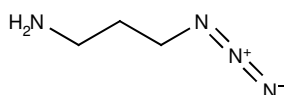


Azido-Propylamine

Cat. No.	Amount
CLK-AZ115-100	100 mg
CLK-AZ115-1000	1 g



Molecular Formula: C₃H₈N₄

Molecular Weight: 100.12 g/mol

Storage conditions: store at -20°C

Shelf life: 12 months

Purity: >95%

Appearance: Slightly yellow oil

Solubility: DMSO, DMF, DCM, Chloroform, THF, Water

Product Features and Benefits:

Azido-containing reagent with carbonyl/carboxyl reactivity.

For research use only!