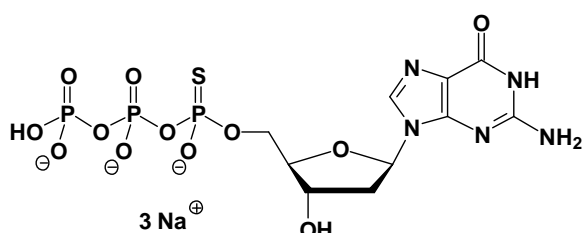


dGTP α S

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

Cat. No.	Amount
NU-424S	100 Units
NU-424L	500 Units



Cat. No.: NU-424

Molecular Formula: C₁₀H₁₃N₅O₁₂P₃S (Anion)

Molecular Weight: 520.22 (Anion)

Purity: > 95%, clear aqueous solution, pH 7.5

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

For research use only!

1 unit = 1 μ l of a 10 mM solution

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.