

N-FMOC-ethylenediamine

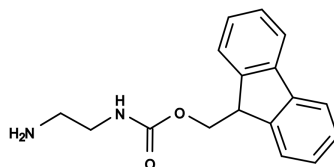
Catalog number: 4016, 4017
Unit size: 5 g, 25 g

Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	282.34
Soluble In	DMF
Chemical Structure	



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.