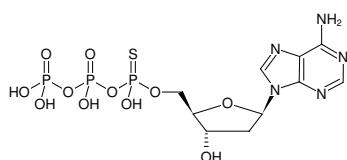


**dATPαS**2'-Deoxyadenosine-5'-(α-thio)-triphosphate, Sodium salt; (Mixture of R_p and S_p isomers)

Cat. No.	Amount
NU-426S	100 µl (10 mM)
NU-426L	5 x 100 µl (10 mM)



Structural formula of dATPαS

For research use only!**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:** C₁₀H₁₆N₅O₁₁P₃S (free acid)**Molecular Weight:** 507.24 g/mol (free acid)**Exact Mass:** 506.98 g/mol (free acid)**CAS#:** 64145-28-4**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)**Selected References:**Gaur *et al.* (1993) Enzymatic RNA synthesis with deoxynucleoside 5'-O-(1-thiotriphosphates). *FEBS Lett.* **315** (1):56.Abbotts *et al.* (1988) Studies on the mechanism of Escherichia coli DNA polymerase I large fragment. Effect of template sequence and substrate variation on termination of synthesis. *J. Biol. Chem.* **263** (29):15094.Nakamaye *et al.* (1988) Direct sequencing of polymerase chain reaction amplified DNA fragments through the incorporation of deoxynucleoside alpha-thiotriphosphates. *Nucleic Acids Res.* **16** (21):9947.Nyren *et al.* (1997) Detection of single-base changes using a bioluminometric primer extension assay. *Anal. Biochem.* **244** (2):367.