

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 »](#)

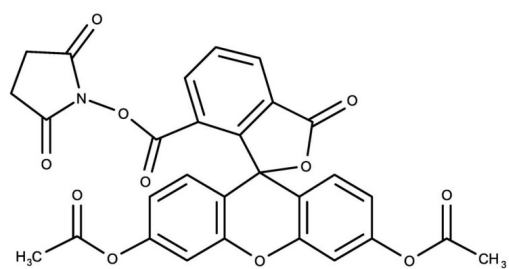
Ordering Information

[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. |

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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