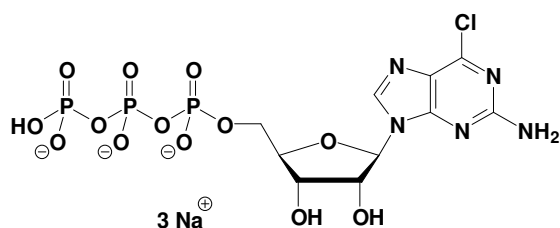


## 6Cl-GTP

6-Chloro-Guanosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-143S	50 Units
NU-143L	250 Units



**Cat. No.:** NU-143

**Molecular Formula:**  $C_{10}H_{12}N_5O_{13}P_3Cl$  (Anion)

**Molecular Weight:** 538.60 (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  $\mu$ l of a 10 mM solution

**For research use only!**