

Metal Fluor™ Zn-520, AM

Catalog Number: 21263

Unit Size: 1 mg

Product Details

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

Chemical Properties

Appearance	Solid off-white
Molecular Weight	721.62
Soluble In	DMSO

Spectral Properties

Excitation Wavelength	493 nm
Emission Wavelength	516 nm

Applications

Metal Fluor™ Zn-520 is designed for detection of higher zinc ion concentrations that are present in synaptic vesicles and released in response to electrical stimulation or excitotoxic agonists (0.05-50 μ M) with minimal interfering calcium sensitivity. Metal Fluor™ Zn-520 has shown great fluorescence enhancement upon binding zinc ion (>250 folds). This cell permeant AM-ester is useful for detecting intracellular zinc ion levels.