

5(6)-FAM cadaverine

PRODUCT INFORMATION SHEET

Catalog number: 127 Unit size: 100 mg

Product Details		
Storage Conditions	Freeze (<-15 °C), Minimize light exposure	
Expiration Date	12 months upon receiving	
Chemical Properties		
Appearance	Orange solid	
Molecular Weight	574.50	
Soluble In	DMSO	
Chemical Structure HO + CHORE + O + O + O + O + O + O + O + O + O +		
Spectral Properties		

Excitation Wavelength	493 nm
Emission Wavelength	517 nm

Applications

FAM cadaverine is an excellent building block for developing fluoresceinated bioconjugates. It is also a fluorescent transglutaminase substrate to label proteins by transaminidation.