

Product Information

CFDA

Catalog Number	Product Name	Unit Size
C031	CFDA	100 mg

Storage upon receipt:

- -20°C
- Protect from light

Product Description

5(6)-CFDA 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, this ester yields carboxyfluorescein. As compared with fluorescein, carboxyfluorescein contains extra negative charges and is therefore better retained in cells.

Specifications:

Chemical name: 5-(and-6)-Carboxyfluorescein Diacetate

Molecular Formula: C₂₅H₁₆O₉

Molecular Weight: 460.40

CAS Number: 124387-19-5