

Calcein violet, AM

Table 1. Contents and storage information.

Material	Amount	Storage	Stability
Calcein violet AM	20 vials, 25 µg each	<ul style="list-style-type: none"> • ≤-20°C • Desiccate * • Protect from light 	When stored as directed, product should remain stable for at least 1 year.

* Calcein violet AM may hydrolyze if exposed to moisture

Number of labelings: At the recommended reagent concentrations and volumes, this kit contains sufficient material to perform approximately 1,000 assays using flow cytometry.

Approximate fluorescence excitation/emission maxima: Calcein violet AM: 400/452 nm

Introduction

Cell vitality as measured by intracellular esterase activity is a recognized parameter of cell health. Live cells are distinguished by the presence of ubiquitous intracellular esterase activity, determined by the enzymatic conversion of the virtually non-fluorescent cell-permeant calcein violet AM to the intensely fluorescent calcein violet (ex/em 400/452 nm, Figure 1). Calcein violet AM is an optimal dye for this application; utilizing the violet laser allows other laser lines to be used with more conventional markers.

Principle of the Method

The acetoxymethyl (AM) ester derivatives of fluorescent indicators 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 is retained in cells to a much greater extent than its parent compound. The calcein violet AM ester is colorless and non fluorescent until hydrolyzed. The polyanionic dye calcein violet is well retained within live cells, producing an intense uniform violet fluorescence in live cells (ex/em 400/452 nm, Figure 2).

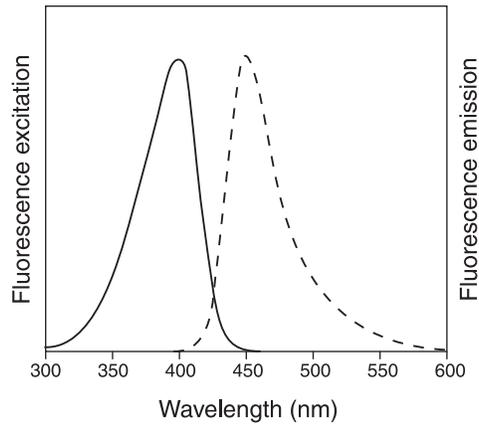


Figure 1. Fluorescence excitation and emission spectra for the Calcein violet stain in PBS, pH 7.2.

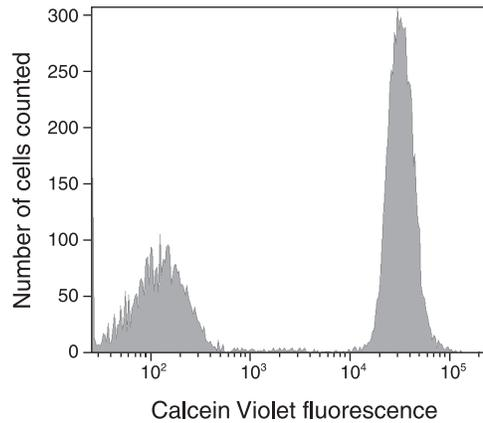


Figure 2. A mixture of heat-killed and untreated Jurkat cells (human leukemia T-cell) was stained according to the protocol in the Calcein violet, AM kit. Cells were analyzed using a flow cytometer equipped with a 405 nm laser and a 450/50 bandpass filter.

Before You Begin

Material Required but Not Provided

- Dimethylsulfoxide (DMSO): reconstitute calcein violet AM using high-quality, anhydrous DMSO.

Working with the Calcein Violet AM Stock Solution

Allow the vial to warm to room temperature before opening. Calcein violet AM is susceptible to hydrolysis when exposed to moisture. Once prepared, DMSO stock solutions of calcein violet AM should preferably be used within a short time period for one set of experiments, while aqueous working solutions containing calcein violet AM should be prepared immediately prior to use and used within one day.

Caution

Hazards posed by these stains have not been fully investigated. DMSO is known to facilitate entry of organic molecules into tissue. These reagents should be handled using equipment and practices appropriate for the hazards posed by such material. Please dispose of reagents in compliance with all pertaining local regulations.

