

## Image-iT™ LIVE Lysosomal and Nuclear Labeling Kit (I34202)

### Quick Facts

#### Storage upon receipt:

- $\leq -20^{\circ}\text{C}$
- Protect from light
- Desiccate
- Avoid freeze-thaw cycles

#### Ex/Em:

- 577/590 nm (LysoTracker® Red DND-99 dye)
- 350/461 nm (Hoechst 33342 dye)

**Number of assays:** 500

### Introduction

The Image-iT™ LIVE Lysosomal and Nuclear Labeling Kit provides two stains — red-fluorescent LysoTracker® Red DND-99 dye and blue-fluorescent Hoechst 33342 dye — for highly selective staining of lysosomes and the nucleus, respectively, in live green-fluorescent protein (GFP)–transfected cells. When used according to the sample protocol, cell-permeant LysoTracker Red DND-99 dye provides highly selective lysosomal staining with minimal background. A significant amount of specific staining is retained after formaldehyde fixation, although some cytoplasmic background staining may be seen. Hoechst 33342 dye, a cell-permeant nucleic acid stain that is selective for DNA and spectrally similar to DAPI, is UV excitable and emits blue fluorescence when bound to DNA. This dye should not interfere with GFP fluorescence and is retained after fixation and permeabilization.

It is not recommended that the dyes be combined into one staining solution. The dyes should be used in separate labeling steps, with Hoechst 33342 staining first.

### Materials

#### Kit Contents

- **LysoTracker Red DND-99 dye** (Component A), 50  $\mu\text{L}$  at 1.0 mM in dimethylsulfoxide (DMSO)
- **Hoechst 33342 dye** (Component B), 3 vials, each containing 400  $\mu\text{L}$  at 1.0 mM in water

#### Storage and Handling of Stock Solutions

Upon receipt, the kit should be stored upright, desiccated, and protected from light at  $\leq -20^{\circ}\text{C}$ . Avoid freeze-thaw cycles. Vials should be allowed to warm to room temperature before opening. When stored properly, components should be stable for at least 6 months.

**Note:** Handle stock solutions containing DMSO with care. DMSO (found in the LysoTracker Red DND-99 dye solution) is readily absorbed through the skin. Vials should be tightly closed and stored with desiccant.

#### Spectral Characteristics

LysoTracker Red DND-99 dye has excitation/emission maxima of approximately 577/590 nm; Hoechst 33342 dye has excitation/emission maxima of approximately 350/461 nm. Cells labeled with LysoTracker Red DND-99 dye and Hoechst 33342 dye can be imaged using standard filter sets.

#### Materials Recommended but Not Provided

Hank's balanced salt solution (HBSS, available from Gibco (14025-092)).

### Experimental Protocol

#### Labeling Live Eukaryotic Cells

This is a general procedure for labeling live, cultured cells that are adhering to coverslips. The protocol has been optimized using HBSS for HeLa cells transfected with GFP. LysoTracker Red DND-99 and Hoechst 33342 dyes must be used in separate staining steps. Recommended times and concentrations may vary in different model systems and may require optimization.

**1.1 Prepare labeling solutions.** Dilute the 1.0 mM solution of Hoechst 33342 dye (Component B) into HBSS or the appropriate medium for the cell line being used; a recommended concentration for Hoechst 33342 dye is 2.0  $\mu\text{g}/\text{mL}$ . Dilute the 1.0 mM of LysoTracker Red DND-99 dye into a separate volume of HBSS; a recommended concentration for LysoTracker Red DND-99 dye is 100 nM. Note: the two dyes cannot be combined in one staining solution.

**1.2 Stain cells with Hoechst 33342 dye.** Apply a sufficient amount of 2.0  $\mu\text{g}/\text{mL}$  Hoechst 33342 labeling solution (prepared above) to cover cells adhering to coverslip(s). Incubate for 5 minutes at 37°C.

**1.3 Wash cells.** When labeling is complete, remove the labeling solution and wash cells twice in room-temperature HBSS or suitable buffer.

**1.4 Stain cells with LysoTracker Red.** Apply a sufficient amount of 100 nM LysoTracker Red DND-99 labeling solution (prepared above) to cover cells adhering to coverslip(s). Incubate for 1 minute at room temperature. Longer label times may result in unwanted cytoplasmic background, perhaps due to an alkalinizing effect on the lysosomes.

**1.5 Wash and view.** Quickly remove labeling solution and gently wash the cells. Unless the cells will be fixed, samples are ready to be mounted in HBSS or suitable buffer for imaging.

**1.6 (Optional) Fix cells.** Wash the stained cells gently with HBSS or suitable buffer and fix with 4% formaldehyde for 15 minutes at 37°C. Wash in HBSS or suitable buffer and add any additional counterstains. Fixation may cause some cytoplasmic background.

---

**Product List** *Current prices may be obtained from our Web site or from our Customer Service Department.*

Cat #	Product Name	Unit Size
I34202	Image-iT™ LIVE Lysosomal and Nuclear Labeling Kit *counterstains for GFP-expressing cells* .....	1 kit

---

**Contact Information**

Further information on Molecular Probes products, including product bibliographies, is available from your local distributor or directly from Molecular Probes. Customers in Europe, Africa and the Middle East should contact our office in Paisley, United Kingdom. All others should contact our Technical Service Department in Eugene, Oregon.

Please visit our website — [probes.invitrogen.com](http://probes.invitrogen.com) — for the most up-to-date information.

**Molecular Probes, Inc.**

29851 Willow Creek Road, Eugene, OR 97402  
Phone: (541) 465-8300 • Fax: (541) 335-0504

**Customer Service:** 6:00 am to 4:30 pm (Pacific Time)

Phone: (541) 335-0338 • Fax: (541) 335-0305 • [probesorder@invitrogen.com](mailto:probesorder@invitrogen.com)

**Toll-Free Ordering for USA:**

Order Phone: (800) 438-2209 • Order Fax: (800) 438-0228

**Technical Service:** 8:00 am to 4:00 pm (Pacific Time)

Phone: (541) 335-0353 • Toll-Free (800) 438-2209  
Fax: (541) 335-0238 • [probestech@invitrogen.com](mailto:probestech@invitrogen.com)

**Invitrogen European Headquarters**

Invitrogen, Ltd.  
3 Fountain Drive  
Inchinnan Business Park  
Paisley PA4 9RF, UK  
Phone: +44 (0) 141 814 6100 • Fax: +44 (0) 141 814 6260  
Email: [euroinfo@invitrogen.com](mailto:euroinfo@invitrogen.com)  
Technical Services: [eurotech@invitrogen.com](mailto:eurotech@invitrogen.com)

Molecular Probes products are high-quality reagents and materials intended for research purposes only. These products must be used by, or directly under the supervision of, a technically qualified individual experienced in handling potentially hazardous chemicals. Please read the Material Safety Data Sheet provided for each product; other regulatory considerations may apply.

**Limited Use Label License**

For research use only. Not intended for any animal or human therapeutic or diagnostic use. The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) to not transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. Invitrogen Corporation will not assert a claim against the buyer of infringement of the above patents based upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Invitrogen is willing to accept return of the product with a full refund. For information on purchasing a license to this product for purposes other than research, contact Molecular Probes, Inc., Business Development, 29851 Willow Creek Road, Eugene, OR 97402. Tel: (541) 465-8300. Fax: (541) 335-0504.

Several Molecular Probes products and product applications are covered by U.S. and foreign patents and patents pending. All names containing the designation ® are registered with the U.S. Patent and Trademark Office.

Copyright 2005, Molecular Probes, Inc. All rights reserved. This information is subject to change without notice.