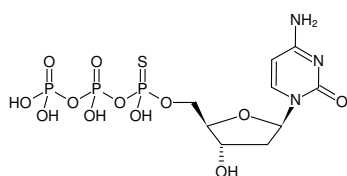


**dCTPaS**2'-Deoxycytidine-5'-( $\alpha$ -thio)-triphosphate, Sodium salt; (Mixture of R<sub>p</sub> and S<sub>p</sub> isomers)

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



Structural formula of dCTPaS

**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:** C<sub>9</sub>H<sub>16</sub>N<sub>3</sub>O<sub>12</sub>P<sub>3</sub>S (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:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5  $\pm$  0.5**Spectroscopic Properties:**  $\lambda_{\max}$  271 nm,  $\epsilon$  9.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (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.