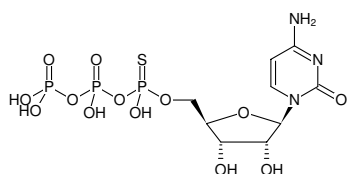


**CTP α S**Cytidine-5'-(α -thio)-triphosphate, Sodium salt; (Mixture of R_p and S_p isomers)

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

Structural formula of CTP α S**For general laboratory use.****Shipping:** shipped on gel packs**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:** \geq 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 100 mM - 110 mM**pH:** 7.5 \pm 0.5**Spectroscopic Properties:** λ_{max} 271 nm, ϵ 9.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)