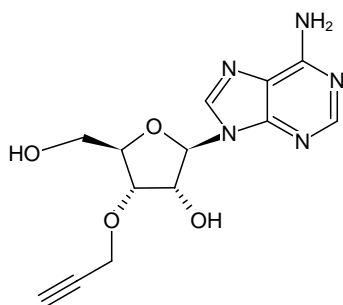


**3'-(O-Propargyl)-adenosine**

Cat. No.	Amount
CLK-RP-3301-10	10 mg
CLK-RP-3301-100	100 mg



Structural formula of 3'-(O-Propargyl)-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:** C₁₃H₁₅N₅O₄**Molecular Weight:** 305.29 g/mol**Purity:** > 98 % (HPLC)**Form:** white crystalline powder**Solubility:** methanol