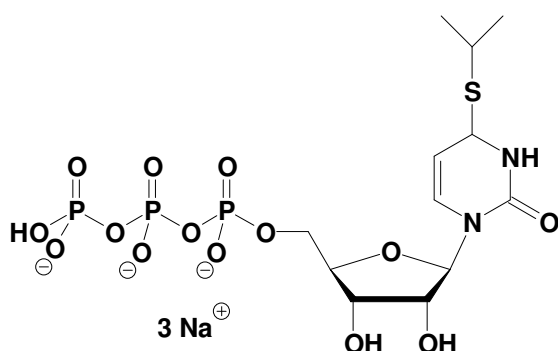


4-Thio-*i*-propyl-UTP

4-Thio-*iso*-propyl-uridine-5'-triphosphate

Cat. No.	Amount
NU-1189S	20 Units
NU-1189L	100 Units



Cat. No.: NU-1189

Molecular Formula: $C_{12}H_{20}N_2O_{14}P_3S$ (Anion)

Molecular Weight: 541.27 (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 μ l of a 10 mM solution

For research use only!