

## 5-CFDA, AM [5-Carboxyfluorescein Diacetate, Acetoxymethyl Ester]

Catalog Number: 21910

Unit Size: 1 mg

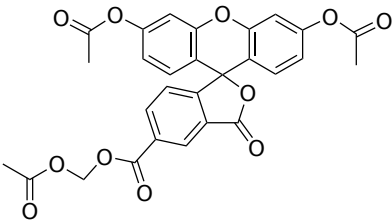
### Product Details

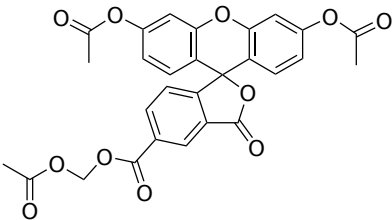
---

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

### Chemical Properties

---

|                    |  |
|--------------------|--|
| Appearance         | Solid white  |
| Molecular Weight   | 532.46   |
| Soluble In         | DMSO   |
| Chemical Structure |  |



### Spectral Properties

---

|                       |     |
|-----------------------|-----|
| Excitation Wavelength | N/A |
| Emission Wavelength   | N/A |

### Applications

---

5-Carboxyfluorescein diacetate, acetoxymethyl ester (CFDA, AM) is a cell-permeant esterase substrate that can serve as a viability probe that measures both enzymatic activity, which is required to activate its fluorescence, and cell-membrane integrity, which is required for intracellular retention of their fluorescent product. Upon hydrolysis by intracellular esterases, CFDA, AM yields the 5-carboxyfluorescein product that gives bright green fluorescence when excited with a blue light or the common 488 nm Argon laser. It is often used to monitor cell viability or label live cells for cell analysis.