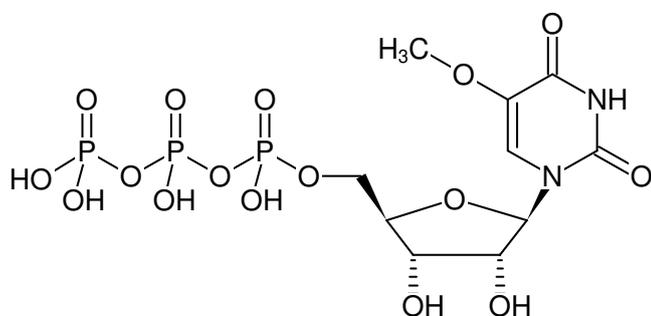


**5-Methoxy-UTP**

5moUTP

5-Methoxyuridine-5'-triphosphate, Sodium Salt

Cat. No.	Amount
NU-972S	10 µl (100 mM)
NU-972L	5 x 10 µl (100 mM)



Structural formula of 5-Methoxy-UTP

For research use only!**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:** C₁₀H₁₇N₂O₁₆P₃ (free acid)**Molecular Weight:** 514.17 g/mol (free acid)**Exact Mass:** 513.98 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 100 mM - 110 mM**pH:** 7.5 ± 0.5**Solubility:** water**Spectroscopic Properties:** λ_{max} 278 nm, ε 8.3 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Related Products:**

HighYield T7 RNA Synthesis Kit, #RNT-101

Selected References:An *et al.* (2019) Long-term efficacy and safety of mRNA therapy in two murine models of methylmalonic acidemia. *EBioMedicine* **45**:519.Novakowski *et al.* (2019) Delivery of mRNA to platelets using lipid nanoparticles. *Sci. Rep.* **9** (1):552.An *et al.* (2017) Systemic Messenger RNA Therapy as a Treatment for Methylmalonic Acidemia. *Cell Rep.* **21** (12):3548.Li *et al.* (2016) Effects of Chemically Modified Messenger RNA on Protein Expression *Bioconjugate Chem.* **27**:849.