

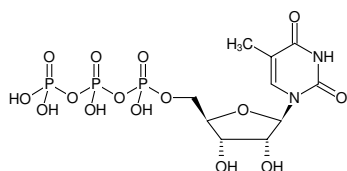
**5-Methyl-UTP**

m5UTP

5-Me-UTP

5-Methyluridine-5'-O-triphosphate, Sodium salt

Cat. No.	Amount
NU-880S	10 µl (100 mM)
NU-880L	5 x 10 µl (100 mM)



Structural formula of 5-Methyl-UTP

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: 12 months after date of delivery**Molecular Formula:** C₁₀H₁₇N₂O₁₅P₃ (free acid)**Molecular Weight:** 498.17 g/mol (free acid)**Exact Mass:** 497.98 g/mol (free acid)**CAS#:** 23198-01-8 (free acid), 1801968-79-5 (Tetrasodium salt)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 100 mM - 110 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 267 nm, ε 9.65 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)