

bBBr [Dibromobimane] *CAS 68654-25-1*

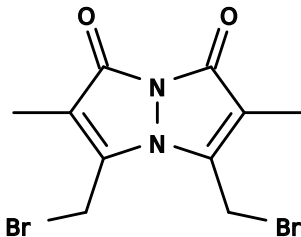
Catalog Number: 634

Unit Size: 5 mg

Product Details

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

Chemical Properties

Appearance	Solid yellow
Molecular Weight	350.01
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	394 nm
Emission Wavelength	490 nm

Applications

bBBr is one of the most popular thiol-reactive fluorescent tags, and is widely used to detect various thiol-containing biomolecules such as glutathione in cells. It is fluorogenic upon reacting with thiol-containing molecules. Dibromobimane is a unique fluorogenic crosslinking reagent for proteins because it is unlikely to fluoresce until both of its alkylating groups have reacted. It has been used to crosslink thiols in myosin, hemoglobin, Escherichia coli lactose permease and mitochondrial ATPase. Dibromobimane is also used to probe for the proximity of dual-cysteine mutagenesis sites in ArsA ATPase and to crosslink thiols in actin, myosin and P-glycoprotein.