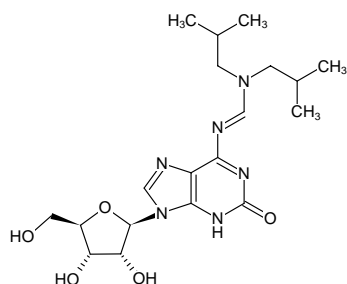


**Isoguanosine (N/N-diisobutylformamide)**

Cat. No.	Amount
N-PM-2643-100	100 mg
N-PM-2643-250	250 mg



Structural formula of Isoguanosine (N/N-diisobutylformamide)

For research use only!**Shipping:** shipped at ambient temperature**Storage Conditions:** store at -20 °C**Shelf Life:** 30 months**Molecular Formula:** C₁₉H₃₀N₆O₅**Molecular Weight:** 422.48 g/mol**Purity:** ≥ 98 % (HPLC)**Form:** white crystalline powder**Solubility:** methanol