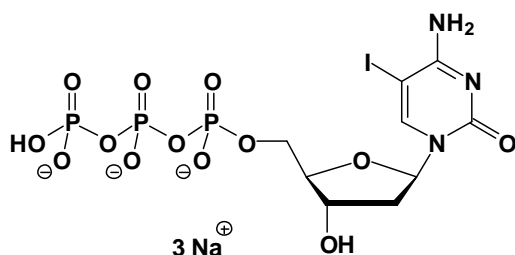


5I-dCTP

5-Iodo-2'-deoxy-cytidine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-128S	50 Units
NU-128L	250 Units



Cat. No.: NU-128

Molecular Formula: C₉H₁₂N₃O₁₃P₃I (Anion)

Molecular Weight: 590.03 (Anion)

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

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