

mBBR [Monobromobimane] *CAS 71418-44-5*

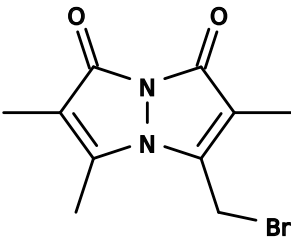
Catalog Number: 633

Unit Size: 25 mg

Product Details

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

Chemical Properties

| | |
|--------------------|--|
| Appearance | Solid light yellow |
| Molecular Weight | 271.11 |
| Soluble In | DMSO |
| Chemical Structure |  |

Spectral Properties

| | |
|-----------------------|--------|
| Excitation Wavelength | 395 nm |
| Emission Wavelength | 490 nm |

Applications

mBBR 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. mBBR is used for the determination of the redox status of low molecular weight and protein thiols in biological systems. MBBR-based in situ derivatization results in maximal recovery of both free, reduced low molecular weight and bromobimane accessible protein thiols. The quantitation of the corresponding adducts of protein thiols is achieved by fluorescence spectroscopy following protein precipitation.