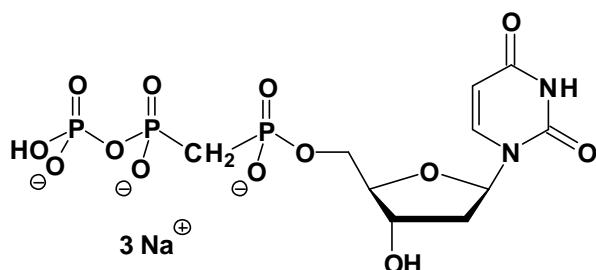


dUpCpp (dUMPCPP)

2'-Deoxy-uridine-5'-[(α,β)-methylene]triphosphate, Sodium salt

Cat. No.	Amount
NU-429S	50 Units
NU-429L	250 Units



Cat. No.: NU-429

Molecular Formula: $C_{10}H_{14}N_2O_{13}P_3$ (Anion)

Molecular Weight: 463.14 (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 6 months after date of delivery.

For research use only!

1 unit = 1 μ l of a 10 mM solution