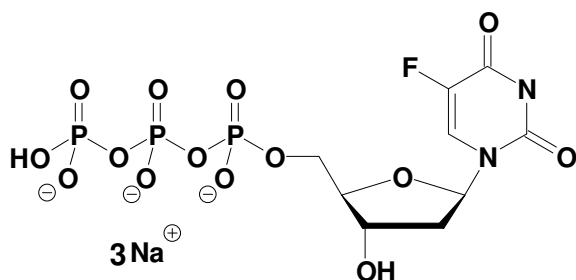


5F-dUTP

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

Cat. No.	Amount
NU-154S	50 Units
NU-154L	250 Units



Cat. No.: NU-154

Molecular Formula: C₉H₁₀N₂O₁₄P₃F (Anion)

Molecular Weight: 482.10 (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