

## Cal-520® amine

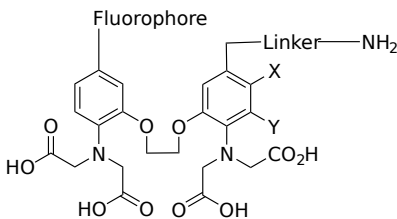
Catalog Number: 20612

Unit Size: 100 ug

### Product Details

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

### Chemical Properties

Appearance	Solid orange
Molecular Weight	909.68
Soluble In	DMSO
Chemical Structure	

### Spectral Properties

Excitation Wavelength	492 nm
Emission Wavelength	515 nm

### Applications

Cal-520® amine 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 a carboxy group via activated esters or EDC coupling 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.