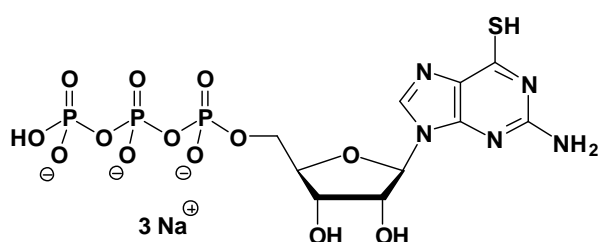


6-Thio-GTP

6-Thio-guanosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1106S	150 Units
NU-1106L	750 Units



Cat. No.: NU-1106

Molecular Formula: $C_{10}H_{13}N_5O_{13}P_3S$ (Anion)

Molecular Weight: 536.22 (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.

For research use only!

1 unit = 1 μ l of a 10 mM solution

Selected References:

Du *et al.* (2007) Enhanced cardiac allograft survival by Vav1-Rac signalling blockade in a mouse model. *Transplant Immunology* **18 (1)**: 53

Pestova *et al.* (2006) Specific functional interactions of nucleotides at key -3 and +4 positions flanking the initiation codon with components of the mammalian 48S translation initiation complex. *Genes&Development* **20**:624.

Tiede *et al.* (2003) CD28-dependent Rac1 activation is the molecular target of azathioprine in primary human CD4 + T lymphocytes. *J. Clin. Invest.* **111**:1133.

Fishback *et al.* (1984) Interaction of 6-mercapto-GTP with bovine brain tubulin - equilibrium aspects. *J. Biol. Chem.* **259 (3)**:1968.

Poland *et al.* (1997) Entrapment of 6-thiophosphoryl-IMP in the active site of crystalline adenylosuccinate synthetase from *Escherichia coli*. *J. Biol. Chem.* **272 (24)**:15200.

Yarbrough *et al.* (1985) Kinetics of interaction of 2-amino-6-mercapto-9-beta-ribofuranosylpurine 5'-triphosphate with bovine brain tubulin. *Biochemistry-US* **24 (7)**:1708.