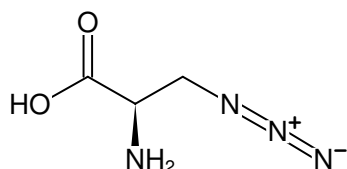


**3-Azido-D-alanine HCl**

(R)-2-Amino-3-azidopropanoic acid hydrochloride

Cat. No.	Amount
CLK-AA004-10	10 mg
CLK-AA004-100	100 mg
CLK-AA004-500	500 mg



Structural formula of 3-Azido-D-alanine HCl

For general laboratory use.**Shipping:** shipped on gel packs**Storage Conditions:** store at 4 °C**Additional Storage Conditions:** store dry**Shelf Life:** 12 months after date of delivery**Molecular Formula:** C₃H₆N₄O₂**Molecular Weight:** 130.11 g/mol**Exact Mass:** 130.05 g/mol**CAS#:** 105928-88-9**Purity:** mass identification conforms (ESI-MS)**Form:** powder**Color:** white