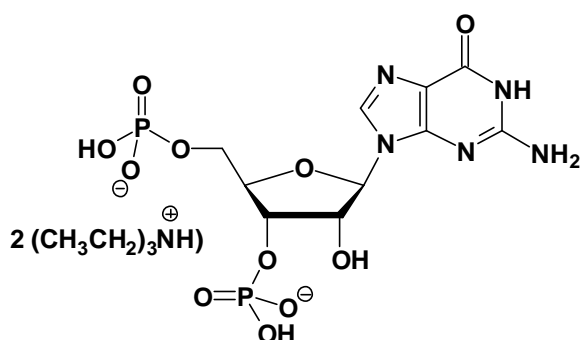


## Guanosine-3',5'-bisphosphate (pGp)

Guanosine-3',5'-bisphosphate, Triethylammonium salt

Cat. No.	Amount
NU-1107S	100 Units
NU-1107L	500 Units



**Cat. No.:** NU-1107

**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.

### For research use only!

1 unit = 1 µl of a 10 mM solution

### Selected References:

Kvint *et al.* (2000) Emergency derepression: stringency allows RNA polymerase to override negative control by an active repressor. *Mol. Microbiol.* **35** (2):435.

Acharya *et al.* (1999) The transmission of the electronic character of guanin-9-yl drives the sugar-phosphate backbone torsions in guanosine 3',5'-bisphosphate. *Angew. Chem. Int. Edit.* **38** (24):3645.

Ishikawa *et al.* (1996) Crystal structure of ribonuclease T1 carboxymethylated at Glu58 in complex with 2'-GMP. *Biochemistry-US* **35** (25):8329.

Atgie *et al.* (1993) Specific decrease of mitochondrial thermogenic capacity in brown adipose-tissue of obese SHR/N-CP rats. *Am. J. Physiol.* **265** (6):c1674 Part 1.

Lenz *et al.* (1993) 3-dimensional structure of the ternary complex between Ribonuclease-T1, guanosine-3',5'-bisphosphate and inorganic-phosphate at 0.19 nm resolution. *Eur. J. Biochem.* **211** (1-2):311.

Lenz *et al.* (1991) X-ray-analysis of cubic-crystals of the complex formed between Ribonuclease-T1 and guanosine-3',5'- bisphosphate. *Acta Crystallogr. B* **47**:521.

Plesner (1984) Guanosine 3',5'-diphosphate is found in the medium of eukaryotic cells during G1 to S-phase transition. *H-S Z. Physiol. Chem.* **365** (6):608.