



FxCycle™ Violet Stain

Catalog no. F10347

Table 1. Contents and storage information.

Material	Amount	Storage*	Stability
FxCycle™ Violet stain	5 × 100 μg	2–6°C Desiccate Protect from light	When stored as directed the product is stable for at least 1 year.

Number of assays: Sufficient material is supplied for 500 assays based on the protocol below.

Approximate fluorescence excitation/emission maxima: FxCycle™ Violet stain: 358/461 nm, bound to DNA.

Introduction

Analysis of nucleic acids is a common application of flow cytometry. Measurement of DNA content allows the study of cell populations in various phases of the cell cycle as well as analysis of DNA ploidy. In a given population, cells are distributed among three major phases of cell cycle: G0/G1 phase (one set of paired chromosomes per cell), S phase (DNA synthesis with variable amount of DNA), and G2/M phase (two sets of paired chromosomes per cell, prior to cell division).¹⁻⁴ DNA content can be measured using fluorescent DNA stains that exhibit emission signals proportional to the DNA mass. Flow cytometric analysis of these stained populations is then used to produce a frequency histogram that reveals the various cell cycle phases.

Univariate DNA content analysis is an established assay method and is widely used for studies in oncology, cell biology, and molecular biology. Using flow cytometry, multicolor cell cycle studies are possible, and it is advantageous to analyze DNA content on alternative lasers to preserve the common 488 nm laser for other markers. Well suited for the popular violet laser line, FxCycle™ Violet stain (4', 6-diamidino-2-phenylindole, dihydrochloride) can also be used with UV excitation. With DNA content measurement on the violet laser, other parameters such as cyclins, cyclin-dependent kinases, cell cycle checkpoints, nuclear proteins, and proliferation markers can be measured on the familiar 488 nm laser. FxCycle™ Violet stain preferentially stains dsDNA; it appears to associate with AT clusters in the minor groove.⁵ Binding of FxCycle™ Violet stain to dsDNA produces a ~20-fold fluorescence enhancement.⁶

For long-term storage, the dye is supplied as five vials of 100 μg solid dye and stored at 2–6°C; individual vial stock solution in deionized water is also stored at 2-6°C for convenience.

Materials Required but Not Provided

- · Deionized water
- Reagents for fixing cells such as alcohol or formaldehyde
- Reagents for permeabilizing cells such as Triton® X-100
- Buffer such as phosphate buffered saline (PBS)

Caution

The hazards posed by FxCycle™ Violet stain has not been fully investigated. The stain is a known mutagen and may cause sensitization by inhalation and skin contact, and is irritating to eyes, respiratory system, and skin. Do not breath dust. In case of contact with eyes, rinse immediately with plenty of water and seek medical advice. Wear suitable protective clothing, safety glasses, and gloves. Avoid contact with skin and eyes. Dispose of the reagents in compliance with all pertaining local regulations.

Preparing Stock Solution

To make a 1 mg/mL stock solution of FxCycle $^{\!\scriptscriptstyle{\text{\tiny{TM}}}}$ Violet stain, add 100 μL deionized water to one vial of the stain. Mix well. Store this solution at 2–6°C, **protected from light**. When stored as directed, this FxCycle[™] Violet stock solution is stable for at least six months.

Spectral Characteristics

The fluorescence excitation and emission spectra of the FxCycle™ Violet stain are shown in Figure 1. The spectra were obtained from samples of the stain bound to DNA with fluorescence excitation and emission maxima of 358/461 nm respectively. FxCycle™ Violet stain may be used with the violet 405 nm excitation laser commonly found on flow cytometers, as well as UV excitation.

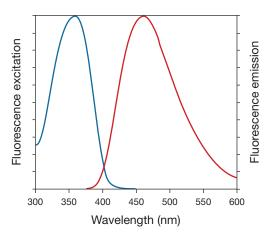


Figure 1. Fluorescence excitation and emission spectra of FxCycle™ Violet stain bound to dsDNA.

Experimental Protocol

The following staining procedure was developed using the Jurkat T-cell leukemia cell line, but can be adapted for any cell type. Fixative, permeabilization reagent, cell density, cell type variations, and other factors may influence staining.

In initial experiments, try a range of stain concentrations to determine the concentration that yields optimal staining for the given cell type and experimental conditions. All fixative should be removed from cells before proceeding with cell staining, however staining with FxCycle[™] Violet stain may be done concurrent with the addition of a permeabilization reagent if desired. For a given experiment, each flow cytometry sample should contain the same number of cells, as sample-to-sample variation in cell number leads to significant differences in fluorescence signal.

If FxCycle™ Violet stain is used in combination with other stains for multicolor applications, apply the other stain(s) to the sample first, following all manufacturers instructions, including wash steps. FxCycle™ Violet stain should be the last stain applied to the sample, and do not wash samples prior to flow cytometric analysis.

