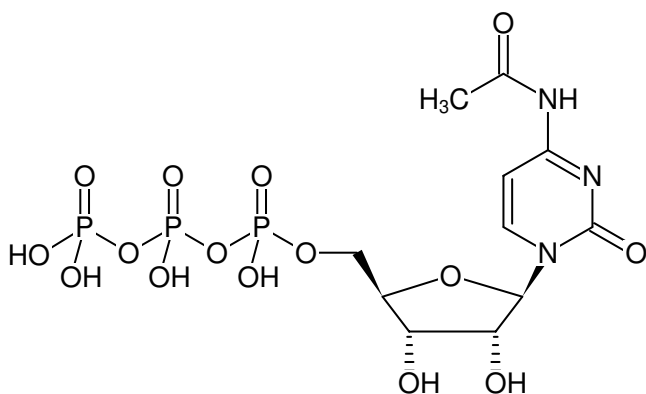


**N⁴-Acetyl-CTP**ac⁴CTPN⁴-Acetyl-cytidine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-988S	10 µl (100 mM)
NU-988L	5 x 10 µl (100 mM)

Related Products:

HighYield T7 RNA Synthesis Kit, #RNT-101

Selected References:Arango *et al.* (2018) Acetylation of Cytidine in mRNA Promotes Translation Efficiency. *Cell* **175** (7):1872.Sinclair *et al.* (2017) Profiling Cytidine Acetylation with Specific Affinity and Reactivity. *ACS Chem. Neurosci.* **12** (12):2922.Structural formula of N⁴-Acetyl-CTP**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:** C₁₁H₁₈N₃O₁₅P₃ (free acid)**Molecular Weight:** 525.19 g/mol (free acid)**Exact Mass:** 525.00 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 100 mM - 110 mM**pH:** 7.0 ±0.5**Spectroscopic Properties:** λ_{max} 243 nm, ε 11.2 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)