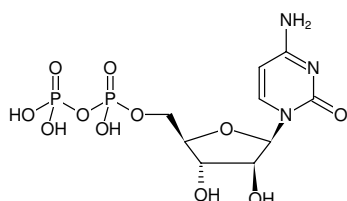



ara-Cytidine-5'-diphosphate (ara-CDP)

Cytarabine diphosphate
Cytidine-arabinofuranoside-5'-diphosphate, Sodium salt

Cat. No.	Amount
NU-1171S	50 µl (10 mM)
NU-1171L	5 x 50 µl (10 mM)



Structural formula of *ara*-Cytidine-5'-diphosphate (*ara*-CDP)

For general laboratory use.

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: 403.17 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 7.0)