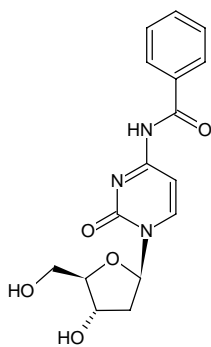


**2'-Deoxycytidine (N-Bz)**

Cat. No.	Amount
N-PM-1102-5G	5 g
N-PM-1102-10G	10 g



Structural formula of 2'-Deoxycytidine (N-Bz)

**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>16</sub>H<sub>17</sub>N<sub>3</sub>O<sub>5</sub>**Molecular Weight:** 331.32 g/mol**Purity:** ≥ 98 % (HPLC)**Form:** white crystalline powder**Solubility:** methanol