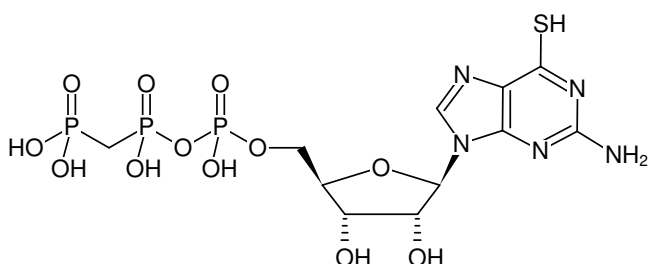


**6-Thio-GppCp**

(6-Thio-GMPPCP)

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

Cat. No.	Amount
NU-418S	50 μ l (10 mM)
NU-418L	5 x 50 μ l (10 mM)



Structural formula of 6-Thio-GppCp

For research use only!**Shipping:** shipped on blue ice**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:** C₁₁H₁₈N₅O₁₂P₃S (free acid)**Molecular Weight:** 537.27 g/mol (free acid)**Exact Mass:** 537.99 g/mol (free acid)**Purity:** \geq 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 \pm 0.5**Spectroscopic Properties:** λ_{\max} 257 nm, ϵ 8.8 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5), λ_{\max} 342 nm, ϵ 24.8 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)