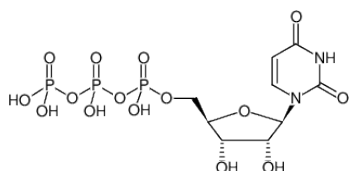


**UTP - Solid**

Uridine 5'-triphosphate, Trisodium salt trihydrate

Cat. No.	Amount
NU-1051-100	100 mg
NU-1051-100G	100 g
NU-1051-10G	10 g
NU-1051-1G	1 g

Quality Control Specifications: in vitro transcription (T7 RNA polymerase): visible RNA fragments after 5 min incubation, Proteases: not detectable



Structural formula of UTP - Solid

For general laboratory use.**Shipping:** Shipped on gel packs**Storage Conditions:** Store at -20 °C**Shelf Life:** 12 months**Molecular Formula:** C₉H₁₅N₂O₁₅P₃ (free acid)**Molecular Weight:** 484.14 g/mol (free acid)**CAS#:** 19817-92-6**Purity:** ≥ 90 % (HPLC)**Form:** Powder**Color:** White**Solubility:** Soluble in water. Clear solution.**Spectroscopic Properties:** λ_{max} 262 nm, ε 9.8 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.0)