

N-FMOC-ethylenediamine hydrochloride *CAS 166410-32-8*

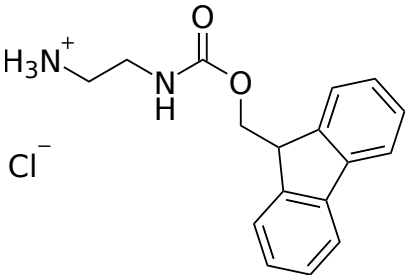
Catalog Number: 4006, 4007

Unit Size: 5 g, 25 g

Product Details

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

Chemical Properties

Appearance	Solid white
Molecular Weight	318.80
Soluble In	DMF
Chemical Structure	

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

N-FMOC-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.