

iFluor® 350 acid

Catalog Number: 2900

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	N/A
Soluble In	DMSO

Spectral Properties

Excitation Wavelength	345 nm
Emission Wavelength	450 nm

Applications

iFluor® 350 is a blue-fluorescent dye that exhibits spectral properties similar to AMCA and Alexa Fluor® 350. It offers significant advantages over traditional fluorophores, including enhanced brightness, improved water solubility, and greater pH resistance across a wide molar range (pH 3-11). With an excitation peak at 345 nm and maximum emission at 450 nm, iFluor® 350 is ideally suited for use with UV excitation sources (e.g., the 350 nm laser). Its exceptional photostability, coupled with its moderate brightness, makes iFluor® 350 conjugates an excellent choice for direct imaging of high-abundance targets.

iFluor® 350 Key Features

- **High performance:** iFluor® 350 shows brighter fluorescence than Alexa Fluor® 350 and AMCA
- **Ex/Em maxima:** 345/450 nm
- **Quantum yield:** 0.95
- **Extinction coefficient:** 20,000 cm⁻¹M⁻¹
- **Spectrally similar dyes:** AMCA, Marina Blue, Alexa Fluor® 350

iFluor® 350 acid is a non-reactive compound that can be employed as a reference standard in studies utilizing iFluor® 350 conjugates. It is also suitable for use as a control in confocal microscopy, immunocytochemistry (ICC), high-content screening (HCS), flow cytometry, and live cell imaging applications. Furthermore, it can be utilized in the synthesis of activated esters and STP and can be coupled to hydrazines, hydroxylamines, or amines in aqueous solutions using water-soluble carbodiimides (e.g., EDAC). This allows for the conjugation of the dye to amino-containing molecules, such as proteins, antibodies, amine-modified oligonucleotides, and peptides.