

# LysoTracker® and LysoSensor™ Probes

Table 1 Contents and storage

Material	Amount	Concentration	Storage	Stability	
LysoTracker <sup>®</sup> and LysoSensor <sup>™</sup> dyes	20 vials, each containing 50 µL	1 mM stock	<ul><li> ≤-20°C</li><li>Desiccate</li><li>Protect from light</li></ul>	When stored as directed, products	
LysoTracker <sup>®</sup> Deep Red	5 vials, each containing 50 μL	solution in anhydrous DMSO	<ul> <li>Avoid freeze-thaw cycles</li> <li>Do not store in a frost-free freezer</li> <li>Store in single-use aliquots, if possible</li> </ul>	are stable for at least 6 months*	
LysoSensor™ Yellow/Blue dextran	5 mg, lyophilized solid	NA	• ≤-20°C • Desiccate • Protect from light	When stored as directed, product is stable at least 1 year	

<sup>\*</sup> If refreezing after use, seal the vial tightly.

Approximate fluorescence excitation and emission, in nm: See Table 2, page 2.

## Introduction

LysoTracker® Probes

Weakly basic amines selectively accumulate in cellular compartments with low internal pH and can be used to investigate the biosynthesis and pathogenesis of lysosomes.<sup>1,2</sup> The LysoTracker® probes are fluorescent acidotropic probes for labeling and tracking acidic organelles in live cells.<sup>4,5</sup> These probes have several important features, including high selectivity for acidic organelles and effective labeling of live cells at nanomolar concentrations. Furthermore, the LysoTracker® probes are available in several fluorescent colors (Table 2, page 2), making them especially suitable for multicolor applications.

The LysoTracker® probes, which consist of a fluorophore linked to a weak base that is only partially protonated at neutral pH, are freely permeant to cell membranes and typically concentrate in spherical organelles. Their mechanism of retention has not been firmly established but is likely to involve protonation and retention in the membranes of the organelles, although staining is generally not reversed by subsequent treatment of the cells with weakly basic cell-permeant compounds. Note that in LysoTracker® dye–stained cells, the lysosomal fluorescence may constitute only a small portion of total cellular fluorescence, making it difficult to quantitate the number of lysosomes by flow cytometry or fluorometry.

MAN0001924 | MP 07525 Revision: 2.0

**Table 2** Summary of our LysoTracker<sup>®</sup> and LysoSensor<sup>™</sup> probes.

Cat. no.	Probe	Abs * (nm)	Em * (nm)	pK <sub>a</sub>
L7525	LysoTracker® Blue DND-22	373	422	NA
L12490	LysoTracker® Blue-White DPX	380	†	NA
L7526	LysoTracker® Green DND-26	504	511	NA
L12491	LysoTracker® Yellow-HCK-123	465	535	NA
L7528	LysoTracker® Red DND-99	577	590	NA
L12492	LysoTracker® Deep Red	647	668	NA
L7533	LysoSensor™ Blue DND-167	373	425	5.1
L7535	LysoSensor <sup>™</sup> Green DND-189	443	505	5.2
L7534	LysoSensor <sup>™</sup> Green DND-153	442	505	7.5
L7545	LysoSensor™ Yellow/Blue DND-160 (PDMP0)	329, 384 ‡	440, 540 ‡	4.2
L22460	LysoSensor™ Yellow/Blue dextran	335, 381 ‡	452, 521 ‡	3.9

<sup>\*</sup> Absorption (Abs) and fluorescence emission (Em) maxima, determined in aqueous buffer or methanol; values may vary somewhat in cellular environments. † Emission is extremely sensitive to environment; stained lysosomes appear blue-white, although the emission maximum in methanol is 576 nm. ‡ Dual-absorption and dual-emission maxima, sensitive to pH (see Figure 1, page 3).

