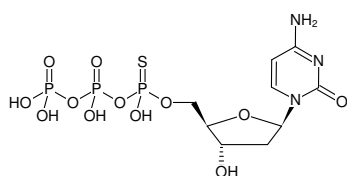


**dCTPaS**

2'-Deoxycytidine-5'-(α -thio)-triphosphate, Sodium salt; (1 : 1 Mixture of R_p and S_p isomers)

Cat. No.	Amount
NU-425S	100 μ l (10 mM)
NU-425L	5 x 100 μ l (10 mM)



Structural formula of dCTPaS

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: 12 months after date of delivery

Molecular Formula: C₉H₁₆N₃O₁₂P₃S (free acid)

Molecular Weight: 483.22 g/mol (free acid)

Exact Mass: 482.97 g/mol (free acid)

CAS#: 64145-29-5

Purity: \geq 95 % (HPLC)

Form: clear aqueous solution

Concentration: 10 mM - 11 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{\max} 271 nm, ϵ 9.0 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.