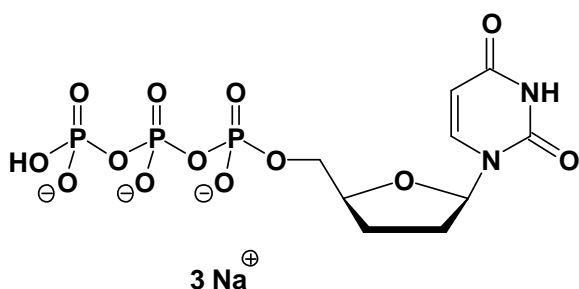


ddUTP

2',3'-Dideoxy-uridine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1021S	100 Units
NU-1021L	500 Units



Cat. No.: NU-1021

Molecular Formula: C₉H₁₂N₂O₁₃P₃ (Anion)

Molecular Weight: 449.12 (Anion)

Purity: at least 99%, clear aqueous solution, pH 8,4 - 8,6

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C.

For research use only!

1 unit = 1 μl of a 10 mM solution