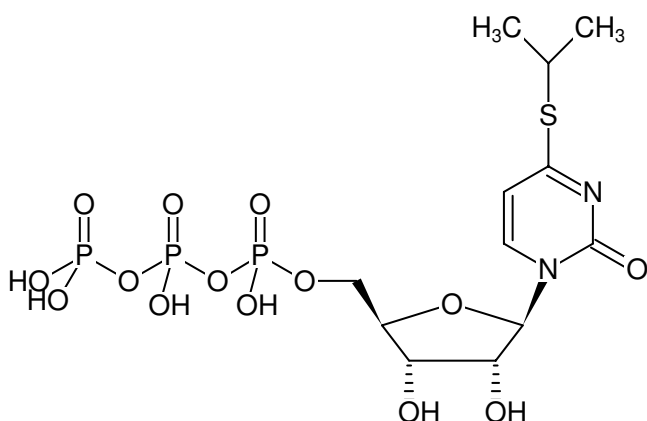




4-Thio-*i*-propyl-UTP

4-Thio-*iso*-propyl-uridine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1189S	20 µl (10 mM)
NU-1189L	5 x 20 µl (10 mM)



Structural formula of 4-Thio-*i*-propyl-UTP

For research use only!

Shipping: shipped on blue ice

Storage Conditions: store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: C₁₂H₂₁N₂O₁₄P₃S (free acid)

Molecular Weight: 542.28 g/mol (free acid)

Exact Mass: 541.99 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: clear aqueous solution

Concentration: 10 mM - 11 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{max} 303 nm, ε 14.1 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)