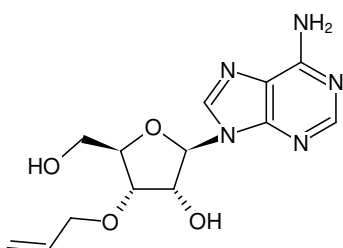


**3'-(O-Allyl)-adenosine**

Cat. No.	Amount
N-RP-3111-10	10 mg
N-RP-3111-100	100 mg



Structural formula of 3'-(O-Allyl)-adenosine

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:** C₁₃H₁₇N₅O₄**Molecular Weight:** 307.31 g/mol**CAS#:** 400820-99-7**Purity:** ≥ 98 % (HPLC)**Form:** white crystalline powder**Solubility:** methanol