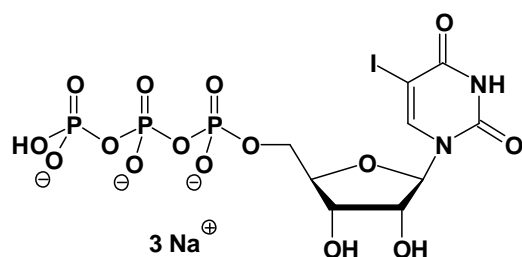


5I-UTP

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

Cat. No.	Amount
NU-119S	50 Units
NU-119L	250 Units



Selected References:

Gilles *et al.* (2007) Regulatory Mechanism Differ in UMP Kinase from Gram-negative and Gram-positive Bacteria. *J. Biol. Chem.* **282** (10):7242.

Cat. No.: NU-119

Molecular Formula: C₉H₁₁N₂O₁₅P₃I (Anion)

Molecular Weight: 607.01 (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