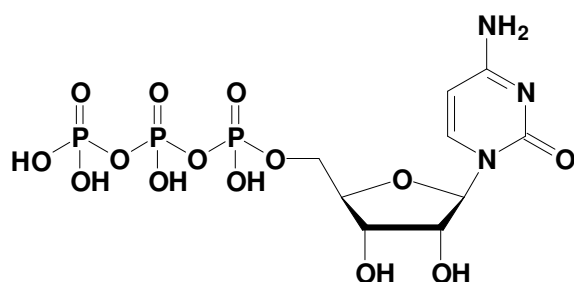


CTP, solution

Cytidine 5'-triphosphate, sodium salt

100 mM

Cat. No.	Amount
NU-1011	1 ml (100 μ mol)



Description

Ultrapure CTP supplied as clear aqueous solution (pH 8.0).

Selected References:

Spangler *et al.* (2011) Interaction of the diguanylate cyclase YdeH of *Escherichia coli* with 2',(3')-substituted purine and pyrimidine nucleotides. *J. Pharmacol. Exp. Ther.* **336** (1):234.

For *in vitro* use only

Quality guaranteed for 12 months

Store at -20°C, short term (up to one week) exposure to ambient temperature possible

Concentration

100 mM +/-2%

Form

clear aqueous solution, pH 8.0 +/-0.2 (4°C)

Purity

>99%

Molecular Formula

C₉H₁₃N₃O₁₄P₃ (Anion)

Molecular Weight

480.16 (Anion)

Absorbance

absorbance max: 271 nm (pH 7)

ϵ at absorbance max: 8.9 mmol⁻¹ cm⁻¹

Quality Control Specifications

in vitro transcription:

suitable

contamination with bacterial and human DNA:

not detectable

activity of DNase, Protease or Phosphatase:

not detectable