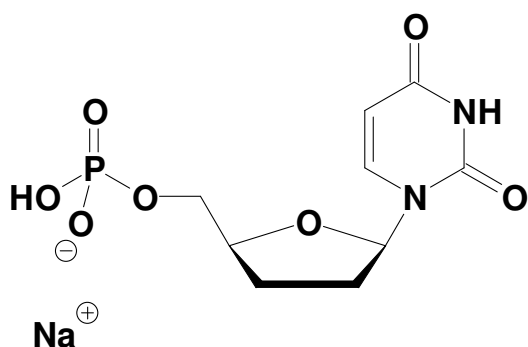


ddUMP

2',3'-Dideoxy-Uridine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-1169S	50 Units
NU-1169L	250 Units



Cat. No.: NU-1169

Molecular Formula: C₉H₁₂N₂O₇P (Anion)

Molecular Weight: 291.17 (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.

* 1 unit = 1 µl of a 10 mM solution

For research use only!