

Acetoxymethyl (AM) and Acetate Esters

Table 1. Contents and storage information.

| Product | Amount | Concentration | Storage | Stability |
|--------------------------------------|--------|---------------|--|--|
| <i>Please refer to product label</i> | | | <ul style="list-style-type: none"> • $\leq -20^{\circ}\text{C}$ • Desiccated • Protected from light <p style="text-align: center;">or</p> <ul style="list-style-type: none"> • $\leq -20^{\circ}\text{C}$ • Desiccated <p style="text-align: center;"><i>(see the label)</i></p> | When stored as directed, product is stable for 6 months. |

Introduction

The acetoxymethyl (AM) ester derivatives of fluorescent indicators and chelators make up one of the most useful groups of compounds for the study of live cells. Modification of carboxylic acids with AM ester groups results in an uncharged molecule that can permeate cell membranes. Once inside the cell, the lipophilic blocking groups are cleaved by nonspecific esterases, resulting in a charged form that leaks out of cells far more slowly than its parent compound. Frequently, hydrolysis of the esterified groups is essential for binding of the target ion. In some cases (e.g., calcein AM), the AM ester is colorless and nonfluorescent until hydrolyzed. This property is useful in diagnosing spontaneous hydrolysis during storage. Acetate groups, used on many fluorescent indicators, are analogous to AM ester groups and should be treated similarly.

Guidelines for Use

Preparing the Stock Solutions

AM or acetate esters should be reconstituted only as required using high-quality, anhydrous dimethylsulfoxide (DMSO). Reagent-grade DMSO should be stored well sealed under argon or nitrogen, and desiccated; desiccant beads (e.g., molecular sieves) can be used for short-term storage. Dissolution of the pure AM or acetate esters in DMSO may be slow (particularly in the 1 mg sizes). Once prepared, DMSO stock solutions of AM or acetate esters should preferably be used within a short time period for one series of experiments. DMSO stock solutions must be kept anhydrous, since the solvent will readily take up moisture, leading to decomposition of the dye. These stock solutions should be stored well sealed, frozen, and desiccated. Many of our AM and acetate esters are available in small aliquots. Use of the prealiquoted product is strongly recommended.

It is advisable to keep the AM ester or acetate ester in as concentrated a stock as possible so that minimal amounts (ideally $\leq 0.1\%$) of DMSO are present in the loading solution.

Using Pluronic® F-127 Dispersing Agent

Since some AM esters (particularly SBF1, AM and PBFI, AM) are relatively insoluble in aqueous solutions, the low-toxicity dispersing agent Pluronic® F-127 is often used to facilitate cell loading. This nonionic detergent can be made up to a final concentration of 20% (w/v) in DMSO, and this solution can be used to prepare the dye stock. Gentle warming (~40°C) may assist in getting the detergent into DMSO. Pluronic® F-127 may decrease the stability of AM esters, so it should only be added to working stocks.

For the convenience of our customers, Molecular Probes offers Pluronic® F-127 in three forms: 1 mL of a 20% (w/v) solution in DMSO (P3000), 30 mL of a 0.2 µm–filtered 10% (w/v) solution in water (P6866) and 2 g solid (P6867).

Whether Pluronic® is used or not, it is advisable to keep the AM ester or acetate ester in as concentrated a stock as possible so that minimal amounts (ideally ≤ 0.1%) of DMSO are present in the loading solution.

Loading of Cells with AM or Acetate Esters

This is intended as an introduction only. Specific protocols for any particular dye and cell type should be obtained from the literature. As a rule, AM and acetate esters are used at a final working concentration of between 1 and 10 µM. Higher concentrations of weakly fluorescent indicators such as Fura Red™ and quin-2 may be required. The AM or acetate ester concentration should be kept as low as possible to reduce potential artifacts from overloading, including incomplete hydrolysis, compartmentalization, and toxic effects of hydrolysis by-products such as formaldehyde or acetic acid. Generally, loading times of between 15 minutes and 1 hour are sufficient, although probes such as SBF1, AM and PBFI, AM may require 1–4 hours. Loading may be done at a temperature that is optimal for the cells, although some investigators have reported greater degrees of compartmentalization at physiological temperatures than at room temperature. In addition to assisting in dye uptake, Pluronic® F-127 may help in reducing compartmentalization. To keep extracellular hydrolysis of the AM and acetate esters to a minimum, it is recommended that a loading buffer free of primary and secondary amines such as PBS be used. Cells should be washed in dye-free buffer after loading.

Assessing Dye Responsiveness

The following protocol is optional and can be used to test the responsiveness of the dye (for example following extended storage) or as a means to calibrate the ion response of the dye. Following the calibration or testing of the dye using this protocol, it should be disposed of and not loaded into cells.

We generally recommend that the separately available salt or free acid form of an indicator be used for calibrating the ion response. However, the following protocol for AM or acetate ester hydrolysis provides a less-preferred alternative and may also be useful to assess spontaneous hydrolysis during storage. *This procedure is not always successful for AM esters, probably because of the formation of formaldehyde in the reaction.*

- 1.1 Dissolve a small amount of the AM or acetate ester (e.g., 50 µg calcein AM) in 50 µL dioxane, DMSO, or other water-miscible solvent.
- 1.2 Add an equal volume of methanol.
- 1.3 Add 25 µL of 2 M KOH/water. If the dye is not in solution at this point, add more methanol.
- 1.4 Wait one hour.
- 1.5 Adjust pH to ~7 with HCl.

1.6 Test for fluorescent response. For example, to test a calcium indicator, dilute 5 μ L of the dye solution into 100 μ L of water, add this separately to high-calcium buffer and to low-calcium buffer.

1.7 If the dye does not respond properly, add more KOH/methanol to the dye solution and repeat steps 1.6 and 1.7.

Related Products

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

| Cat # | Product Name | Unit Size |
|--------|--|-------------------|
| P6867 | Pluronic® F-127 *low UV absorbance* | 2 g |
| P3000 | Pluronic® F-127 *20% solution in DMSO* | 1 mL |
| P6866 | Pluronic® F-127 *10% solution in water* *0.2 μ m filtered* | 30 mL |
| P10020 | PowerLoad™ 100X concentrate | 5 mL |
| P36400 | Probenecid, water soluble | 10 \times 77 mg |

Contact 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
probetech@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

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.

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 No. 223: Labeling and Detection Technology

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) 335-0354.

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 2008, Molecular Probes, Inc. All rights reserved. This information is subject to change without notice.