

# CFSE

## Amine dye

Non-toxic cell permeable fluorescent dye that reacts with amine compounds. Readily accumulates inside viable cells. Hardly leaks from cells when covalently attached to intracellular proteins. Hydrolyzed CFSE emits fluorescence (Ex(max): 495nm; Em(max): 519nm). Therefore, CFSE is utilized for the long-term observation of cell activities by fluorescent microscopy as well as for viable cell detection.

**Citations:** 12

[View Online »](#)

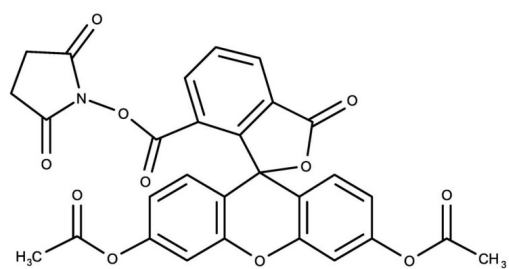
## Ordering Information

[Order Online »](#)

ALX-610-030-M025	25mg
------------------	------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	5(6)-(N-Succinimidylloxycarbonyl)-3',6',0,0'-diacetylfluorescein, 5(6)-Carboxyfluorescein diacetate, N-succinimidyl ester-
Appearance	White or slightly yellow powder.
CAS	150347-59-4
Formula	$C_{29}H_{19}NO_{11}$
MW	557.5
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO, dimethyl formamide or methanol.



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)