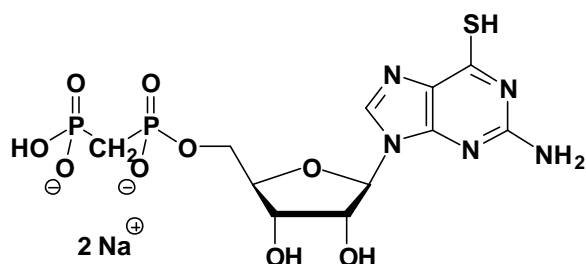


6-Thio-GpCp (6-Thio-GMPCP)

6-Thio-guanosine-5'-[(α,β)-methylene]diphosphate, Sodium salt

Cat. No.	Amount
NU-417S	30 Units
NU-417L	150 Units



Cat. No.: NU-417

Molecular Formula: C₁₁H₁₅N₅O₉P₂S (Anion)

Molecular Weight: 455.27 (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 μ l of a 10 mM solution