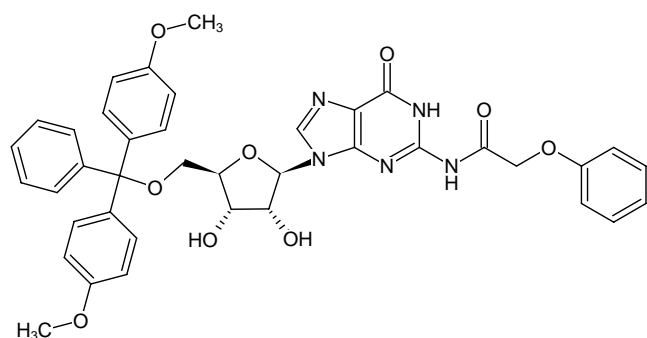


**5'-DMT-guanosine (N-PAC)**

Cat. No.	Amount
N-PM-2205-5G	5 g
N-PM-2205-10G	10 g



Structural formula of 5'-DMT-guanosine (N-PAC)

For research use only!**Shipping:** shipped at ambient temperature**Storage Conditions:** store at -20 °C**Shelf Life:** 30 months**Molecular Formula:** C₃₉H₃₇N₅O₉**Molecular Weight:** 719.74 g/mol**Purity:** ≥ 98 % (HPLC)**Form:** white crystalline powder**Solubility:** acetonitrile