

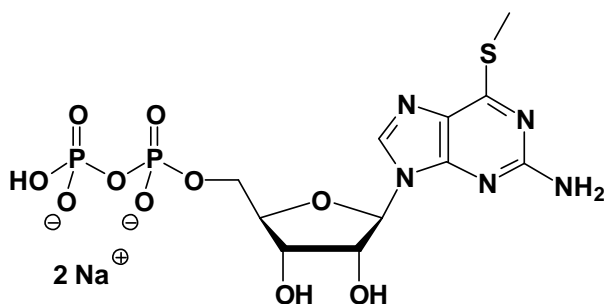
6-Methylthio-GDP

6-Methylthioguanosine-5'-diphosphate, Sodium salt

Cat. No.	Amount
NU-1129S	50 Units
NU-1129L	250 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-1129

Molecular Formula: $C_{11}H_{15}N_5O_{10}P_2S$ (Anion)

Molecular Weight: 471.27 (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 μ l of a 10 mM solution