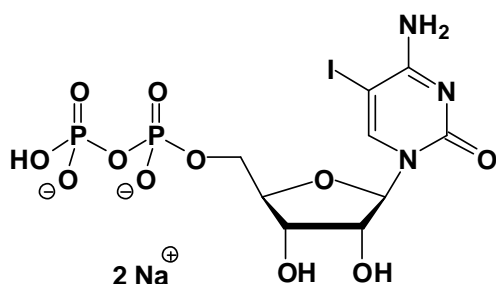


5I-CDP

5-Iodo-cytidine-5'-diphosphate, Sodium salt

Cat. No.	Amount
NU-138S	30 Units
NU-138L	150 Units



Cat. No.: NU-138

Molecular Formula: C₉H₁₂N₃O₁₁P₂I (Anion)

Molecular Weight: 527.05 (Anion)

Purity: > 95%, clear aqueous solution, pH 7.5

Spectroscopic properties: λ_{max} 293 nm; ε 5700

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

For research use only!

1 unit = 1 μl of a 10 mM solution