

mBCl [Monochlorobimane] *CAS 76421-73-3*

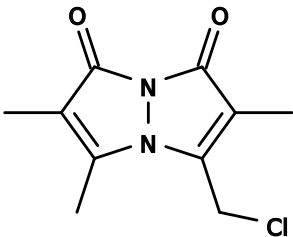
Catalog Number: 635

Unit Size: 10 mg

Product Details

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

Chemical Properties

Appearance	Solid yellow
Molecular Weight	226.66
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	389 nm
Emission Wavelength	483 nm

Applications

Monochlorobimane (mBCl) is essentially nonfluorescent until reacted with a thiol group. It readily reacts with several low molecular weight thiols, including cysteine, glutathione, N-acetylcysteine, mercaptopurine, peptides and plasma thiols. The glutathione conjugate of mBCl has absorption/emission maxima ~394/490 nm.