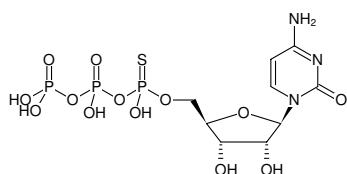


**CTPαS**Cytidine-5'-(α-thio)-triphosphate, Sodium salt; (1 : 1 Mixture of R_p and S_p isomers)

Cat. No.	Amount
NU-436S	25 µl (100 mM)
NU-436L	5 x 25 µl (100 mM)



Structural formula of CTPαS

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:** C₉H₁₆N₃O₁₃P₃S (free acid)**Molecular Weight:** 499.21 g/mol (free acid)**Exact Mass:** 498.96 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 100 mM - 110 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 271 nm, ε 9.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)