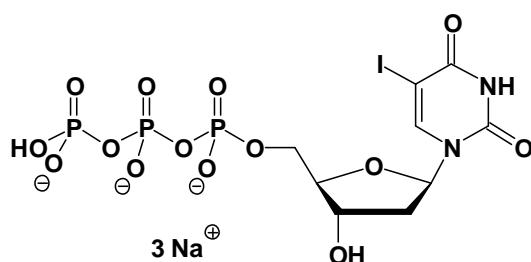


5I-dUTP

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

Cat. No.	Amount
NU-120S	50 Units
NU-120L	250 Units



Cat. No.: NU-120

Molecular Formula: C₉H₁₁N₂O₁₄P₃I (Anion)

Molecular Weight: 591.01 (Anion)

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

Spectroscopic properties:

$\epsilon_{\text{max}} = 7700 \text{ M}^{-1} \text{ cm}^{-1}$ at 287 nm (Tris-HCl 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