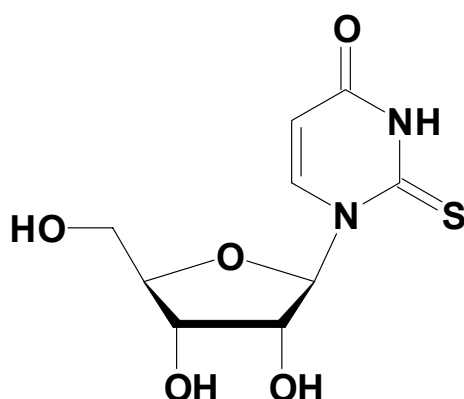


2-Thio-uridine

Cat. No.	Amount
N-1062-20	20 mg
N-1062-100	100 mg

Selected References:

Nawrot *et al.* (1989) The tRNA "WOBBLE POSITION" Uridines. III. 1 the Synthesis of 5-[S-Methoxycarbonyl (Hydroxy)Methyl] Uridine and its 2-Thio Analogue. *Nucleosides, Nucleotides and Nucleic Acids* **8**:1499.



CAS number: 20235-78-3

Molecular formula: C₉H₁₂N₂O₅S

Molecular weight: 260.27 g/mol

Purity: ≥ 98% (TLC)

Appearance: white to off-white powder

Storage conditions: store at -20°C

Shelf life: 12 months

For research use only!