

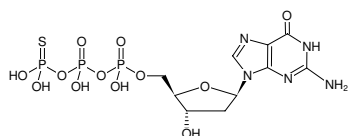
**dGTPyS**

2'-dGua-5'-PPP[S]

2'-Deoxyguanosine-5'-(γ-thio)-triphosphate, Lithium salt

2'-Deoxyguanosine-5'-(3-thio)-triphosphate, 2'-Deoxyguanosine-5'-(3-thiotriphosphate)

Cat. No.	Amount
NU-266	5 mg



Structural formula of dGTPyS

The limited stability of dGTPyS makes it impossible to keep it stocked and to offer it as catalog product that is available immediately. Therefore, dGTPyS is freshly prepared per order. For this reason the price is higher than for stable nucleotides which can be synthesized and kept in stock.

For research use only!**Shipping:** shipped on dry ice**Storage Conditions:** store at -20 °C**Shelf Life:** 6 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:** ≥ 90 % (HPLC), contains < 10 % GDP**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ_{max} 252 nm, ε 13.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)