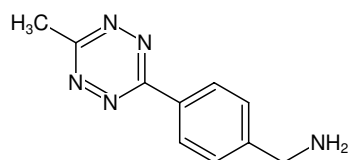


**6-Methyl-Tetrazine-Amine (HCl salt)**

(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:** C<sub>10</sub>H<sub>11</sub>N<sub>5</sub>**Molecular Weight:** 201.23 g/mol**CAS#:** 1345955-28-3**Purity:** > 95 % (HPLC)**Form:** pink solid**Solubility:** DMSO, DMF