

ER-Tracker™ Dyes for Live-Cell Endoplasmic Reticulum Labeling

ER-Tracker™ Blue-White DPX
ER-Tracker™ Green (glibenclamide BODIPY® FL)
ER-Tracker™ Red (glibenclamide BODIPY® TR)

Quick Facts

Storage upon receipt:

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

Ex/Em:

- 374/430–640 nm (ER-Tracker™ Blue-White DPX dye)
- 504/511 nm (ER-Tracker™ Green dye)
- 587/615 nm (ER-Tracker™ Red dye)

Introduction

ER-Tracker™ dyes are cell-permeant, live-cell stains that are highly selective for the endoplasmic reticulum (ER). These dyes rarely stain mitochondria, unlike the conventional ER stain DiOC₆(3), and staining at low concentrations does not appear to be toxic to cells. When cells are stained using the optimized protocol provided, staining patterns are partially retained after treatment with formaldehyde.

ER-Tracker™ Blue-White DPX (E12353) is a member of our Dapoxyl™ (DPX) dye family. DPX dyes have long emission wavelengths, high extinction coefficients, high quantum yields, and large Stokes shifts,¹ and are environment-sensitive probes— with increasing solvent polarity, the fluorescence maxima shift to longer wavelengths and the quantum yield decreases.

ER-Tracker™ Green and ER-Tracker™ Red dyes are the drug conjugates glibenclamide BODIPY® FL and glibenclamide BODIPY® TR, respectively. Glibenclamide (glyburide), a drug taken daily by millions of diabetic patients to correct hyperglycemia, has been used to investigate pancreatic β -cell activity and insulin secretion^{2,3} and to study myocardial cell function and

heart arrhythmia.^{4,5} Glibenclamide binds to the sulphonylurea receptors of ATP-sensitive K⁺ channels,⁶ which are prominent on ER; the pharmacological activity of glibenclamide could potentially affect ER function. Variable expression of sulphonylurea receptors in some specialized cell types may result in non-ER labeling.

Materials

Contents

ER-Tracker™ Blue-White DPX dye (MW = 580.53) is packaged as a set of 20 separate vials, each containing 50 μL of a 1 mM stock solution in high-quality, anhydrous dimethylsulfoxide (DMSO). ER-Tracker™ Green dye (MW = 783.1) and ER-Tracker™ Red dye (MW = 915.23) are each provided as 100 μg of lyophilized material.

Storage and Handling

Upon receipt, these products should be stored with desiccant at $\leq -20^{\circ}\text{C}$. AVOID REPEATED FREEZING AND THAWING; AVOID STORING THESE PRODUCTS IN A FROST-FREE FREEZER. When stored properly, these reagents are stable for at least six months.

Note: DMSO stock solutions should be handled with particular caution as DMSO is known to facilitate the entry of organic molecules into tissues. Dispose of these reagents in compliance with all pertaining local regulations.

Materials Required but not Provided

High-quality, anhydrous DMSO (for ER-Tracker™ Green and ER-Tracker™ Red dyes)

Spectral Characteristics

Spectral properties of the ER-Tracker™ dyes are listed in Table 1.

Table 1. Spectral characteristics of ER-Tracker™ dyes.

Dye	Ex/Em *	Filter
ER-Tracker™ Blue-White DPX dye	374/430–640 †	DAPI or UV longpass
ER-Tracker™ Green dye	504/511	FITC
ER-Tracker™ Red dye	587/615	TRITC

* Excitation (Ex) and emission (Em) maxima, in nm, measured in methanol.
† Dapoxyl™ dyes have highly environment-sensitive fluorescence, resulting in variable peak emission values.

Experimental Protocol

This protocol was optimized using bovine pulmonary artery endothelial cells and has been confirmed in other common cell lines. Recommendations for experimental protocols should be used as a starting point, and optimal labeling conditions for each cell type should be determined empirically.

Reagent Preparation

ER-Tracker™ Blue-White DPX is supplied as aliquots of a 1 mM stock solution in DMSO. Allow each vial to warm to room temperature before use, then briefly centrifuge to deposit the DMSO solution at the bottom of the vial.

ER-Tracker™ Green and ER-Tracker™ Red dyes are supplied as 100 µg of lyophilized material. Prepare a 1 mM stock solution of the appropriate dye: for ER-Tracker™ Green dye, dissolve the contents of the vial in 128 µL of DMSO; for ER-Tracker™ Red dye, dissolve the contents of the vial in 110 µL of DMSO. It is recommended that the 1 mM solution then be separated into aliquots and stored frozen with desiccant.

Cell Preparation and Staining

1.1 Prepare staining solution. Dilute the 1 mM stock solution to the final working concentration. We recommend working

concentrations of 100 nM–1 µM for ER-Tracker™ Blue-White DPX and ~1 µM for ER-Tracker™ Green and ER-Tracker™ Red dyes. To minimize potential labeling artifacts, use the lowest dye concentrations possible. Best results are obtained when staining is performed in Hank's Balanced Salt Solution with calcium and magnesium (HBSS/Ca/Mg, Gibco cat. #14025-092) at 37°C/5% CO₂.

1.2 Stain the cells. For adherent cells, remove the medium from the culture dish, rinse with HBSS, and add prewarmed staining solution. Incubate the cells for approximately 15–30 minutes at 37°C. Replace the staining solution with fresh probe-free medium and view the cells using a fluorescence microscope. If the stained cells are to be fixed, refer to the fixation steps below for the appropriate dye.

Fixation and Permeabilization for ER-Tracker™ Blue-White DPX

2.1 Fix and permeabilize cells. ER-Tracker™ Blue-White DPX signal is only partially retained after formaldehyde fixation. Fix stained cells with 4% formaldehyde for 10–20 minutes at 37°C. If additional staining will be performed, cells can be permeabilized with 0.2% Triton® X-100 for 10 minutes.

2.2 Wash and view cells. After cells are fixed, perform two 5-minute washes in any suitable buffer and view.

Fixation for ER-Tracker™ Green and ER-Tracker™ Red Dyes

3.1 Fix cells. If stained cells are to be fixed, fixation is recommended using 4% formaldehyde for 2 minutes at 37°C.

3.2 Wash and view cells. After fixation, perform two 5-minute washes in any suitable buffer prior to mounting, viewing, or further staining. Permeabilization is not recommended; signal is not retained after permeabilization with Triton® X-100.

References

1. Photochem Photobiol 66, 424 (1997);
2. Diabetes 52, 767 (2003);
3. Biochem Pharmacol 67, 1437 (2004);
4. Eur J Anaesthesiol 18, 811 (2001);
5. Turk J Med Sci 30, 517 (2000);
6. Br J Pharmacol 136, 995 (2002).

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

Cat #	Product Name	Unit Size
E12353	ER-Tracker™ Blue-White DPX *for live-cell imaging* *1 mM solution in DMSO*	20 x 50 µL
E34250	ER-Tracker™ Red (BODIPY® TR glibenclamide) *for live-cell imaging*	100 µg
E34251	ER-Tracker™ Green (BODIPY® FL glibenclamide) *for live-cell imaging*	100 µg

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 — 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

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

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
Technical Services: 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.