

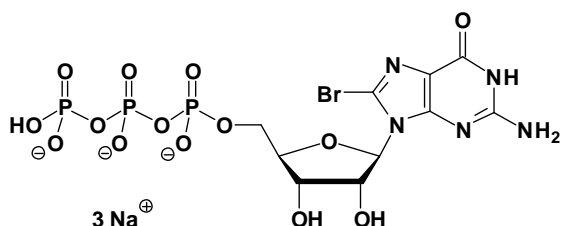
## 8Br-GTP

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

Cat. No.	Amount
NU-118S	50 Units
NU-118L	250 Units

### Selected References:

Läppchen *et al.* (2005) GTP Analogue Inhibits Polymerization and GTPase Activity of the Bacterial Protein FtsZ without Affecting Its Eukaryotic Homologue Tubulin. *Biochemistry* **44** (21):7879.



**Cat. No.:** NU-118

**Molecular Formula:**  $C_{10}H_{11}N_5O_{14}P_3Br$  (Anion)

**Molecular Weight:** 598.04 (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  $\mu$ l of a 10 mM solution