

Product Information Sheet

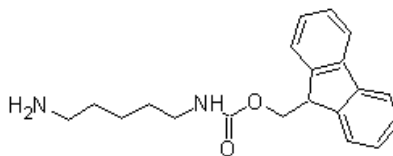
Ordering Information

Product Number: 4019
Product Name: N-FMOC-cadaverine
Unit Size: 25 g
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 324.42

Chemical Structure:



Soluble in: DMF

Excitation Wavelength: NA

Emission Wavelength: NA

Application Notes

N-FMOC-cadaverine is a monoprotected diamine that is widely used as a building block in bioconjugate chemistry.