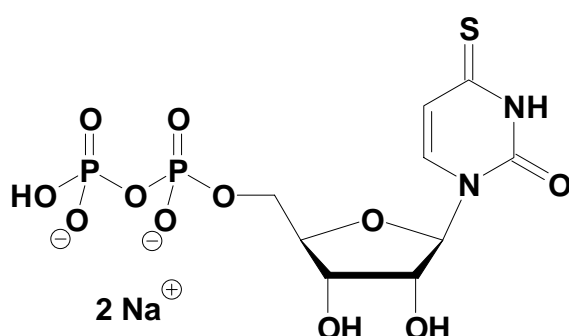


4-Thio-UDP

4-Thio-Uridine-5'-diphosphate, Sodium salt

Cat. No.	Amount
NU-1155S	20 Units
NU-1155L	100 Units



Cat. No.: NU-1155

Molecular Formula: C₉H₁₁N₂O₁₂P₂S (Anion)

Molecular Weight: 433.20 (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.

* 1 unit = 1 µl of a 10 mM solution

For research use only!

Selected References:

Zaher *et al.* (2006) A general RNA-capping ribozyme retains stereochemistry during cap exchange. *J. Am. Chem. Soc.* **128** (42):13894.

Kwon *et al.* (2001) DNA sequencing and genotyping by transcriptional synthesis of chain-terminated RNA ladders and MALDI-TOF mass spectrometry. *Nucleic Acids Res.* **29** (3):e11.

Testa *et al.* (1999) Thermodynamics of RNA-RNA Duplexes with 2- or 4-Thiouridines: Implications for Antisense Design and Targeting a Group I Intron. *Biochemistry* **38**:16655.

Dontsova *et al.* (1994) Stem-loop IV of 5S rRNA lies close to the peptidyltransferase center. *Proc. Natl. Acad. Sci. USA* **91** (10):4125.

Sheng *et al.* (1993) Active site labeling of HIV-1 reverse transcriptase. *Biochemistry* **32** (18):4938.

Khanna *et al.* (1991) Photoaffinity labelling of the pea chloroplast transcriptional complex by nascent RNA in vitro. *Nucleic Acids Res.* **19** (18):4849.

Dissingner *et al.* (1990) Active site labeling of Escherichia coli transcription elongation complexes with 5-[4-azidophenacyl]thio)uridine 5'-triphosphate. *J. Biol. Chem.* **265** (13):7662.

Tanner *et al.* (1988) Binding interactions between yeast tRNA ligase and a precursor transfer ribonucleic acid containing two photoreactive uridine analogues. *Biochemistry* **27** (24):8852.

Bartholomew *et al.* (1986) RNA contacts subunits Ilo and IIc in HeLa RNA polymerase II transcription complexes. *J. Biol. Chem.* **261** (30):14226.

Eshaghpour *et al.* (1979) Specific chemical labeling of DNA fragments. *Nucleic Acids Res.* **7** (6):1485.