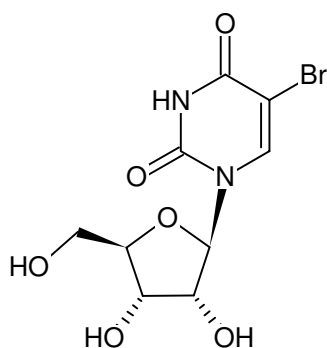


**5-Bromo-uridine (5-BrU)**

Cat. No.	Amount
N-RP-1192-1G	1 g



Structural formula of 5-Bromo-uridine (5-BrU)

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:** C₉H₁₁BrN₂O₆**Molecular Weight:** 323.10 g/mol**CAS#:** 957-75-5**Purity:** ≥ 98 % (HPLC)**Form:** white crystalline powder**Solubility:** water