



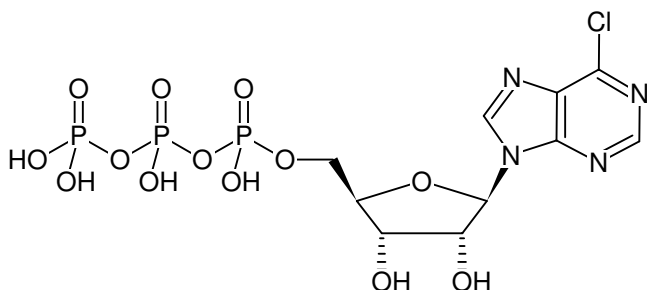
6-Chloropurine-riboside-5'-triphosphate

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

Cat. No.	Amount
NU-1109S	10 µl (100 mM)
NU-1109L	5 x 10 µl (100 mM)

Selected References:

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



Structural formula of 6-Chloropurine-riboside-5'-triphosphate

For research use only!

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: C₁₀H₁₄N₄O₁₃P₃Cl (free acid)

Molecular Weight: 526.61 g/mol (free acid)

Exact Mass: 525.95 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 100 mM - 110 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{max} 263 nm, ε 8.9 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)