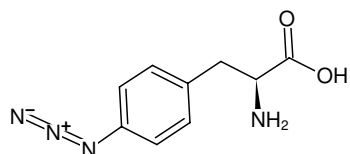


**4-Azido-L-phenylalanine***p*-Azido-Phenylalanine

Cat. No.	Amount
CLK-AA001-10	10 mg
CLK-AA001-100	100 mg
CLK-AA001-500	500 mg



Structural formula of 4-Azido-L-phenylalanine

For general laboratory use.**Shipping:** shipped on gel packs**Storage Conditions:** store at -20 °C**Additional Storage Conditions:** store dry**Shelf Life:** 12 months after date of delivery**Molecular Formula:** C₉H₁₀N₄O₂**Molecular Weight:** 206.20 g/mol**Exact Mass:** 206.08 g/mol**CAS#:** 33173-53-4**Purity:** ≥ 98 % (HPLC)**Form:** powder**Color:** white to off-white