

FilmTracer™ FM® 1-43 Green Biofilm Cell Stain

Catalog no. F10317

Table 1. Contents and storage information.

Material	Amount	Storage	Stability
FilmTracer™ FM® 1-43 green biofilm cell stain	1 mg	 2–25°C Desiccate Protect from light	When stored as directed, the product is stable for at least 6 months.

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

Approximate fluorescence excitation/emission maxima: FilmTracer™ FM® 1-43 green biofilm cell stain: 472/580 nm.

Introduction

Bacterial biofilms present a unique set of challenges for fluorescent staining and subsequent imaging. A typical biofilm not only exhibits heterogeneous thickness throughout the surface, placing stringent restrictions on stain penetration, but also contains regions of widely varying environmental conditions. Evidence suggests that bacterial cells exist in various physiological states within these biofilm microenvironments. Furthermore, biofilms contain many undefined components (e.g., the extracellular polymeric matrix) that differ with species and conditions.

FM[®] dyes are lipophilic styryl compounds used in a wide variety of studies involving plasma membrane and vesiculation. The water-soluble FM® dyes are virtually nonfluorescent in aqueous media, and are believed to insert into the surface membrane where they become intensely fluorescent. FM® 1-43 stain has been used successfully to stain the cell bodies specifically in a complex biofilm mileu, including Pseudomonas aeruginosa, Escherichia coli, Staphylococcus sp., Acidothiobacillus caldus, and Vibrio cholerae.^{1,2}

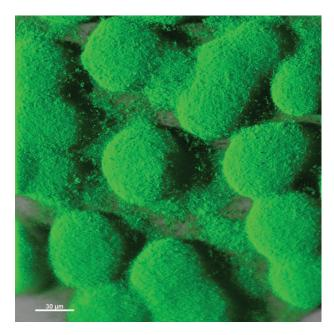


Figure 1. FM® 1-43 applied to a Pseudomonas aeruginosa biofilm. FM® 1-43 appears to bind to the cell membrane, and has been shown to work equally well on Staphylococcus epidermidis, Pseudomonas aeruginosa, and Escherichia coli, exhibiting exceptional cell specificity in each case. The image was obtained using a Leica TCS-SP2 AOBS confocal microscope and a 63×/0.9 NA water immersion objective.

Before You Begin

Materials Required but Not Provided

- Biofilm samples: Biofilms may be grown on coupons in a biofilm reactor, as colony biofilms, in flow-cell system, or in drip-flow reactors. This protocol describes staining biofilms grown on glass coupons in a CDC reactor. For more information on the CDC reactor or other reactor types, refer to the BioSurface Technologies website (www.imt.net/~mitbst/Products.html), contact Center for Biofilm Engineering, Montana State University, Bozeman, Montana, or refer to the standard protocols outlined in the following ASTM methods: ASTM E2647, ASTM E2562, ASTM E2196...
- Fluorescence microscope with appropriate FITC excitation/emission filters
- 0.2 µm filter-sterilized water
- Staining dishes (e.g., 60-mm dish, 6-well plate, etc.)
- Anhydrous dimethylsulfoxide (DMSO)

Preparing Reagents

Add 100 μ L DMSO to the contents of the vial. Mix thoroughly until dissolved. The stock solution in DMSO may be stored at 4°C; do not store the staining solution that is diluted in water. Dispose of dyes according to local regulations.

Caution

FilmTracer™ FM® 1-43 stain is potentially harmful, and may cause irritation by inhalation. Avoid prolonged or repeated exposure. Wear appropriate gloves, protective clothing and eyewear and follow safe laboratory practices. Wash thoroughly after handling. If eye or skin contact occurs, wash affected area with water for 15 minutes and seek medical advice. If inhaled, move individual to fresh air and seek medical advice. If swallowed, seek immediate medical advice.

Use appropriate protective equipment and methods to clean up spilled substances promptly. Absorb spill onto an appropriate material. Collect and dispose of all waste in accordance with applicable laws.

