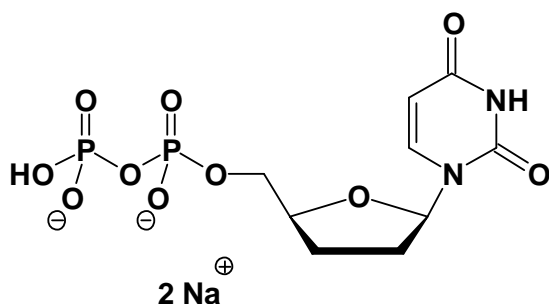


ddUDP

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

Cat. No.	Amount
NU-1022S	50 Units
NU-1022L	250 Units



Cat. No.: NU-1022

Molecular Formula: C₉H₁₂N₂O₁₀P₂ (Anion)

Molecular Weight: 370.14 (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