

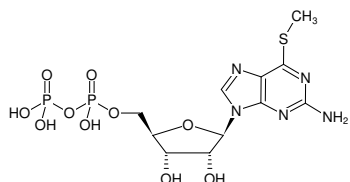
**6-Methylthio-GDP**

6-Methylthioguanosine-5'-diphosphate, Sodium salt

Cat. No.	Amount
NU-1129S	50 µl (10 mM)
NU-1129L	5 x 50 µl (10 mM)

Selected References:

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



Structural formula of 6-Methylthio-GDP

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:** C₁₁H₁₇N₅O₁₀P₂S (free acid)**Molecular Weight:** 473.29 g/mol (free acid)**Exact Mass:** 473.02 g/mol (free acid)**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} 242 nm, ε 12.8 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)