

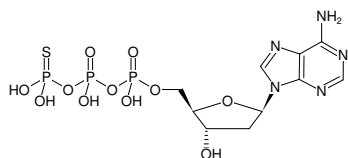
**dATPyS**

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

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

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

Cat. No.	Amount
NU-265	5 mg



Structural formula of dATPyS

For general laboratory use.**Shipping:** shipped on dry ice**Storage Conditions:** store at -20 °C**Shelf Life:** 6 months after date of delivery**Molecular Formula:** C₁₀H₁₆N₅O₁₁P₃S (free acid)**Molecular Weight:** 507.24 g/mol (free acid)**Exact Mass:** 506.98 g/mol (free acid)**Purity:** ≥ 90 % (HPLC), contains < 10 % ADP**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ_{max} 259 nm, ε 15.4 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)