

2',7'-

Dichlorodihydrofluorescein diacetate

Peroxynitrite indicator

Cell permeable, sensitive indicator of peroxynitrite formation. After hydrolysis of the diacetate groups by cytosolic esterases or base-catalyzed cleavage of the diacetate groups, DCDHF is oxidized by peroxynitrite to the highly fluorescent product dichlorofluorescein (DCF). Formation of DCF can be monitored by fluorescence spectroscopy (Ex(max): 502nm, Em(max): 523nm), or by absorbance spectroscopy at 500nm ($\epsilon=59,500\text{M}^{-1}\text{cm}^{-1}$). Neither nitric oxide, superoxide nor hydrogen peroxide alone appear to oxidize DCDHF.

Citations: 16

[View Online »](#)

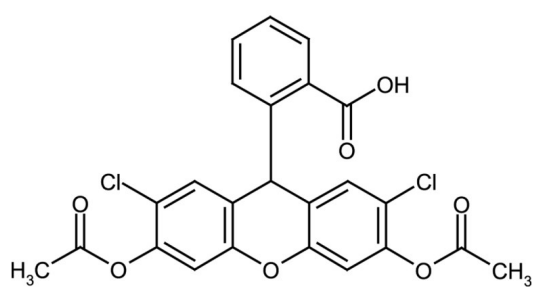
Ordering Information

[Order Online »](#)

ALX-610-022-M050	50mg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solution in organic solvents is stable for at least 6 months if stored at -20°C. For maximum solubility in aqueous buffers, dissolve directly in 0.1M disodium carbonate (7.7mg/ml) and then immediately dilute with PBS, pH 7.2, to the desired concentration.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	DCDHF diacetate, 2',7'-Dichlorofluorescein diacetate
Appearance	White to light orange-brown solid.
CAS	4091-99-0
Formula	$C_{24}H_{16}O_7Cl_2$
MW	487.3
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO (≥25mg/ml), 100% ethanol or dimethyl formamide; sparingly soluble in aqueous buffers.



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-
fr@enzolifesciences.com](mailto:info-fr@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)