5(6)-CDCFDA (ultra pure)

ROS Detector and pH indicator

5(6)-CDCFDA is an amine—reactive fluorescein diacetate (FDA) derivative used to prepare a variety of FDA conjugates. 5(6)-CDCFDA is a non-fluorescent molecule that diffuse into cells and is hydrolyzed by intracellular non-specific esterases, yielding a fluorescent probe. The probe accumulates exclusively in cells with intact cell membranes. 5(6)-CFCFDA does not stain dead cells. 5(6)-CDCFDA labeling can be used for monitoring cells by flow cytometry or fluorescence microscopy. The probe is also frequently used to monitor reactive oxygen species (ROS) generation in live cells. Wavelength Maxima: Excitation 504nm, Emission 529nm

Citations: 5

View Online »

Ordering Information

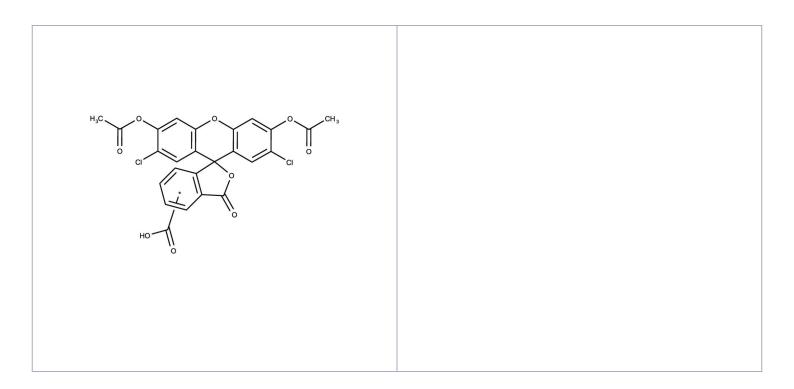
Order Online »

ENZ-52103

100mg

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability Stable for at least one year after receipt when stored as recommended.

Handling Protect from light. Keep cool and dry.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 5(6)-Carboxy-2',7'-dichlorofluorescein diacetate

CAS 127770-45-0

Formula $C_{25}H_{14}Cl_2O_9$

MW 529.3

Purity ≥98% (HPLC)

Solubility Soluble in DMSO.

Technical Info / Product NotesThis product is a member of the CELLESTIAL[®] product

line, reagents and assay kits comprising fluorescent molecular probes that have been extensively benchmarked for live cell analysis applications.

CELLESTIAL® reagents and kits are optimal for use in demanding imaging applications, such as confocal

microscopy, flow cytometry and HCS, where consistency

and reproducibility are required.

Last modified: May 29, 2024

