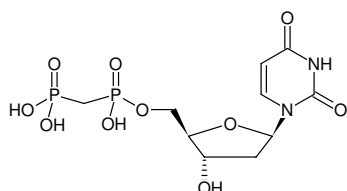


**dUpCp**

(dUMPCP)

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

Cat. No.	Amount
NU-428S	100 μ l (10 mM)
NU-428L	5 x 100 μ l (10 mM)



Structural formula of dUpCp

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:** C₁₀H₁₆N₂O₁₀P₂ (free acid)**Molecular Weight:** 386.19 g/mol (free acid)**Exact Mass:** 386.03 g/mol (free acid)**Purity:** \geq 95 % (HPLC)**Form:** clear aqueous solution**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)