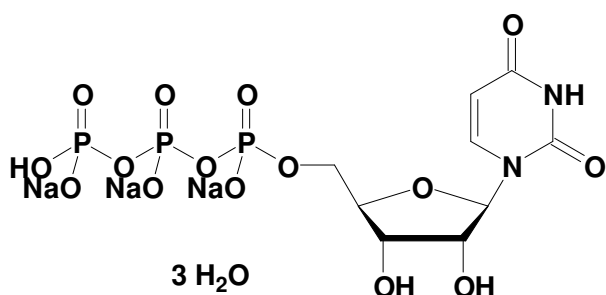


UTP, lyophilized

Uridine 5'-triphosphate, sodium salt

Cat. No.	Amount
NU-1013-100	100 mg, lyophilized
NU-1013-1G	1 g, lyophilized
NU-1013-10G	10 g, lyophilized
NU-1013-100G	100 g, lyophilized



Molecular Formula: C₉H₁₂N₂Na₃O₁₅P₃ × 3H₂O

Molecular Weight: 604.12

Absorbance

absorbance max: 262 nm (pH 7)

ε at absorbance max: 9.8 mmol⁻¹ cm⁻¹

Purity: >90%

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C.

For research use only!