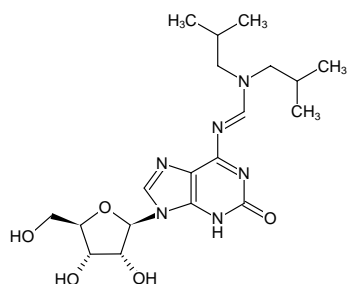


**Isoguanosine (N/N-diisobutylformamide)**

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 after date of delivery**Molecular Formula:** C<sub>19</sub>H<sub>30</sub>N<sub>6</sub>O<sub>5</sub>**Molecular Weight:** 422.48 g/mol**Purity:** ≥ 98 % (HPLC)**Form:** white crystalline powder**Solubility:** methanol