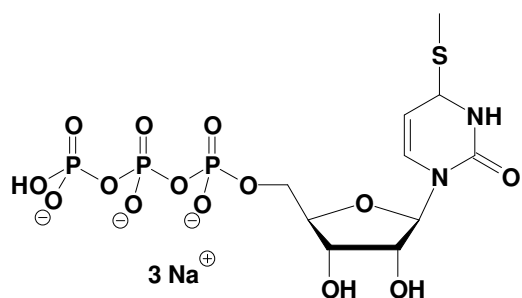


## 4-Thio-methyl-UTP

### 4-Thio-methyl-uridine-5'-triphosphate

Cat. No.	Amount
NU-1187S	20 Units
NU-1187L	100 Units



**Cat. No.:** NU-1187

**Molecular Formula:**  $C_{10}H_{16}N_2O_{14}P_3S$  (Anion)

**Molecular Weight:** 513.22 (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  $\mu$ l of a 10 mM solution

**For research use only!**