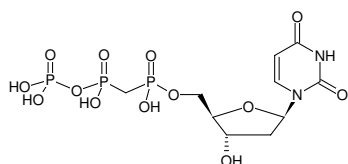


**dUpCpp**

(dUMPCPP)

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

Cat. No.	Amount
NU-429S	50 μ l (10 mM)
NU-429L	5 x 50 μ l (10 mM)



Structural formula of dUpCpp

For research use only!**Shipping:** shipped on blue ice**Storage Conditions:** store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery**Molecular Formula:** C₁₀H₁₇N₂O₁₃P₃ (free acid)**Molecular Weight:** 466.17 g/mol (free acid)**Exact Mass:** 465.99 g/mol (free acid)**CAS#:** 170428-84-9**Purity:** \geq 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 \pm 0.5**Spectroscopic Properties:** λ_{\max} 262 nm, ϵ 10.2 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Applications:**Catalytically active complex^[1]**Selected References:**

[1] Guranowski *et al.* (2006) Methylene analogues of adenosine 5-tetraphosphate. Their chemical synthesis and recognition by human and plant mononucleoside tetraphosphates and dinucleotide tetraphosphates. *FEBS Journal* **273**:829.

Xia *et al.* (2011) Structural insights into complete metal ion coordination from ternary complexes of B family RB69 DNA polymerase. *Biochemistry*. **50** (42): 9114.