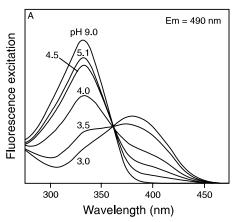
LysoSensor<sup>™</sup> pH Indicators

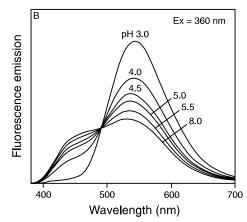
For researchers studying the dynamic aspects of lysosome biogenesis and function in live cells, we offer LysoSensor probes—fluorescent pH indicators that partition into acidic organelles. The LysoSensor dyes are acidotropic probes that appear to accumulate in acidic organelles as the result of protonation. This protonation also relieves the fluorescence quenching of the dye by its weak base side chain, resulting in an increase in fluorescence intensity. Thus, the LysoSensor reagents exhibit a pH-dependent increase in fluorescence intensity upon acidification, in contrast to the LysoTracker probes, which exhibit fluorescence that is largely independent of pH.

Molecular Probes offers five LysoSensor<sup>™</sup> reagents that differ in color and pK<sub>a</sub> (Table 2). Because these probes may localize in the membranes of organelles, it is probable that the actual pK<sub>a</sub> values in cellular environments will differ from the values listed in Table 2 and that only qualitative and semiquantitative comparisons of organelle pH will be possible. The blue and green fluorescent LysoSensor™ probes are available with optimal pH sensitivity in either the acidic or neutral range (p $K_a \sim 5.2$  or  $\sim 7.5$ ). Because of their low pK<sub>a</sub> values, LysoSensor<sup>™</sup> Blue DND-167 and LysoSensor<sup>™</sup> Green DND-189 are almost nonfluorescent except when inside acidic compartments, whereas LysoSensor™ Green DND-153 is brightly fluorescent at neutral pH. LysoSensor™ Yellow/Blue DND-160 (PDMPO) is unique in that it exhibits both dual-excitation and dual-emission spectral peaks that are pH-dependent (Figure 1). Nevertheless, this LysoSensor™ only exhibits the pH-dependent dual-emission spectra in living cells. In acidic organelles LysoSensor<sup>™</sup> Yellow/Blue DND-160 (PDMPO) has predominantly yellow fluorescence, and in less acidic organelles it has blue fluorescence. Dual-emission measurements may permit ratio imaging of the pH in acidic organelles such as lysosomes or the acrosomes of spermatozoa. LysoSensor<sup>™</sup> Yellow/Blue dextran allows loading of the cells by endocytosis. This conjugate should prove useful for studying the endocytic pathway. The pK<sub>a</sub> is somewhat lower than the p $K_a$  of the free LysoSensor<sup>m</sup> Yellow/Blue dye.

These probes can be used singly (or potentially in combination) to investigate the acidification of lysosomes and alterations of lysosomal function or trafficking that occur in cells. For example, lysosomes in some tumor cells have a lower pH than normal lysosomes, while other tumor cells contain lysosomes with higher pH. In addition, cystic fibrosis and other diseases result in defects in the acidification of some intracellular organelles, and the LysoSensor™ probes may prove useful in studying these aberrations. As in LysoTracker® dye—stained cells, the lysosomal fluorescence in LysoSensor™ dye—stained cells may constitute only a small portion of total cellular fluorescence, making it difficult to quantitate the number of lysosomes or their pH by flow cytometry or fluorometry.

Figure 1 The pH-dependent spectral response of LysoSensor™ Yellow/Blue DND-160 (PDMPO, Cat. no. L7545). (A) fluorescence excitation spectra and (B) fluorescence emission spectra.





## Guidelines for Use

Before opening, allow the vial to warm to room temperature and then briefly centrifuge the vial in a microcentrifuge to deposit the DMSO solution at the bottom of the vial.

The concentration of probe for optimal staining will vary depending on the application. Here we suggest some initial conditions to use as a guideline. The staining conditions may need to be modified depending upon the particular cell type and the permeability of the cells or tissues to the probe, among other factors.

LysoTracker® and LysoSensor™ Dyes

1.1 Dilute the 1 mM probe stock solution to the final working concentration in the growth medium or buffer of choice. For the LysoTracker® probes, we recommend working concentrations of 50–75 nM and for the LysoSensor $^{\text{TM}}$  probes at least 1  $\mu$ M. To reduce potential artifacts from overloading, the concentration of dye should be kept as low as possible.

