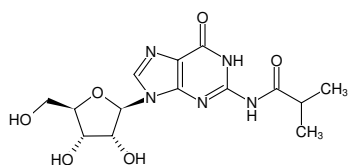


**Guanosine (N-iBu)**

Cat. No.	Amount
N-PM-2103-5G	5 g



Structural formula of Guanosine (N-iBu)

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:** C₁₄H₁₉N₅O₆**Molecular Weight:** 353.33 g/mol**CAS#:** 64350-24-9**Purity:** ≥ 98 % (HPLC)**Form:** white crystalline powder**Solubility:** methanol