

Product Information

Monobromobimane (mBBr)

Catalog Number	Unit Size
C038	25 mg

Storage upon receipt:

- -20°C
- Protect from light

Product Description

Monobromobimane (mBBr) is essentially nonfluorescent, and freely passively diffuse across the plasma membrane into the cytoplasm, where they form blue-fluorescent adducts with intracellular glutathione and thiol containing proteins. The glutathione conjugate of monobromobimane has absorption/emission maxima ~394/490 nm. mBBr has been reported to be used as effective tracer of gap-junctional communication.

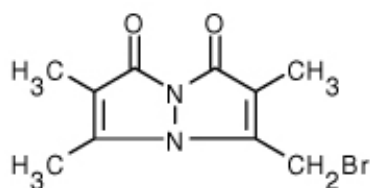
Specifications

Chemical Name: monobromobimane (mBBr)

Molecular Formula: C₁₀H₁₁BrN₂O₂

Molecular Weight: 271.11

CAS Number: 71418-44-5



Application

This reactive probe.