

## MitoProbe™ DiIC<sub>1</sub>(5) Assay Kit for Flow Cytometry (M34151)

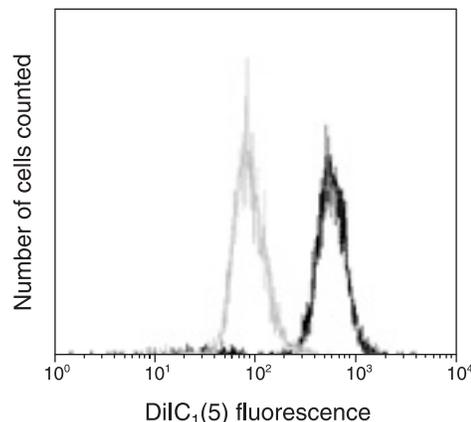
### Quick Facts

#### Storage upon receipt:

- 2–6°C
- Protect from light
- Do not freeze

**Ex/Em:** 638/658 nm

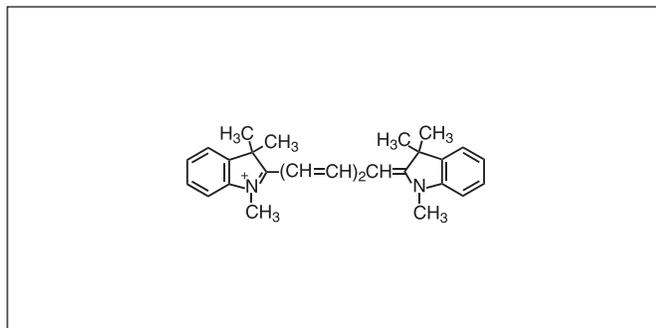
**Number of Assays:** 100, based on labeling volumes of 1.0 mL



**Figure 2.** Decrease in DiIC<sub>1</sub>(5) fluorescence with the addition of CCCP. Jurkat cells were stained with 50 nM DiIC<sub>1</sub>(5) alone (black line) or in the presence of 50 μM CCCP (gray line).

### Introduction

Cationic cyanine dyes have been shown to accumulate in cells in response to membrane potential<sup>1</sup> and membrane potential changes have been studied in association with apoptosis.<sup>2-3</sup> The MitoProbe™ DiIC<sub>1</sub>(5) Assay Kit provides solutions of the cyanine dye DiIC<sub>1</sub>(5) (1,1',3,3,3',3'-hexamethylindodicarbocyanine iodide, Figure 1) and CCCP (carbonyl cyanide 3-chlorophenylhydrazone), for the study of mitochondrial membrane potential. DiIC<sub>1</sub>(5) penetrates the cytosol of eukaryotic cells. At concentrations below 100 nM, the dye accumulates primarily in mitochondria with active membrane potentials. DiIC<sub>1</sub>(5) stain intensity decreases when cells are treated with reagents that disrupt mitochondrial membrane potential, such as CCCP (Figure 2).



**Figure 1.** Structure of DiIC<sub>1</sub>(5), molecular weight: 510.46.

Cells stained with DiIC<sub>1</sub>(5) can be visualized by flow cytometry with red excitation and far red emission. The reagent can be paired with other reagents, such as blue-excited propidium iodide and annexin V–Alexa Fluor® 488 dye (Vybrant® Apoptosis Assay Kit #2, V13241), for multiparametric study of vitality and apoptosis (Figure 3). Combining DiIC<sub>1</sub>(5) dye with an annexin V conjugate results in superior resolution of subpopulations when compared to results obtained from other commonly used dyes.

### Materials

#### Kit Contents

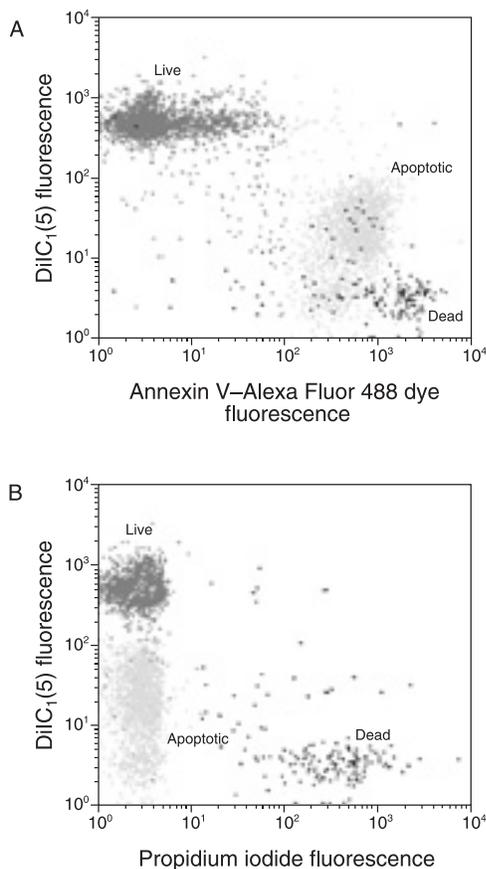
- DiIC<sub>1</sub>(5), 625 μL of 10 μM DiIC<sub>1</sub>(5) in DMSO
- CCCP, 125 μL of 50 mM CCCP in DMSO

#### Storage and Handling

Upon receipt, components should be stored at 2–6°C. DO NOT FREEZE. Before opening, each vial must be at room temperature. When stored properly, both the DiIC<sub>1</sub>(5) and CCCP solutions should be stable for at least twelve months.

