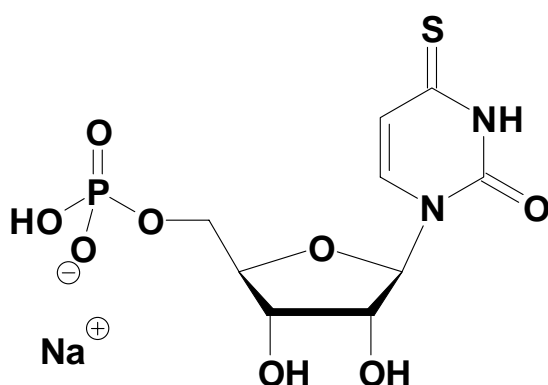


4-Thio-UMP

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

Cat. No.	Amount
NU-1154S	50 Units
NU-1154L	250 Units



Cat. No.: NU-1154

Molecular Formula: C₉H₁₁N₂O₉PS (Anion)

Molecular Weight: 354.22 (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:

Sohn *et al.* (2005) Sequential multiple functions of the conserved sequence in sequence-specific termination by T7 RNA polymerase. *Proc. Natl. Acad. Sci. USA* **102** (1):75.

Shen *et al.* (2001) Two site contact of elongating transcripts to phage T7 RNA polymerase at C-terminal regions. *J. Biol. Chem.* **276** (6):4080.

Powell *et al.* (1996) Elongation factor SII contacts the 3'-end of RNA in the RNA polymerase II elongation complex. *J. Biol. Chem.* **1271** (37):22301.

Sheng *et al.* (1993) Wheat germ and yeast RNA polymerase II: photoaffinity labeling by 4-thiouracil 5'-monophosphate positioned uniquely at the 3' end of an enzyme-bound [32P]-containing transcript. *Biochemistry* **32** (9):2248.

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

Kelly *et al.* (1990) Yeast RNA polymerase I. Derivatization of the 190 and 135 subunits by 4-thiouridine monophosphate positioned uniquely at the 3' terminus of an enzyme-bound 32P-containing transcript initiated by a triribonucleotide primer on synthetic single-stranded DNA. *J. Biol. Chem.* **265** (14):7787.

Shostak *et al.* (1990) Direct spectrophotometric assays for orotate phosphoribosyltransferase and orotidylate decarboxylase. *Anal. Biochem.* **191** (2):365.