

## iFluor® 430 acid

Catalog Number: 2950

Unit Size: 5 mg

### Product Details

---

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

### Chemical Properties

---

Appearance	Solid yellow
Molecular Weight	490.53
Soluble In	DMSO

### Spectral Properties

---

Excitation Wavelength	433 nm
Emission Wavelength	497 nm

### Applications

---

iFluor® 430 is a blue-fluorescent dye that is excited by the 405 nm laser line and emits green fluorescence with a peak at 498 nm. It offers brighter signal and stronger absorption than Alexa Fluor® 430, with improved aqueous solubility and excellent stability across a wide pH range (pH 3 to 11). These properties make iFluor® 430 ideal for demanding multicolor applications like flow cytometry and confocal microscopy, where spectral separation and photostability are critical. iFluor® 430 acid is the non-reactive version of the dye and serves as a reference standard or fluorescence control in assays using iFluor® 430 conjugates. It is suitable for use in ICC, confocal microscopy, flow cytometry, HCS, and live-cell imaging. 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, enabling efficient conjugation to proteins, antibodies, peptides, and amine-modified oligonucleotides.