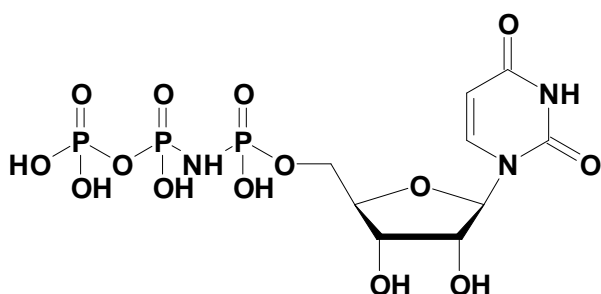


UpNHpp (UMPNPP)

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

Cat. No.	Amount
NU-930S	50 Units
NU-930L	250 Units



Cat. No.: NU-930

Molecular Formula: $C_9H_{16}N_3O_{14}P_3$ (free acid)

Molecular Weight: 483.15 (free acid)

Purity: > 95%, clear aqueous solution, pH 7.5

Spectroscopic Properties: λ_{max} 262nm;
 ϵ 10,200 $M^{-1}cm^{-1}$ Tris HCl pH 7.5

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