General Guidelines

For optimal DNA content cell cycle analysis, follow these general guidelines:

- Eliminate cell clumps and aggregates from the cell suspension before staining
- Validate flow cytometry instrument performance on the day of use
- · Use linear amplification for DNA content
- Use low flow rate for acquisition
- Collect adequate number of events for the intended application
- Cells must be fixed before staining with FxCycle[™] Violet stain for DNA content cell cycle
- Do not wash cells after staining with FxCycle[™] Violet stain

Staining Procedure

- **1.1** Harvest the cell sample(s).
- **1.2** Fix cells according to your preferred protocol.
- 1.3 Wash cells. Remove all fixative from cells before proceeding with the cell staining.
- 1.4 Using an appropriate buffer, adjust the sample cell concentration to be 1×10^6 cells/mL.
- 1.5 Prepare flow cytometry samples each containing 1 mL cell suspension. Optional permeabilization reagent may be added.
- **1.6** Add 1 μL of FxCycle™ Violet stain to each flow cytometry sample and mix well.
- 1.7 Incubate flow cytometry tubes for 30 minutes at room temperature or 2–6°C, protected from light.
- 1.8 Analyze samples without washing in a flow cytometer, using 405 nm excitation and emission collected in a 450/50 bandpass or equivalent.

Example of results obtained using FxCycle™ Violet stain for DNA content cell cycle analysis is shown in Figure 2.

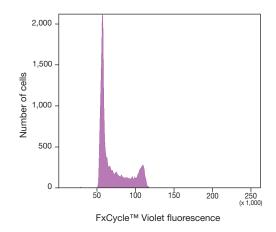


Figure 2. Histogram of HL-60 promyeloblast cells stained with FxCycle™ Violet stain showing DNA content distribution. HL-60 cells were fixed overnight with alcohol, washed, and then resuspended in 0.1% Triton® X-100/PBS/1% BSA before staining with FxCycle™ Violet stain for 30 minutes at room temperature. G0/G1 and G2/M phase histogram peaks are separated by the S-phase distribution. Analysis was performed at 405 nm excitation with a 450/50 bandpass filter.

References

1. Current Protocols in Cytometry, 7.0.1–7.27.7 (2004); 2. Practical Flow Cytometry, 4th Ed., Shapiro HM, Ed. (2003); 3. Methods Mol Biol 281, 301 (2004); 4. Cytometry A 58, 21 (2004); 5. Biochemistry 26, 4545 (1987); 6. Biochem Biophys Res Commun 170, 270 (1990).

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

Cat. no.	Product Name	Unit Size			
F10347	FxCycle™ Violet Stain *for flow cytometry* *500 assays* *DAPI*	1 set			
F10348	FxCycle™ Far Red Stain *for flow cytometry* *500 assays*	1 set			
Related Products					
L10120	LIVE/DEAD® Fixable Far Red Dead Cell Stain Kit *for 633 or 635 nm excitation* *200 assays*	1 kit			
L23101	LIVE/DEAD® Fixable Green Dead Cell Stain Kit *for 488 nm excitation* *200 assays*	1 kit			
L23102	LIVE/DEAD® Fixable Red Dead Cell Stain Kit *for 488 nm excitation* *200 assays*	1 kit			
L23105	LIVE/DEAD® Fixable Blue Dead Cell Stain Kit *for UV excitation* *200 assays*	1 kit			
L34955	LIVE/DEAD® Fixable Violet Dead Cell Stain Kit *for 405 nm excitation* *200 assays*	1 kit			
L34957	LIVE/DEAD® Fixable Aqua Dead Cell Stain Kit *for 405 nm excitation* *200 assays*	1 kit			
GAS-003	Fixation and Permeabilization, $1 \times 5 \text{ ml}$	50 tests			
GAS-004	Fixation and Permeabilization, $4 \times 5 \text{ ml}$	200 tests			
GAS001S-10	0 Fixation Medium - Bulk, (MEDIUM A), 1 × 100 ml Fixation Medium	1000 tests			
GAS002S-10	0 Permeabilization Medium - Bulk, (MEDIUM B), 1 \times 100 ml	1000 tests			
10010-049	Phosphate Buffered Saline (PBS) 7.2 (1X), liquid	\times 500 mL			
20012-050	Phosphate Buffered Saline (PBS) 7.4 (1X), liquid	\times 500 mL			
14025-092	Hanks' Balanced Salt Solution (HBSS) (1X), liquid, contains calcium and magnesium, but no phenol red	500 mL			
14170-112	Hanks' Balanced Salt Solution (HBSS) (1X), liquid, contains no calcium chloride, magnesium chloride, or magnesium sulfate	500 mL			
14175-095	Hanks' Balanced Salt Solution (HBSS) (1X), liquid, contains no calcium chloride, magnesium chloride, magnesium sulfate,				
	or phenol red	500 mL			
24020-117	Hanks' Balanced Salt Solution (HBSS) (1X), liquid, contains calcium and magnesium	500 mL			
12091-039	RNAse A (20 mg/ml)	25 mL			

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 an extension of the product of tfrom 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 transfer (a) this product or its components to a third party 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 transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party 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 transfer (b) its components or (c) materials made using this product or its components to a third party or otherwise transfer (b) its components or (c) materials made using the components of the component of the compone erwise 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 2009, Molecular Probes, Inc. All rights reserved. This information is subject to change without notice.