

MM 4-64 Fixable Membrane Stain

Catalog Number: 21488

Unit Size: 5x100 ug

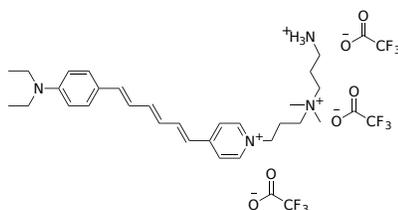
Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	24 months upon receiving

Chemical Properties

Appearance	Solid purple
Molecular Weight	788.75
Soluble In	Water

Chemical Structure



Spectral Properties

Excitation Wavelength	506 nm
Emission Wavelength	725 nm

Applications

MM4-64 Fixable Membrane Stain is a lipophilic probe with low fluorescence in aqueous solution that intensifies upon binding to the plasma membrane's outer leaflet, enabling precise membrane staining. Its rapid, reversible binding is compatible with aldehyde fixation and is effective in identifying active neurons and studying activity-dependent vesicle cycling.