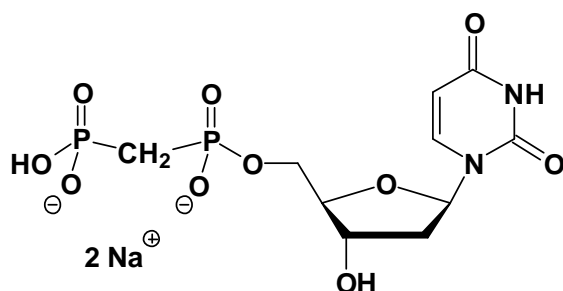


dUpCp (dUMPCP)

2'-Deoxy-uridine-5'-[(α,β)-methyleno]diphosphate, Sodium salt

Cat. No.	Amount
NU-428S	100 Units
NU-428L	500 Units



Cat. No.: NU-428

Molecular Formula: C₁₀H₁₄N₂O₁₀P₂ (Anion)

Molecular Weight: 384.17 (Anion)

Purity: > 95%, clear aqueous solution, pH 7.5

Spectroscopic properties: λ_{\max} 262 nm; ϵ 10200

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 μ l of a 10 mM solution