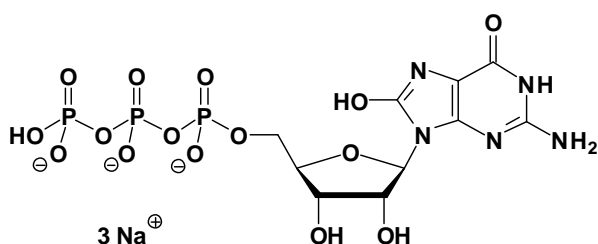


## 8-Oxo-GTP

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

Cat. No.	Amount
NU-1116S	50 Units
NU-1116L	250 Units



**Cat. No.:** NU-1116

**Molecular Formula:** C<sub>10</sub>H<sub>12</sub>N<sub>5</sub>O<sub>15</sub>P<sub>3</sub> (Anion)

**Molecular Weight:** 535.15 (Anion)

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

**Spectroscopic properties:** λ<sub>max</sub> 245nm; ε 12300

### 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:

Tanaka *et al.* (2005) Novel reaction mechanism of GTP cyclohydrolase I. High-resolution X-ray crystallography of *Thermus thermophilus* HB8 enzyme complexed with a transition state analogue, the 8-oxoguanine derivative. *J. Biochem. (Tokyo)* **138** (3):263.

Nakabeppu (2001) Molecular genetics and structural biology of human MutT homolog, MTH1. *Mutat Res.* **477** (1-2):59.

Kobayashi *et al.* (1998) Potential of *Escherichia coli* GTP cyclohydrolase II for hydrolyzing 8-oxo-dGTP, a mutagenic substrate for DNA synthesis. *J. Biol. Chem.* **273** (41):26394.

Kuryavyi *et al.* (1989) Kinetics of inhibition of *Escherichia coli* RNA polymerase-catalyzed synthesis of the dinucleotide pppApU by 8-oxo-GTP and 8-BrGTP on the A1 promoter of Bacteriophage- $\tau$  delta-d111 DNA using a limited set of substrates. *Mol. Biol.* **23** (3):648.