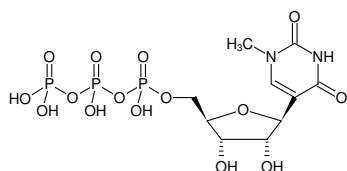


**N¹-Methylpseudo-UTP**me¹ψ-UTPN¹-Methylpseudouridine-5'-triphosphate, Sodium Salt

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

Structural formula of N¹-Methylpseudo-UTP**For research use only!****Shipping:** shipped on blue ice**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.