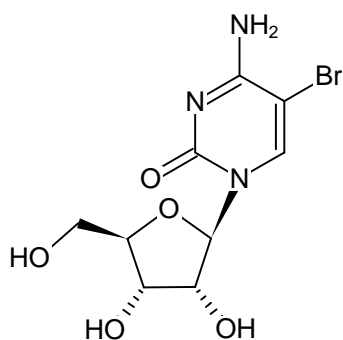


**5-Bromo-cytidine**

Cat. No.	Amount
N-RP-1215-1G	1 g
N-RP-1215-5G	5 g



Structural formula of 5-Bromo-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:** C₉H₁₂BrN₃O₅**Molecular Weight:** 322.11 g/mol**Purity:** ≥ 98 % (HPLC)**Form:** white crystalline powder**Solubility:** Water