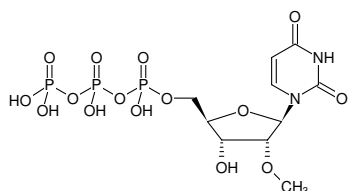


**2'OMe-UTP**

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

Cat. No.	Amount
NU-1212S	50 µl (100 mM)
NU-1212L	5 x 50 µl (100 mM)



Structural formula of 2'OMe-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:** C<sub>10</sub>H<sub>17</sub>N<sub>2</sub>O<sub>15</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 498.17 g/mol (free acid)**Exact Mass:** 497.98 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**Spectroscopic Properties:** λ<sub>max</sub> 262 nm, ε 10.1 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)**Selected References:**Lauridsen *et al.* (2012) Enzymatic recognition of 2'-modified ribonucleoside 5'-triphosphates: towards the evolution of versatile aptamers. *ChemBiochem.* **13** (1):19.Xiao *et al.* (2012) Engineering of Targeted Nanoparticles for Cancer Therapy Using Internalizing Aptamers Isolated by Cell-Uptake Selection. *ACS Nano* **6** (1):696.Keefe *et al.* (2008) SELEX with modified nucleotides. *Curr. Opin. Chem. Biol.* **12** (4):448.Burmeister *et al.* (2006) 2'-Deoxy purine, 2'-O-methyl pyrimidine (dRmY) aptamers as candidate therapeutics. *Oligonucleotides* **16** (4):337.Burmeister *et al.* (2005) Direct in vitro selection of a 2'-O-methyl aptamer to VEGF. *Chem. Biol.* **12** (1):25.