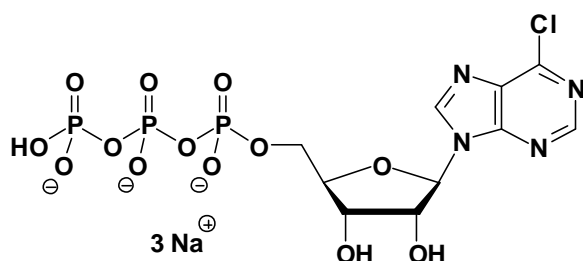


6-Chlorpurine-riboside-5'-triphosphate

6-Chlorpurine-riboside-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1109S	100 Units
NU-1109L	500 Units



Cat. No.: NU-1109

Molecular Formula: $C_{10}H_{11}N_4O_{13}P_3Cl$ (Anion)

Molecular Weight: 523.59 (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 $< -20^{\circ}C$. 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 μ l of a 10 mM solution

Selected References:

Patzelt *et al.* (1974) Study of (Na+K)-ATPase with 6-chlorpurine triphosphate, 6-mercaptopurine triphosphate and dinitrophenyl-6-thiopurine triphosphate. *H-S Z. Physiol. Chem.* **355** (10):1237.