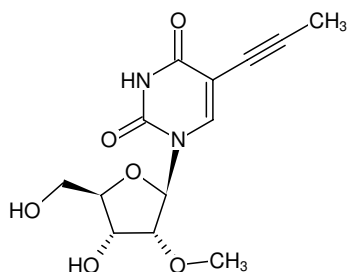


**2'-(O-Methyl)-5-propynyl-uridine**

Cat. No.	Amount
N-RP-1856-10	10 mg
N-RP-1856-100	100 mg



Structural formula of 2'-(O-Methyl)-5-propynyl-uridine

**For research use only!****Shipping:** shipped at ambient temperature**Storage Conditions:** store at -20 °C**Shelf Life:** 30 months after date of delivery**Molecular Formula:** C<sub>13</sub>H<sub>16</sub>N<sub>2</sub>O<sub>6</sub>**Molecular Weight:** 296.28 g/mol**CAS#:** 179817-95-9**Purity:** ≥ 98 % (HPLC)**Form:** white crystalline powder**Solubility:** methanol