Experimental Protocols

Guidelines for Staining

- We recommend performing staining in water as the phosphates in buffers may interfere with fluorescent staining.
- If you need to stain and image multiple samples, do not stain more than two samples at a time. Evidence suggests that, in many cases, stain might be drawn from cells over time as they sit in water. Stagger staining, so that samples are stained, rinsed, and imaged following the same schedule. Image immediately following rinsing.
- For imaging biofilm on CDC reactor coupons, use glass coupons only. In particular, avoid polycarbonate coupons for imaging purposes as polycarbonate is autofluorescent, and the rough surface interferes with imaging.
- If you follow the procedure below, you do not need to use fixatives on the biofilm.

Staining Procedure

The protocol below describes how to stain biofilms grown on glass CDC reactor coupons. For any other growth surface, you may need to adjust the staining volumes.

- 1.1. Prepare staining solutions by diluting 10 µL stock solution into 990 µL DMSO, followed by diluting 100 µL into 0.9 mL filter-sterilized water. We recommend a final staining concentration of $1 \mu g/mL$.
- 1.2. Add 200 μL (or appropriate volume) of staining solution onto the biofilm sample. Add the stain very gently so as not to disturb the biofilm. It is important to immediately add the stain before the biofilm dries.
- **1.3.** Incubate the sample for up to 30 minutes at room temperature, **protected from light**.
- 1.4. Rinse the sample gently with filter-sterilized water. Remove all excess stain and rinse water from the base of the support material.
- 1.5. For best results with reactor coupons, place coupon in a 60-mm dish, fill the dish with filtersterilized water to cover the coupon surface by 1-3 mm, and observe on the microscope using a 40X 0.7NA 3.3 mm WD water objective or a 63X 0.9NA 2.2 mm WD water immersion objective.

References

1. Microscopy Today 16, 18 (2008); 2. J Bacteriol 190, 311 (2008).

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

Cat. no.	Product Name	Unit Size
F10317	FilmTracer™ FM® 1-43 green biofilm cell stain	1 mg
Related Produ	icts	
F10318	FilmTracer™ SYPRO® Ruby biofilm matrix stain.	200 mL
F10319	FilmTracer™ calcein red-orange biofilm stain	0 × 50 μg
F10320	FilmTracer™ calcein violet biofilm stain	0 × 25 μg
F10322	FilmTracer™ calcein green biofilm stain	0 × 50 μg
F35355	FM® 1-43FX *fixable analog of FM® 1-43 membrane stain*	\times 100 μg
L10316	FilmTracer™ LIVE/DEAD™ Biofilm Viability Kit	1 kit
S34854	SYTO® 9 green fluorescent nucleic acid stain	100 μL

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 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 composite the purchase of the purchased amount of the product and composite the purchase of the purchased amount of the product and composite the purchase of the purchased amount of the product and composite the purchased amount of the purchased amount of the product and composite the purchased amount of the purchased amountnents 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 (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 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 this product or its components to a third party or otherwise transfer (c) materials made using this product or its component (c) materials made using the component (c) materials erwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyermay 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. For products that are subject to multiple limited use label licenses, the most restrictive terms apply. Invitrogen Corporation will not assert a claim against the buyer of infringement of patents that are owned or controlled by Invitrogen Corporation and/or Molecular Probes, Inc. which cover this product 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 $\ensuremath{^\circ}$ 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.

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 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. For products that are subject to multiple limited use label licenses, the most restrictive terms apply. Invitrogen Corporation will not assert a claim against the buyer of infringement of patents that are owned or controlled by Invitrogen Corporation and/or Molecular Probes, Inc. which cover this product 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 the product of the proneither this product. If the purchaser is not will ling to accept the manufacture of such product. If the purchaser is not will ling to accept the purchaser is not will ling to accept the purchaser in the purchaser is not will ling to accept the purchaser in the purchaser is not will ling to accept the purchaser in the purchaser is not will line to accept the purchaser in the purchaser is not will line to accept the purchaser in the purchaser in the purchaser is not will line to accept the purchaser in the pulimitations 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.