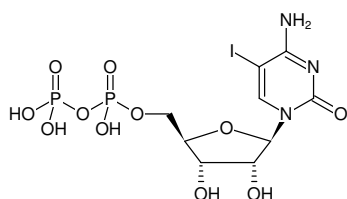


**5-Iodo-CDP**

(5I-CDP)

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

Cat. No.	Amount
NU-138S	30 µl (10 mM)
NU-138L	5 x 30 µl (10 mM)



Structural formula of 5-Iodo-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:** C₉H₁₄N₃O₁₁P₂I (free acid)**Molecular Weight:** 529.07 g/mol (free acid)**Exact Mass:** 528.91 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} 293 nm, ε 5.7 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)