

N-BOC-ethylenediamine *CAS 57260-73-8*

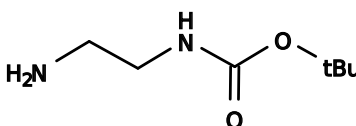
Catalog Number: 4013 4014,
Unit Size: 5 g 25 g

Product Details

Storage Conditions	Freeze (< -15 °C),
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Liquid colorless
Molecular Weight	160.21
Soluble In	DMF
Chemical Structure	



Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

N-BOC-ethylenediamine is a monoprotected diamine that is widely used as a building block in bioconjugate chemistry. It has been used in both oligonucleotide and peptide synthesis.