

**dATPyS**

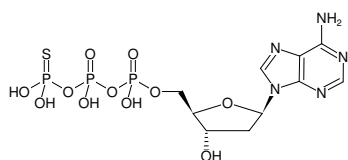
2'-dAdo-5'-PPP[S]

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

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



Structural formula of dATPyS

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: 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:** white to off-white solid**Spectroscopic Properties:** λ_{max} 259 nm, ε 15.4 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)