

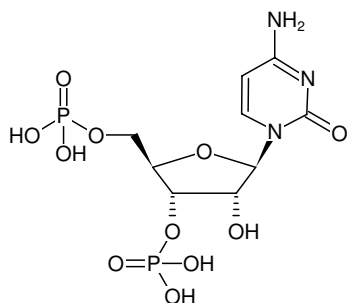


Cytidine-3',5'-bisphosphate

3',5'-pCp

Cytidine-3',5'-bisphosphate, Sodium salt

Cat. No.	Amount
NU-549S	100 µl (10 mM)
NU-549L	5 x 100 µl (10 mM)



Structural formula of Cytidine-3',5'-bisphosphate

For general laboratory use.

Please centrifuge briefly before opening (volume ≤2 ml).

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: C₉H₁₅N₃O₁₁P₂ (free acid)

Molecular Weight: 403.18 g/mol (free acid)

Exact Mass: 403.02 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 10 mM - 11 mM

pH: 7.5 ±0.5

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