

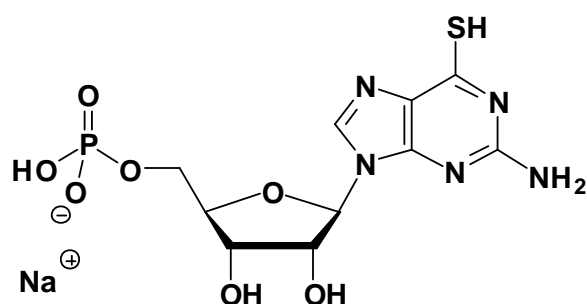
6-Thio-GMP

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

Cat. No.	Amount
NU-1121S	150 Units
NU-1121L	750 Units

Selected References:

Courtois *et. al.* (1969) Interaction entre pigments et acides nucleiques
Étude sur l'interaction de l'Actinomycine D avec les nucléosides par
Dichroïsme Circulaire. *FEBS Lett.* **3**:330.



Cat. No.: NU-1121

Molecular Formula: C₁₀H₁₃N₅O₇PS (Anion)

Molecular Weight: 378.02 (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