

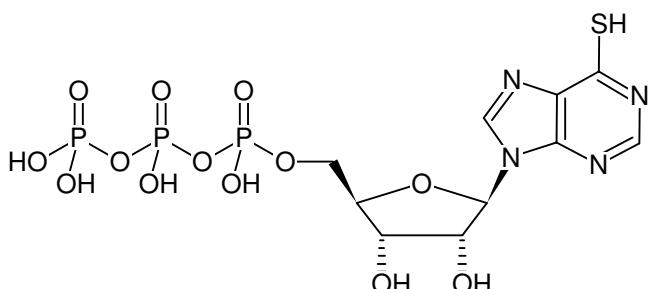


6-Mercaptopurine-riboside-5'-triphosphate

6-Thio-Inosine-5'-triphosphate

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

Cat. No.	Amount
NU-1110S	100 µl (10 mM)
NU-1110L	5 x 100 µl (10 mM)



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

For general laboratory use.

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: C₁₀H₁₅N₄O₁₃P₃S (free acid)

Molecular Weight: 524.22 g/mol (free acid)

Exact Mass: 523.96 g/mol (free acid)

CAS#: 27652-34-2

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 10 mM - 11 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{max} 322 nm, ε 27.3 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Applications:

Competitive inhibitor of T4 RNA ligase^[1]

Kinetic parameters with thiopurine methyltransferase^[2]

Interaction with (sodium-potassium ion)-dependent ATPase^[3]

Selected References:

[1] Labeikyte *et al.* (1995) Structural features of nucleotide binding in T4 RNA ligase determined from inhibition by structural analogs of ATP. *Biologija* **3**:13.

[2] Deininger *et al.* (1994) Purine substrates for human thiopurine methyltransferase. *Biochemical Pharmacology* **48**:2135.

[3] Erdmann *et al.* (1976) The cardiac glycoside receptor: its properties and its correlation to nucleotide binding sites, phosphointermediate, and Na-K-dependent ATPase activity. *Recent Adv. Stud. Cardiac Struct. Metab.* **9**:329.

Bierau *et al.* (2007) Pharmacogenetic significance of inosine triphosphatase. *Pharmacogenomics* **8** (9):1221.

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.