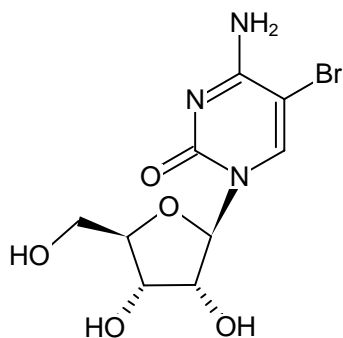


**5-Bromo-cytidine**

| 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