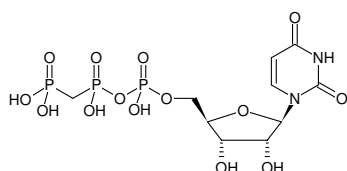


**UppCp**

(UMPPCP)

Uridine-5'-[(β,γ)-methylene]triphosphate, Sodium salt

Cat. No.	Amount
NU-255	5 mg



Structural formula of UppCp

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:** C₁₀H₁₇N₂O₁₄P₃ (free acid)**Molecular Weight:** 482.17 g/mol (free acid)**Exact Mass:** 481.99 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** white to off-white solid**Spectroscopic Properties:** λ_{\max} 262 nm, ϵ 10.2 L mmol⁻¹ cm⁻¹
(Tris-HCl pH 7.5)