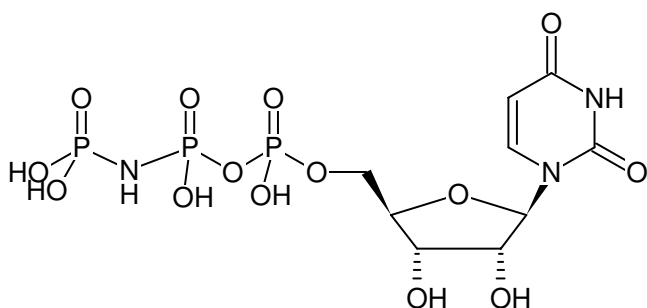


**UppNHp - Solid**

(UMPPNP)

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

Cat. No.	Amount
NU-415-5	5 mg



Structural formula of UppNHp - Solid

For general laboratory use.**Shipping:** shipped on gel packs**Storage Conditions:** store at -20 °C

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

Shelf Life: 6 months after date of delivery**Molecular Formula:** C₉H₁₆N₃O₁₄P₃ (free acid)**Molecular Weight:** 483.15 g/mol (free acid)**Exact Mass:** 482.98 g/mol (free acid)**CAS#:** 82145-58-2**Purity:** ≥ 95 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ_{\max} 262 nm, ϵ 10.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)