
CellVue™ NIR815 Cell Labeling Kit

Catalog Number: 88-0874

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

Product Information

Contents: CellVue™ NIR815 Cell Labeling Kit



Catalog Number: 88-0874



Formulation: 1 mM dye stock in ethanol
Temperature Limitation: Store at 2-8°C Light sensitive material.



Batch Code: Refer to vial



Use By: Refer to vial

Description

CellVue® dyes are lipophilic dyes that can be used to label the cell membrane for the purpose of identifying and tracking labeled cells. Cell labeling is rapid and stable and can be combined with fluorescently labeled antibodies and other markers of cellular function for flow cytometric analysis and fluorescent microscopy. Mini CellVue® Kits are supplied with one vial of dye stock (1 mM in ethanol) and one vial of labeling vehicle (Diluent C).

CellVue® NIR815 is a near-infrared fluorescent cell labeling reagent. It is optimally excited at 786 nm and has a peak emission of 814 nm. CellVue® NIR815 has been reported to be useful for short-term in vivo cell tracking studies using a Pearl™ Imager or Odyssey Infrared Imaging System (Li-COR Biosciences, Inc., Lincoln, NE).

Applications Reported

CellVue® NIR815 Cell Labeling Kit has been reported for use in microscopy, immunocytochemistry, and cell labeling.

References

Roy EJ, Sivaguru M, Fried G, Gray BD, Kranz DM. Imaging membrane intercalating near infrared dyes to track multiple cell populations. *J Immunol Methods*. 2009 Aug 31;348(1-2):18-29.

Thomas DL, Kranz DM, Roy EJ. Experimental manipulations of afferent immune responses influence efferent immune responses to brain tumors. *Cancer Immunol Immunother*. 2008 Sep;57(9):1323-33.

Related Products

00-4501 CellVue Diluent C (For CellVue™ Cell Labeling Kits)

Legal

CellVue® is a registered trademark of PTI Research, Inc. Patent <http://www.mtarget.com/mm5/pdfs/patents.pdf>

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

CellVue™ Labeling Kit Protocol

Protocols: CellVue™ Kit General Cell Labeling

Experimental Procedure

Note: All procedures should be carried out at room temperature. No azide or other metabolic inhibitors should be present during labeling.

1. Wash cells one time with serum-free media to remove serum proteins and lipids that may interfere with cell labeling.
2. Carefully discard the supernatant, leaving no more than 25 μ L of residual media on the pellet.
3. Gently resuspend cells at 2×10^7 cells/mL with Diluent C. Do not vortex, pipette several times to be sure cells are in a single-cell suspension.
4. Immediately before labeling, prepare a 2X working solution of dye in Diluent C. For a final labeling concentration of 2 μ M, prepare a 4 μ M solution by adding 4 μ L of the 1 mM dye stock to 1 mL of Diluent C. To ensure uniform labeling, never add dye stock directly to cell suspension.
5. Rapidly add 1 mL of cells to 1 mL of 2X working dye solution and immediately mix the sample by pipetting to ensure uniform labeling. The final cell concentration should be approximately 1×10^7 cells/mL and the final dye concentration should be 2 μ M.
6. Incubate the cells for 2–5 minutes, with periodic mixing. Longer incubations will result in brighter labeling, but may begin to adversely affect cell viability.
7. Stop labeling by adding an equal volume of serum and let sit for 1 minute.
8. Centrifuge cells and discard supernatant.
9. Wash cells three times with complete media (do not use Diluent C). To reduce any effect of residual dye bound to the tube, cells may be transferred to a fresh tube.
10. Count, culture or transfer cells as desired.

Note: This protocol has been tested for in vitro or ex vivo labeling of cells. Because labeling occurs as a function of the dye partitioning into the cell membrane, concentrations of dye that are too high or labeling times that are too long will result in a loss of membrane integrity and poor cell recovery. Labeling with CellVue™ dyes should be done prior to staining with monoclonal antibodies. Labeling conditions and concentrations should be carefully determined by each investigator to ensure optimal performance in the assay of interest.

Additional Diluent C can be purchased separately. Refer to Cat. No. 00-4501.

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