


eBioscience™ CFSE


Catalog Number: 65-0850


Also known as: 5-(and 6)-Carboxyfluorescein diacetate succinimidyl ester, CFDA SE

For Research Use Only. Not for use in diagnostic procedures.

Product Information

Contents: eBioscience™ CFSE
 **Catalog Number:** 65-0850

Formulation: lyophilized
 **Temperature Limitation:** Store at -20°C to -80°C. Protect from light and moisture. It is recommended to use the reconstituted dye within 6 months and to avoid freeze-thawing.

 **Batch Code:** Refer to vial

 **Use By:** Refer to vial

Description

CFSE is widely used for cell tracking and proliferation studies. It has also been used in CTL assays and cell motility studies. CFSE readily crosses intact cell membranes. Once inside the cells, intracellular esterases cleave the acetate groups to yield the fluorescent carboxyfluorescein molecule. The succinimidyl ester group reacts with primary amines, crosslinking the dye to intracellular proteins. Cell division can be measured as successive halving of the fluorescence intensity of CFSE. Cells labeled with CFSE may be fixed and permeabilized for analysis of intracellular targets using standard formaldehyde-containing fixatives and saponin-based permeabilization buffers, such as the Foxp3/Transcription Factor Staining Buffer Set (cat. 00-5523) or the IC Fixation Buffer (cat. 00-8222) and Permeabilization Buffer (10X) (cat. 00-8333).

CFSE has a molecular weight of 557.47. After the acetate groups are cleaved, it has a peak excitation of 494 nm and peak emission of 521 nm. Each vial of CFSE may be reconstituted to a stock concentration of 10 mM with 90 µL of anhydrous DMSO; once reconstituted it should be used within 6 months and protected from light and stored at -20°C with desiccant; avoid freeze-thawing.

Applications Reported

CFSE has been reported for use in flow cytometric analysis and fluorescent microscopy.

Applications Tested

CFSE [5-(and 6)-carboxyfluorescein diacetate succinimidyl ester] has been tested by flow cytometric analysis of stimulated mouse splenocytes. Cells can be labeled with 0.5 to 20 µM CFSE. It is recommended that the concentration used for labeling cells be carefully determined by each investigator for optimal performance in the assay of interest.

References

Budda SA, Girton A, Henderson JG, Zenewicz LA. Transcription Factor HIF-1a Controls Expression of the Cytokine IL-22 in CD4 T Cells. *J Immunol.* 2016 Oct 1;197(7):2646-52. (CFSE, FC, PubMed)

Parish CR, Glidden MH, Quah BJ, Warren HS. Use of the intracellular fluorescent dye CFSE to monitor lymphocyte migration and proliferation. *Curr Protoc Immunol.* 2009 Feb;Chapter 4:Unit4.9.

Ingulli E. Tracing tolerance and immunity in vivo by CFSE-labeling of administered cells. *Methods Mol Biol.* 2007;380:365-76.

Miller MJ, Wei SH, Parker I, Cahalan MD. Two-photon imaging of lymphocyte motility and antigen response in intact lymph node. *Science.* 2002 Jun 7;296(5574):1869-73.

Lyons AB, Parish CR. Determination of lymphocyte division by flow cytometry. *J Immunol Methods.* 1994 May 2;171(1):131-7.

Not for further distribution without written consent.

Copyright © 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

Tel: 888.999.1371 or 858.642.2058 • Fax: 858.642.2046 • thermofisher.com/ebioscience •

info@ebioscience.com

eBioscience™ CFSE

Catalog Number: 65-0850

Also known as: 5-(and 6)-Carboxyfluorescein diacetate succinimidyl ester, CFDA SE

For Research Use Only. Not for use in diagnostic procedures.

Related Products

45-0900 CD90.1 (Thy-1.1) Monoclonal Antibody (HIS51), PerCP-Cyanine5.5, eBioscience™ TDS DISABLED: ABMAINT SKU (HIS51)

48-0081 CD8a Monoclonal Antibody (53-6.7), eFluor 450, eBioscience™ TDS DISABLED: ABMAINT SKU (53-6.7)

65-0840 eBioscience™ Cell Proliferation Dye eFluor™ 670

65-0842 eBioscience™ Cell Proliferation Dye eFluor™ 450

Not for further distribution without written consent.

Copyright © 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

Tel: 888.999.1371 or 858.642.2058 • Fax: 858.642.2046 • thermofisher.com/ebioscience •

info@ebioscience.com

CFSE Cell Labeling Protocol

Protocol: CFSE Cell Labeling

Materials Needed

- PBS
- Flow Cytometry Staining Buffer (cat. 00-4222)

Experimental Procedure

Note: Reconstitute one vial of CFSE to a stock concentration of 10 mM with 90 μ L of anhydrous DMSO. Once reconstituted the dye should be used within 6 months, protected from light, and stored with desiccant at less than or equal to -20°C . Avoid freeze-thawing.

1. Prepare a single-cell suspension of cells to be labeled.
2. Wash cells two times with PBS to remove any serum.
3. Resuspend cells at $5\text{-}10 \times 10^6/\text{mL}$ of PBS (pre-warmed to room temperature).
4. Add CFSE to the desired final concentration (e.g., for a final concentration of $1 \mu\text{M}$, add $0.2 \mu\text{L}$ of a 5 mM stock solution per mL of cells).
5. Mix immediately and incubate for 10 minutes at room temperature in the dark.
6. Stop labeling by adding 4-5 volumes of cold complete media (containing $\geq 10\%$ serum) and incubate on ice for 5 minutes.
7. Wash cells 3 times with complete media.
8. Culture or transfer cells, as desired.

Note: The concentration of CFSE, incubation time, and temperature can be modified to achieve the desired staining intensity. However, very high labeling can lead to compensation issues and may also interfere with cellular functions. Thus, it is highly recommended that each investigator determine the optimal concentration for the assay of interest.

Documentation and support

Customer and technical support

Visit thermofisher.com/support for the latest in services and support, including:

- Worldwide contact telephone numbers
- Product support, including:
 - Product FAQs
 - Software, patches, and updates
- Order and web support
- Product documentation, including:
 - User guides, manuals, and protocols
 - Certificates of Analysis
 - Safety Data Sheets (SDSs; also known as MSDSs)

Note: For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at thermofisher.com/support.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

Corporate entity: Life Technologies | Carlsbad, CA 92008 USA | Toll Free in USA 1.800.955.6288

©2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. All other trademarks are properties of their respective owners.

For support visit thermofisher.com/support or email techsupport@lifetech.com

thermofisher.com

23 January 2017

ThermoFisher
SCIENTIFIC