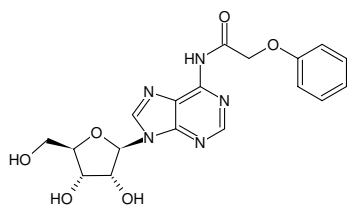


**Adenosine (N-PAC)**

Cat. No.	Amount
N-PM-6001-1G	1 g
N-PM-6001-5G	5 g



Structural formula of Adenosine (N-PAC)

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:** C₁₈H₁₉N₅O₆**Molecular Weight:** 401.37 g/mol**CAS#:** 119824-65-6**Purity:** ≥ 98 % (HPLC)**Form:** white crystalline powder**Solubility:** methanol