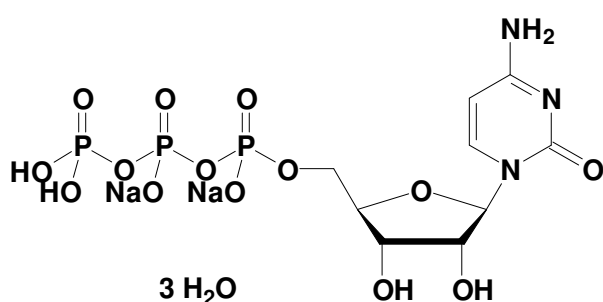


CTP, lyophilized Cytidine 5'-triphosphate, sodium salt

Cat. No.	Amount
NU-1011-100	100 mg, lyophilized
NU-1011-1G	1 g, lyophilized
NU-1011-10G	10 g, lyophilized
NU-1011-100G	100 g, lyophilized



Molecular Formula: $C_9H_{14}N_3Na_2O_{14}P_3 \times 3H_2O$

Molecular Weight: 581.16

Absorbance

absorbance max: 271 nm (pH 7)

ϵ at absorbance max: $8.9 \text{ mmol}^{-1} \text{ cm}^{-1}$

Purity: >95%

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at $< -20^\circ\text{C}$.

For research use only!

Applications:

Physiological role in coronary artery disease^[1]

Physiological role in lipid metabolism^[2]

Physiological role in farnesol induced apoptosis^[3]

Specific Ligands:

CTP synthase^[4]

Phosphocholine cytidyltransferase alpha^[2]

Selected References:

[1] Lui *et al.* (2010) Evaluation of CT perfusion in setting of cerebral ischemia: patterns and pitfalls. *American Journal of Neuroradiology* **31**:1552.

[2] Luoma (2010) Gene activation regresses arteriosclerosis, promotes health, and enhances longevity. *Lipids in health and disease* **9**:67.

[3] Joo *et al.* (2010) Molecular mechanisms involved in farnesol-induced apoptosis. *Cancer letters* **287**:123.

[4] Cabeen *et al.* (2010) A metabolic assembly line in bacteria. *Nature Cell Biology* **12**:731.

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.