**Note:** If the cells are incubated in dye-free medium after staining, we often observe a decrease in fluorescent signal and cell blebbing

1.2 For adherent cells, grow cells on coverslips inside a Petri dish filled with the appropriate culture medium. When cells have reached the desired confluence, remove the medium from the dish and add the prewarmed (37°C) probe-containing medium. Incubate the cells for 30 minutes to 2 hours under growth conditions appropriate for the particular cell type. Then replace the loading solution with fresh medium and observe the cells using a fluorescence microscope fitted with the correct filter set (see Table 2, page 2). If the cells do not appear to be sufficiently stained, we recommend either increasing the labeling concentration or increasing the time allowed for the dye to accumulate in the lysosomes.

**Note:** Kinetic studies on the internalization of the LysoTracker® Green DND-26 and LysoSensor™ Yellow/Blue DND-160 (PDMPO) probes indicate that the rates of uptake of these dyes into living cells can occur within seconds. Unfortunately, these lysosomal probes can exhibit an "alkalizing effect" on the lysosomes, such that longer incubation with these probes can induce an increase in lysosomal pH. We suggest that these probes are useful pH indicators only when they are incubated with cells for 1–5 minutes at 37°C.

1.3 For suspension cells, centrifuge to obtain a cell pellet and aspirate the supernatant. Resuspend the cells gently in prewarmed (37°C) probe-containing medium. Incubate the cells for 30 minutes to 2 hours under growth conditions appropriate for the particular cell type (see note above regarding internalization rate of these probes). Re-pellet the cells by centrifugation and resuspend in fresh prewarmed medium. Observe the cells using a fluorescence microscope fitted with the correct filter set (Table 2, page 2). If the cells do not appear to be sufficiently stained, we recommend either increasing the labeling concentration or increasing the time allowed for the dye to accumulate in the lysosomes.

Alternatively, suspension cells may be attached to coverslips that have been treated with BD Cell-Tak<sup>®</sup> (BD Biosciences) and stained as if they were adherent cells (see step 1.2).

## LysoSensor™ Yellow/Blue Dextran

- **2.1** To prepare a stock solution, reconstitute the lyophilized dextran to 50 mg/mL in phosphate-buffered saline, pH 7.4. Store the stock solution ≤–20°C, protected from light.
- **2.2** Dilute the stock solution to a final working concentration in the growth medium or buffer of choice. We recommend a working concentration of 1–5 mg/mL.
- 2.3 For adherent cells, grow cells on coverslips inside a Petri dish filled with the appropriate culture medium. When cells have reached the desired confluence, remove the medium from the dish and add the prewarmed (37°C) dextran working solution. Incubate the cells for 1–24 hours under growth conditions appropriate for the particular cell type and experiment. Replace the loading solution with fresh medium and observe the cells using a fluorescence microscopy fitted with the correct filter set (Table 2, page 2).
- 2.4 For suspension cells, centrifuge to obtain a cell pellet and aspirate the supernatant. Resuspend the cells gently in pre-warmed (37°C) dextran-containing medium. Incubate the cells for 1–24 hours under growth conditions appropriate for the particular cell type. Re-pellet the cells by centrifugation and resuspend in fresh prewarmed medium. Observe the cells using a fluorescence microscope fitted with the correct filter set (Table 2, page 2).

### References

1. Cell 52, 329 (1988); 2. Lysosomes in Biology and Pathology, J.T. Dingle et al., Eds., North-Holland Publications Co. (1969); 3. J Cell Biol 106, 539 (1988); 4. Cytometry suppl 7, 77 abstract #426B (1994); 5. Mol Biol Cell 5, 113a abstract #653 (1994); 6. J Cell Biol 126, 877 (1994); 7. J Cell Biol 128, 901 (1994); 8. Molecular Aspects of Anticancer Drug Action, S. Neidle and M.J. Waring, Eds., Macmillian (1983) pp. 233–282; 9. J Biol Chem 265, 4775 (1990); 10. Nature 352, 70 (1991).

