

iFluor® 450 acid

Catalog Number: 2905

Unit Size: 5 mg

Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	405.42
Soluble In	DMSO

Spectral Properties

Excitation Wavelength	451 nm
Emission Wavelength	502 nm

Applications

iFluor® 450 is a blue-fluorescent dye that is excited by the 405 nm laser line and emits green fluorescence with a peak at 502 nm (Ex/Em = 451/502 nm). It exhibits stronger absorption and brighter fluorescence than Alexa Fluor® 450, along with excellent photostability, water solubility, and stable performance across a wide pH range (pH 3 to 11). With a moderate Stokes Shift and clean spectral separation, iFluor® 450 is particularly well-suited for multicolor flow cytometry and confocal microscopy. iFluor® 450 acid is the non-reactive form of the dye and serves as a reference standard or fluorescence control in assays utilizing iFluor® 450 conjugates. It is compatible with ICC, HCS, live-cell imaging, and flow cytometry. It can also be chemically modified to generate activated esters or STP derivatives and coupled to hydrazines, hydroxylamines, or amines using water-soluble carbodiimides such as EDAC, allowing efficient labeling of proteins, antibodies, peptides, and amine-modified oligonucleotides.