

Yeast Vacuole Membrane Marker MDY-64

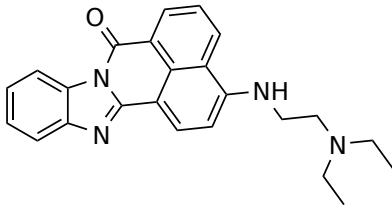
Catalog Number: 24010

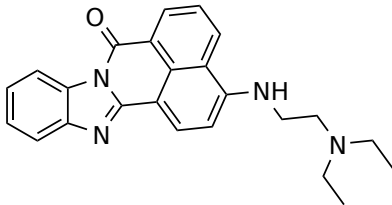
Unit Size: 1 mg

Product Details

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

Chemical Properties

Appearance	Solid orange
Molecular Weight	384.48
Soluble In	DMSO
Chemical Structure	



Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

MDY-64 is a fluorescent dye used as a specific marker for the vacuole membrane in yeast cells, allowing researchers to visualize the vacuole structure and dynamics within a yeast cell under a microscope. It is a stain that specifically binds to the membrane of the yeast vacuole. Its fluorescence can be readily monitored with the FITC filter set. Vacuole morphology can be conveniently monitored by fluorescence microscopy.