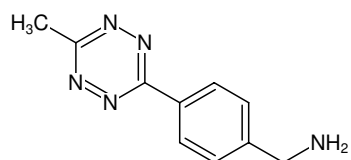


**6-Methyl-Tetrazine-Amine (HCl salt)**

(4-(6-Methyl-1,2,4,5-tetrazin-3-yl)phenyl)methanamine, HCl

Cat. No.	Amount
CLK-1011-25	25 mg
CLK-1011-100	100 mg



Structural formula of 6-Methyl-Tetrazine-Amine (HCl salt)

For research use only!**Shipping:** shipped at ambient temperature**Storage Conditions:** store at -20 °C**Shelf Life:** 12 months after date of delivery**Molecular Formula:** C₁₀H₁₁N₅**Molecular Weight:** 201.23 g/mol**CAS#:** 1345955-28-3**Purity:** > 90 % (HPLC)**Form:** solid**Color:** pink**Solubility:** DMSO, DMF