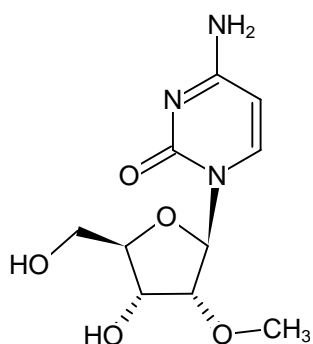


**2'-(O-Methyl)-cytidine**

Cat. No.	Amount
N-RP-1902-1G	1 g
N-RP-1902-10G	10 g



Structural formula of 2'-(O-Methyl)-cytidine

**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>10</sub>H<sub>15</sub>N<sub>3</sub>O<sub>5</sub>**Molecular Weight:** 257.24 g/mol**Purity:** ≥ 98 % (HPLC)**Form:** white crystalline powder**Solubility:** methanol