

Preparation of Dihydrorhod-2 AM

Introduction

The cell-permeant AM ester derivative of rhod-2 (R-1244, R-1245) has a net positive charge. This property promotes its sequestration into mitochondria in some cells, most likely via membrane potential-driven uptake. By reducing rhod-2 AM to the colorless, nonfluorescent dihydrorhod-2 AM, the discrimination between cytosolic and mitochondrially localized dye can be further improved.¹ The AM ester of dihydrorhod-2 exhibits Ca^{2+} -dependent fluorescence only after it is oxidized and its AM esters are cleaved to yield the rhod-2 indicator, processes that occur rapidly in the mitochondrial environment. Reduction of rhod-2 AM (R-1244, R-1245) to dihydrorhod-2 AM can be readily accomplished using the protocol given below.

Protocol

1.1 Dissolve 50 μg of rhod-2 AM (one vial of the set of 20 supplied as product R-1245) in 100 μL of anhydrous dimethylsulfoxide (DMSO).

1.2 Add a small excess of sodium borohydride (NaBH_4) either as a solid or as a methanol solution. The smallest amount of solid NaBH_4 that can be practicably transferred will provide a sufficient excess.

1.3 Incubate for 10 minutes or until the reaction mixture appears colorless, whichever is sooner.

1.4 Use the reaction solution in DMSO (about 0.4 mM dihydrorhod-2 AM) directly for cell loading according to usual AM ester loading protocols (see our product information sheets *Rhod-2* and *X-rhod-1 Calcium Indicators* (MP1244) or *Acetoxymethyl (AM) Esters* (G002)).

1.5 Dihydrorhod-2 AM will spontaneously and quite rapidly revert to the oxidized form, marked by reappearance of color in the stock solution. Therefore experiments using dihydro rhod-2 AM should be carried out as soon as possible after preparation.

References

1. Cell 82, 415 (1995).

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 Leiden, the Netherlands. All others should contact our Technical Assistance Department in Eugene, Oregon.

Please visit our Web site — www.probes.com — for the most up-to-date information

Molecular Probes, Inc.

PO Box 22010, Eugene, OR 97402-0469

Phone: (541) 465-8300 • Fax: (541) 344-6504

Customer Service: 7:00 am to 5:00 pm (Pacific Time)

Phone: (541) 465-8338 • Fax: (541) 344-6504 • order@probes.com

Toll-Free Ordering for USA and Canada:

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

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

Phone: (541) 465-8353 • Fax: (541) 465-4593 • tech@probes.com

Molecular Probes Europe BV

PoortGebouw, Rijnsburgerweg 10

2333 AA Leiden, The Netherlands

Phone: +31-71-5233378 • Fax: +31-71-5233419

Customer Service: 9:00 to 16:30 (Central European Time)

Phone: +31-71-5236850 • Fax: +31-71-5233419

eurorder@probes.nl

Technical Assistance: 9:00 to 16:30 (Central European Time)

Phone: +31-71-5233431 • Fax: +31-71-5241883

eurotech@probes.nl

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.

Several of Molecular Probes' products and product applications are covered by U.S. and foreign patents and patents pending. Our products are not available for resale or other commercial uses without a specific agreement from Molecular Probes, Inc. We welcome inquiries about licensing the use of our dyes, trademarks or technologies. Please submit inquiries by e-mail to busdev@probes.com. All names containing the designation ® are registered with the U.S. Patent and Trademark Office.

Copyright 2001, Molecular Probes, Inc. All rights reserved. This information is subject to change without notice.