

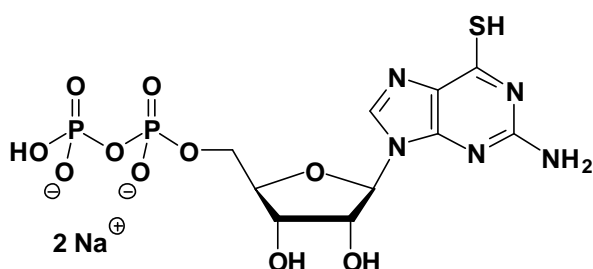
6-Thio-GDP

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

Cat. No.	Amount
NU-1120S	50 Units
NU-1120L	250 Units

Selected References:

Weng *et. al.* (1994) Hydrogen bond interactions of G proteins with the guanine ring moiety of guanine nucleotides. *Protein Sci.* **1**:22.



Cat. No.: NU-1120

Molecular Formula: C₁₀H₁₃N₅O₁₀P₂S (Anion)

Molecular Weight: 457.24 (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 µl of a 10 mM solution