Experimental Protocols

For Flow Cytometry This flow cytometry protocol has been optimized using Jurkat cells (human T-cell leukemia line) at a concentration of 1×10^6 cells/mL. Use of other cell types or other cell concentrations may require optimization of staining. If another staining reaction is to be performed on the sample, the user must determine the optimal staining sequence for the two procedures (See Figure 3).

- 1.1 Allow one vial of calcein violet AM to come to room temperature.
- 1.2 Prepare a 1 mL suspension of cells with $0.1\text{--}5 \times 10^6$ cells/mL for each assay. Cells may be suspended in medium or buffer.
- 1.3 Add 42 μL high-quality, anhydrous DMSO to one vial calcein violet AM to prepare a stock solution. Add 40 μL of this stock solution to 1.25 mL buffer or media to make a working solution of calcein violet AM. Use this working solution within one day.
- 1.4 Add 5 μL of working solution to each mL cell suspension. Mix the sample.
- 1.5 Incubate the cells for 20–30 minutes on ice or at room temperature.
- 1.6 Analyze the stained cells by flow cytometry using violet (~ 405 nm) excitation and violet fluorescence emission (~ 450 nm).

For Fluorescence Microscopy

This microscopy protocol has been optimized using bovine pulmonary artery endothelial cells (BPAEC). Use of other cell types or other cell concentrations may require optimization of staining. If another staining reaction is to be performed on the sample, the user must determine the optimal staining sequence for the two procedures.

- 2.1 Allow one vial of calcein violet AM to come to room temperature.
- 2.2 Add 42 μL high-quality, anhydrous DMSO to one vial calcein violet AM, yielding a 1 mM stock solution. Once prepared, the DMSO stock solution should be used within a short time period for one series of experiments.
- 2.3 Prepare a working solution of dye in buffer which is free of serum. We recommend Hanks Balanced Salt Solution, HBSS (Invitrogen Cat. no. 14025-092) containing stain at a concentration range of 250 nM to 1 μM .
- 2.4 Adherent cells may be stained on coverslips after rinsing once with HBSS to remove residual serum present in the culture medium. Add a sufficient amount of stain solution to adequately cover the adherent cells. Cells in suspension should be pelleted by centrifugation and washed once in HBSS. Incubate for 30 minutes at 37°C .
- 2.5 Observe the samples in staining solution using a fluorescence microscope.

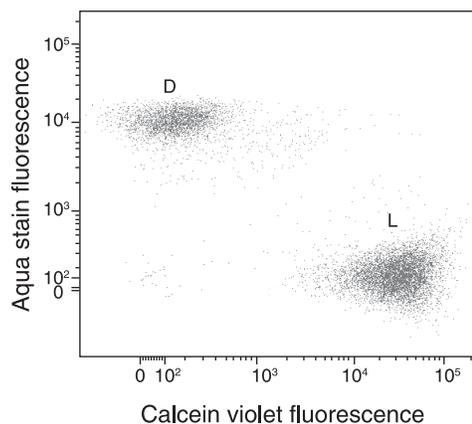


Figure 3. Staining pattern of a mixture of heat-killed and untreated Jurkat cells (human leukemia T-cell). Jurkat cells were stained according to the protocol in the LIVE/DEAD® Violet Viability/Vitality Kit, which combines the calcein violet fluorescence with the dead cell aqua-fluorescent reactive dye. Cells were analyzed using a flow cytometer equipped with a 405 nm laser and 450/50 bandpass for calcein violet labeling live cells (L) and 525/50 bandpass for the aqua-fluorescent reactive dye labeling dead cells (D).

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

Cat. no.	Product Name	Unit Size
C34858	Calcein violet, AM *for flow cytometry* *for 405 nm excitation* *special packaging*	20 × 25 µg

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
 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

For country-specific contact information, visit www.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 2010, Molecular Probes, Inc. All rights reserved. This information is subject to change without notice.