

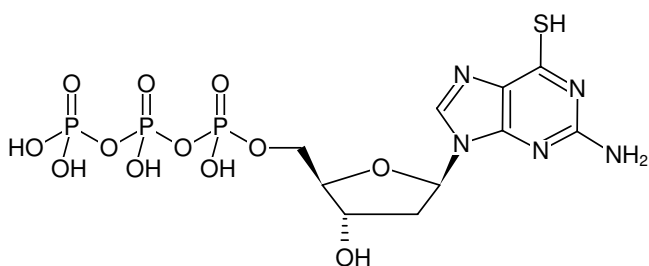
**6-Thio-dGTP**

6-Thio-2'-deoxyguanosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1213S	50 µl (10 mM)
NU-1213L	5 x 50 µl (10 mM)

Selected References:

Tendian *et al.* (2000) Interaction of deoxyguanosine nucleotide analogs with human telomerase. *Mol. Pharmacol.* **57** (4):695.



Structural formula of 6-Thio-dGTP

For general laboratory use.**Shipping:** shipped on gel packs**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:** C₁₀H₁₆N₅O₁₂P₃S (free acid)**Molecular Weight:** 523.24 g/mol (free acid)**Exact Mass:** 522.97 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±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)