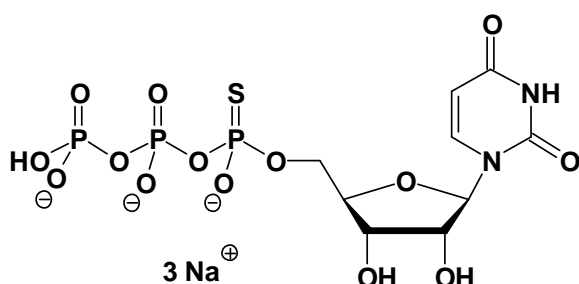


UTP α S

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

Cat. No.	Amount
NU-411S	50 Units
NU-411L	250 Units



Cat. No.: NU-411

Molecular Formula: C₉H₁₂N₂O₁₄P₃S (Anion)

Molecular Weight: 497.18 (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:

Oyelere *et al.* (2002) pKa perturbation in genomic hepatitis delta virus ribozyme catalysis evidenced by nucleotide analogue interference mapping. *Biochemistry-US* **41** (11):3667.

Basu *et al.* (2001) Biochemical detection of monovalent metal ion binding sites within RNA. *Methods* **23** (3):264.

Oyelere *et al.* (2000) Biochemical detection of cytidine protonation within RNA. *J. Am. Chem. Soc.* **122** (42):10259.

Vortler *et al.* (2000) Phosphorothioate modification of RNA for stereochemical and interference analyses. *Method. Enzymol.* **317**:74.

Ryder *et al.* (2000) Chemical probing of RNA by nucleotide analog interference mapping. *Method. Enzymol.* **317**:92.

Burgess *et al.* (2000) Syntheses of nucleoside triphosphates. *Chem. Rev.* **100** (6):2047.

Ryder *et al.* (1999) Nucleotide analog interference mapping. *Methods* **18** (1):38.

Strobel (1999) A chemogenetic approach to RNA function/structure analysis. *Curr. Opin. Struc. Biol.* **9** (3):346.

Boudvillain *et al.* (1998) Defining functional groups, core structural features and inter-domain tertiary contacts essential for group II intron self-splicing: a NAIM analysis. *EMBO J.* **17** (23):7091.

Basu *et al.* (1998) A specific monovalent metal ion integral to the AA platform of the RNA tetraloop receptor. *Nat. Struct. Biol.* **5** (11):986.

Ortoleva-Donnelly *et al.* (1998) Identifying RNA minor groove tertiary contacts by nucleotide analogue interference mapping with N-2-methylguanosine. *Biochemistry-US* **37** (37):12933.

Arabshahi *et al.* (1994) A simplified procedure for synthesizing nucleoside 1-thiotriphosphates - dATP(α)S, dGTP(α)S, UTP- α -S, and dTTP(α)S. *Biochem. Biophys. Res. Co.* **204** (1):150.