

Cal-520® alkyne

Catalog Number: 20614

Unit Size: 100 ug

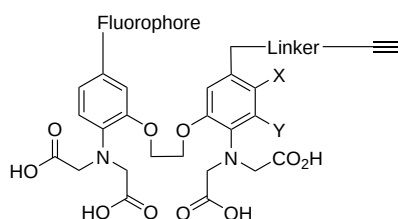
Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	905.77
Soluble In	DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength	492 nm
Emission Wavelength	515 nm

Applications

Cal-520® alkyne is an excellent building block that can be readily used to prepare a calcium-sensitive bioconjugate for monitoring calcium change spatially for a specific target. It can be conjugated to azido-containing biomolecules via the well-known click chemistry. Calcium measurement is critical for numerous biological and neurological investigations. Fluorescent probes showing spectral responses upon binding calcium have enabled researchers to investigate changes in intracellular free calcium concentrations using fluorescence microscopy, flow cytometry, fluorescence spectroscopy, and fluorescence microplate readers. Cal-520® has been proven to be one of the best-known green fluorescent calcium indicators.