

# 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 diffuses 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 »](#)

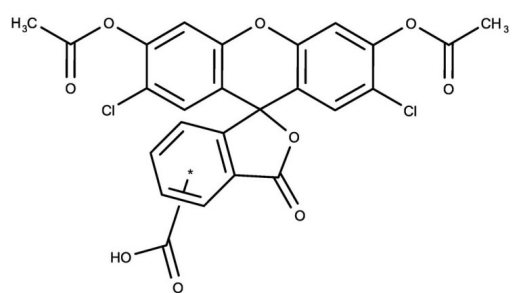
## Ordering Information

[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 Notes	This product is a member of the CELLESTIAL <sup>®</sup> product line, reagents and assay kits comprising fluorescent molecular probes that have been extensively benchmarked for live cell analysis applications. CELLESTIAL <sup>®</sup> 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.
--------------------------------	---



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)