

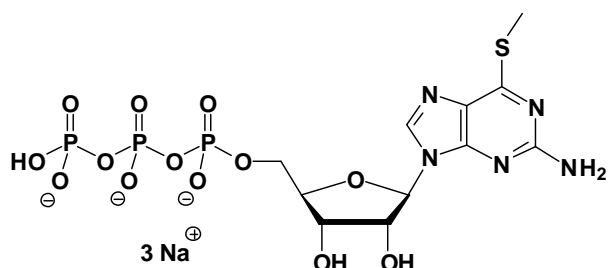
6-Methylthio-GTP

6-Methylthioguanosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1130S	100 Units
NU-1130L	500 Units

Selected References:

Nelson *et al.* (1977) Oxypurine and 6-thiopurine nucleoside triphosphate formation in human erythrocytes. *Adv. Exp. Med. Biol.* **76A**:121.



Cat. No.: NU-1130

Molecular Formula: C₁₁H₁₅N₅O₁₃P₃S (Anion)

Molecular Weight: 550.24 (Anion)

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

Spectroscopic properties: λ_{max} 242 nm;
ε 12800

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 μl of a 10 mM solution