#### Spectral Characteristics

The approximate excitation and emission peaks of DiIC<sub>1</sub>(5) are 638 nm and 658 nm, respectively. Cells labeled with DiIC<sub>1</sub>(5) can be analyzed by flow cytometry using 633 nm excitation and far red emission, and by fluorescence microscopy using standard filters for Alexa Fluor® 633 dye.



**Figure 3.** Camptothecin-treated Jurkat cells stained with DiIC<sub>1</sub>(5) and annexin V–allophycocyanin. Cells were incubated for 3 hours with 10 μM camptothecin at 37°C, 5% CO<sub>2</sub>, then stained with DiIC<sub>1</sub>(5), annexin V–Alexa Fluor<sup>®</sup> 488 dye, and propidium iodide. In Panel A, most of the annexin V–positive cells show reduced DiIC<sub>1</sub>(5) fluorescence. In Panel B, propidium iodide–positive cells show low to very low DiIC<sub>1</sub>(5) fluorescence.

## Experimental Protocol

The following protocol describes introducing DiIC<sub>1</sub>(5) reagent into the cultured cells and analyzing the stained cells by flow cytometry. Suggested initial conditions may require modifications because of differences in cell types and culture conditions. The concentration of probe for optimal staining will vary depending upon the application. A concentration range should be tested starting around 50 nM DiIC<sub>1</sub>(5). CCCP controls should be used to confirm that the DiIC<sub>1</sub>(5) response is sensitive to changes in membrane potential.

### Labeling Cells with DiIC<sub>1</sub>(5)

Before beginning the experiment, ensure that the vials of DiIC<sub>1</sub>(5) and CCCP have equilibrated to room temperature.

**1.1** For each sample, suspend cells in 1 mL warm medium, phosphate-buffered saline, or other buffer at approximately 1 x 10<sup>6</sup> cells/mL.

**1.2** For the control tube, add 1 μL of 50 mM CCCP (supplied with the kit, 50 μM final concentration) and incubate the cells at 37°C for 5 minutes.

**Note:** CCCP can be added simultaneously with DiIC<sub>1</sub>(5). Titration of the CCCP may be required for optimal results with each cell type.

**1.3** Add 5 μL of 10 μM DiIC<sub>1</sub>(5) (supplied with the kit, 50 nM final concentration) and incubate the cells at 37°C, 5% CO<sub>2</sub>, for 15 to 30 minutes. If performing additional labeling, for example with an annexin V conjugate, follow the protocol below, beginning with step 2.1. If no additional staining is to be performed, proceed with step 1.4.

**1.4** OPTIONAL: Wash cells once by adding 2 mL of warm phosphate-buffered saline (PBS) or other buffer to each tube of cells.

**1.5** Pellet the cells by centrifugation.

**1.6** Resuspend by gently flicking the tubes. Add 500 μL PBS (or other suitable buffer) to each tube.

**1.7** Analyze on a flow cytometer with 633 nm excitation using emission filters appropriate for Alexa Fluor<sup>®</sup> 633 dye (Figure 2). Gate on the cells, excluding debris. Using the CCCP-treated sample, perform standard compensation.

### Additional Labeling with an Annexin V Conjugate and Propidium Iodide

It is possible to label the DiIC<sub>1</sub>(5)-stained cells with other markers for apoptosis or viability, as long as the fluorescence emission of the additional label is spectrally resolved from DiIC<sub>1</sub>(5). The example below is a protocol for labeling with an annexin V–Alexa Fluor<sup>®</sup> 488 conjugate and propidium iodide.

**2.1** After step 1.3 (above), wash cells once by adding 2 mL of warm phosphate-buffered saline or other buffer to each tube of cells.

**2.2** Pellet the DiIC<sub>1</sub>(5)-stained cells and resuspend in 100 μL of 1X annexin binding buffer (10 mM HEPES, 140 mM NaCl, and 2.5 mM CaCl<sub>2</sub>, pH 7.4).

**2.3** Add 5 μL annexin V conjugate (e.g. annexin V–Alexa Fluor<sup>®</sup> 488 conjugate, V13241).

**Note:** 5 μL is appropriate for annexin V conjugates from Molecular Probes. Conjugates purchased from other suppliers may require a different volume to be effective.

**2.4** Add 1 μL of a 100 μg/mL propidium iodide solution (V13241, prepared according to instructions accompanying that kit).

**2.5** Incubate the samples at 37°C for 15 minutes. (37°C is important to maintain membrane potential.)

**2.6** Add 400 μL annexin binding buffer.

**2.7** Analyze on a flow cytometer with 488 nm and 633 nm excitation using emission filters appropriate for Alexa Fluor<sup>®</sup> 488, R-phycoerythrin, and Alexa Fluor<sup>®</sup> 633 dye (Figure 3).

---

## References

1. Proc Natl Acad Sci USA 76, 5728 (1979); 2. Meth Cell Biol 63, 467 (2001); 3. Exp Cell Res 214, 323 (1994).

---

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

Cat #	Product Name	Unit Size
M34151	MitoProbe™ DiIC <sub>1</sub> (5) Assay Kit *for flow cytometry* *100 assays* .....	1 kit

---

## 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](http://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](mailto: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](mailto: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](mailto:euroinfo@invitrogen.com)

Technical Services: [eurotech@invitrogen.com](mailto: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.