

N-BOC-cadaverine *CAS 51644-96-3*

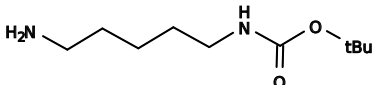
Catalog Number: 4010, 4011

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	202.29
Soluble In	DMF
Chemical Structure	

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

N-BOC-cadaverine 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.