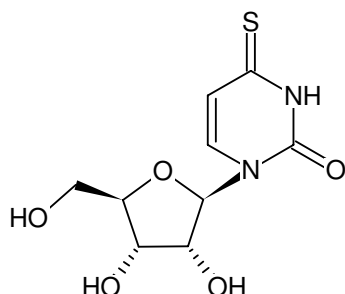




4-Thio-uridine

Cat. No.	Amount
N-1090-100	100 mg
N-1090-250	250 mg



Structural formula of 4-Thio-uridine

For general laboratory use.**Shipping:** shipped at ambient temperature**Storage Conditions:** store at -20 °C**Shelf Life:** 30 months after date of delivery**Molecular Formula:** C₉H₁₂N₂O₅S**Molecular Weight:** 260.27 g/mol**Exact Mass:** 260.04 g/mol**CAS#:** 13957-31-8**Purity:** ≥ 98 % (HPLC)**Form:** solid**Color:** white to off-white**Solubility:** methanol**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.

Dissinger *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.

Favre *et al.* (1986) 4-Thiouridine photosensitized RNA-protein crosslinking in mammalian cells. *Biochem. Biophys. Res. Commun.* **141** (2):847.

Hajnsdorf *et al.* (1987) RNA protein crosslinks introduced into E. coli ribosomes by use of the intrinsic probe 4-thiouridine. *Photochem. Photobiol.* **45** (4):445.