

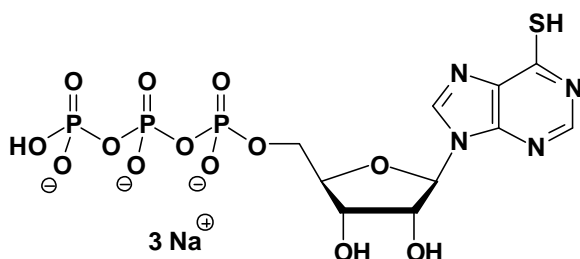
## 6-Mercaptopurine-riboside-5'-triphosphate

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

Cat. No.	Amount
NU-1110S	100 Units
NU-1110L	500 Units

### 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.



**Cat. No.:** NU-1110

**Molecular Formula:**  $C_{10}H_{12}N_4O_{13}P_3S$  (Anion)

**Molecular Weight:** 521.20 (Anion)

**Purity:** > 95%, clear aqueous solution, pH 7.5

**Spectroscopic properties:**  $\lambda_{max}$  322 nm;  $\epsilon$  27300

### 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