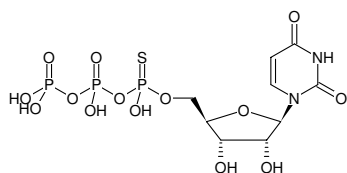


**UTP α S**

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

Cat. No.	Amount
NU-411S	25 μ l (100 mM)
NU-411L	5 x 25 μ l (100 mM)



Structural formula of UTP α S

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: C₉H₁₅N₂O₁₄P₃S (free acid)

Molecular Weight: 500.20 g/mol (free acid)

Exact Mass: 499.95 g/mol (free acid)

CAS#: 71214-29-4

Purity: \geq 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 100 mM - 110 mM

pH: 7.5 \pm 0.5

Spectroscopic Properties: λ_{\max} 262 nm, ϵ 10.0 L mmol⁻¹ cm⁻¹
(Tris-HCl pH 7.5)

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.