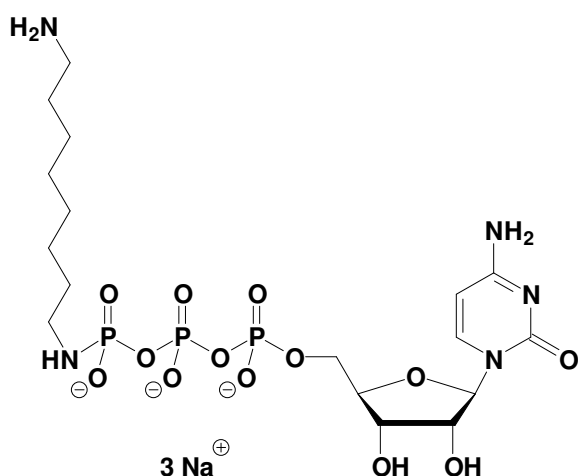


γ -[(8-Aminoethyl)-imido]-CTP

γ -[(8-Aminoethyl)-imido]-Cytosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-844S	5 mg
NU-844L	25 mg



Cat. No.: NU-844

Molecular Formula: C₁₇H₃₁N₅O₁₃P₃ (Anion)

Molecular Weight: 606.38 (Anion)

Purity: > 95%, lyophilized powder

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.

For research use only!