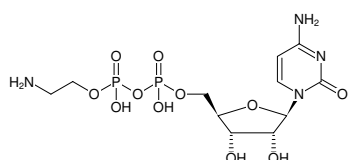




CDP-ethanolamine

Cytidine-5'-diphosphate ethanolamine, free acid
Cytidine-5'-diphosphateethanolamine

Cat. No.	Amount
NU-1208-5	5 mg
NU-1208-50	50 mg



Structural formula of CDP-ethanolamine

For research use only!

Shipping: shipped on blue ice

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: C₁₁H₂₀N₄O₁₁P₂ (free acid)

Molecular Weight: 446.24 g/mol (free acid)

Exact Mass: 446.06 g/mol (free acid)

CAS#: 3036-18-8

Purity: ≥ 95 % (HPLC)

Form: white powder

Spectroscopic Properties: λ_{max} 271 nm, ε 8.9 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Selected References:

Martin *et al.* (1993) 31P-NMR analysis of *Entamoeba histolytica*. Occurrence of high amounts of two inositol phosphates. *Eur. J. Biochem.* **214** (3):711.

Bostwick *et al.* (1989) Purification and characterization of a central cholinergic enhancing factor from rat brain: its identity as phosphoethanolamine. *J. Neurochem.* **53** (2):448.

Zelinsky *et al.* (1982) Phosphatidylethanolamine biosynthesis in isolated hamster heart. *Can. J. Biochem.* **60** (8):817.

Zelinsky *et al.* (1982) Choline regulates phosphatidylethanolamine biosynthesis in isolated hamster heart. *J. Biol. Chem.* **257** (22):13201.

Brunetti *et al.* (1979) Synthesis of phosphatidylcholine and phosphatidylethanolamine at different ages in the rat brain in vitro. *Lipids.* **14** (11):925.

Call *et al.* (1975) Synthesis of ethanolamine phosphoglycerides by human platelets. *J. Lipid Res.* **16** (5):352.

Roberti *et al.* (1975) Enzymic synthesis of 1-alkyl-2-acyl-sn-glycero-3-phosphorylethanolamine through ethanolaminephosphotransferase activity in the neuronal and glial cells of rabbit in vitro. *Lipids.* **10** (3):121.

Chojnacki *et al.* (1966) Enzymatic synthesis of labelled cytidine-5'-diphosphate ethanolamine. *Nature.* **210** (5039):947.

Kennedy *et al.* (1956) The function of cytidine coenzymes in the biosynthesis of phospholipides. *J. Biol. Chem.* **222** (1):193.