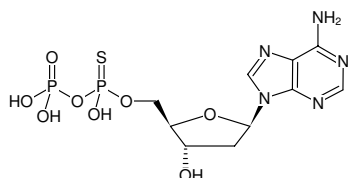


**dADPαS**

2'-Deoxyadenosine-5'-(α-thio)-diphosphate, Triethylammonium salt

Cat. No.	Amount
NU-1165S	20 µl (10 mM)
NU-1165L	5 x 20 µl (10 mM)



Structural formula of dADPαS

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:** C₁₀H₁₅N₅O₈P₂S (free acid)**Molecular Weight:** 427.26 g/mol (free acid)**Exact Mass:** 427.01 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} 259 nm, ε 15.4 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)