## Product List Current prices may be obtained from our website or from our Customer Service Department.

Cat. no.	Product Name	Unit Size
A1301	Acridine orange	1 g
A3568	Acridine orange	10 mL
D1552	N-(3-((2,4-dinitrophenyl)amino)propyl)-N-(3-aminopropyl)methylamine, dihydrochloride	100 mg
L7533	LysoSensor <sup>™</sup> Blue DND-167	_
L7534	LysoSensor <sup>™</sup> Green DND-153	20 × 50 μL
L7535	LysoSensor <sup>™</sup> Green DND-189	
L22460	LysoSensor™ Yellow/Blue dextran, 10,000 MW, anionic, fixable	5 mg
L7545	LysoSensor <sup>™</sup> Yellow/Blue DND-160 (PDMP0) *1 mM solution in DMS0*	20 × 50 μL
L7525	LysoTracker® Blue DND-22	-
L12490	LysoTracker <sup>®</sup> Blue-White DPX	20 × 50 μL
L7526	LysoTracker <sup>®</sup> Green DND-26	20 × 50 μL
L7528	LysoTracker <sup>®</sup> Red DND-99	20 × 50 μL
L12491	LysoTracker® Yellow HCK-123	20 × 50 μL
L12492	LysoTracker® Deep Red	5 × 50 μL

## **Purchaser Notification**

#### Corporate Headquarters

5791 Van Allen Way Carlsbad, CA 92008

USA

Phone: +1 760 603 7200 Fax: +1 760 602 6500

Email: techsupport@lifetech.com

### European Headquarters

Inchinnan Business Park 3 Fountain Drive Paisley PA4 9RF UK Phone: +44 141 814 6100

Phone: +44 141 814 6100
Toll-Free Phone: 0800 269 210
Toll-Free Tech 0800 838 380
Fax: +44 141 814 6260
Tech Fax: +44 141 814 6117

Tech Fax: +44 141 814 6117 Email: euroinfo@invitrogen.com Email Tech: eurotech@invitrogen.com

#### Japanese Headquarters

LOOP-X Bldg. 6F 3-9-15, Kaigan Minato-ku, Tokyo 108-0022 Japan Phone: +81 3 5730 6509

Fax: +81 3 5730 6519 Email: jpinfo@invitrogen.com

Additional international offices are listed at www.lifetechnologies.com

These high-quality reagents and materials must be used by, or directly under the supervision of, a technically qualified individual experienced in handling potentially hazardous chemicals. Read the Safety Data Sheet provided for each product; other regulatory considerations may apply.

#### **Obtaining Support**

For the latest services and support information for all locations, go to www.lifetechnologies.com.

At the website, you can:

- Access worldwide telephone and fax numbers to contact Technical Support and Sales facilities
- Search through frequently asked questions (FAQs)
- Submit a question directly to Technical Support (techsupport@lifetech.com)
- Search for user documents, SDSs, vector maps and sequences, application notes, formulations, handbooks, certificates of
  analysis, citations, and other product support documents
- · Obtain information about customer training
- · Download software updates and patches

#### SDS

Safety Data Sheets (SDSs) are available at www.lifetechnologies.com/sds.

#### Certificate of Analysis

The Certificate of Analysis provides detailed quality control and product qualification information for each product. Certificates of Analysis are available on our website. Go to <a href="https://www.lifetechnologies.com/support">www.lifetechnologies.com/support</a> and search for the Certificate of Analysis by product lot number, which is printed on the product packaging (tube, pouch, or box).

#### Disclaimer

LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

#### 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 <a href="https://www.lifetechnologies.com/termsandconditions">www.lifetechnologies.com/termsandconditions</a>. If you have any questions, please contact Life Technologies at <a href="https://www.lifetechnologies.com/support">www.lifetechnologies.com/support</a>.

#### Limited Use Label License: Research Use Only

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial services of any kind, including, without limitation, reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact **outlicensing@lifetech.com** or Out Licensing, Life Technologies Corporation, 5791 Van Allen Way, Carlsbad, California 92008.

The trademarks mentioned herein are the property of Life Technologies Corporation and/or its affiliate(s) or their respective owners.

©2013 Life Technologies Corporation. All rights reserved.

