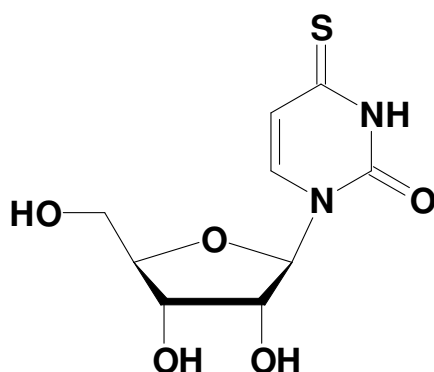


4-Thio-uridine

Cat. No.	Amount
N-1043	100 mg



Cat. No.: N-1043

Molecular Formula: C₉H₁₂N₂O₅S

Molecular Weight: 260.26

Purity: > 98%

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at +4°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

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 labeling of the pea chloroplast transcriptional complex by nascent RNA in vitro. *Nucleic Acids Res.* **19** (18):4849.

Dissing *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.