
Product Information

CDCFDA

Cat. No.	Product Name	Unit Size
C035	CDCFDA	100 mg

Storage upon receipt:

- -20°C
- Protect from light

Product Description

5(6)-CDCFDA is a cell membrane permeant probe with low sensitivity to fluorescence to intracellular pH which can be loaded into the cell where it is hydrolysed by cytosolic esterases to carboxy-dichlorofluorescein. It can be used as an acidic pH sensor with a pKa of 4.8.

Specifications:

Chemical name: 5-(and-6)-Carboxy-2',7'-dichlorofluorescein Diacetate

Molecular Formula: C₂₅H₁₄Cl₂O₉

Molecular Weight: 529.28

CAS Number: 127770-45-0