

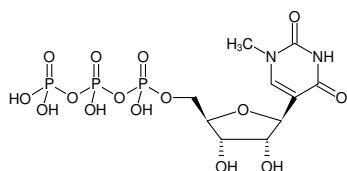


N¹-Methylpseudo-UTP

me¹ψ-UTP

N¹-Methylpseudouridine-5'-triphosphate, Sodium Salt

Cat. No.	Amount
NU-890S	10 µl (100 mM)
NU-890L	5 x 10 µl (100 mM)



Structural formula of N¹-Methylpseudo-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#: 1428903-59-6 (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} 271 nm, ε 8.9 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Related Products:

HighYield T7 RNA Synthesis Kit, #RNT-101

Selected References:

Andies *et al.* (2015) N (1)-methylpseudouridine-incorporated mRNA outperforms pseudouridine-incorporated mRNA by providing enhanced protein expression and reduced immunogenicity in mammalian cell lines and mice. *J. Control. Release* **217**:337.