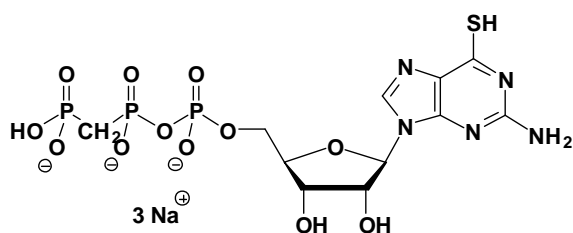


6-Thio-GppCp (6-Thio-GMPPCP)

6-Thio-guanosine-5'-[(β,γ)-methylene]triphosphate, Sodium salt

Cat. No.	Amount
NU-418S	50 Units
NU-418L	250 Units



Cat. No.: NU-418

Molecular Formula: $C_{11}H_{15}N_5O_{12}P_3S$ (Anion)

Molecular Weight: 534.24 (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 $< -20^{\circ}\text{C}$. 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