

FluoroMyelin™ Fluorescent Myelin Stains

FluoroMyelin™ Green Fluorescent Myelin Stain (F34651)

FluoroMyelin™ Red Fluorescent Myelin Stain (F34652)

Quick Facts

Storage upon receipt:

- 4°C
- Protect from light

Ex/Em:

- FluoroMyelin Green fluorescent myelin stain:
479/598 nm
- FluoroMyelin Red fluorescent myelin stain:
558/654 nm

Introduction

The FluoroMyelin™ Red and FluoroMyelin Green fluorescent myelin stains enable quick and selective labeling of myelin in brain cryosections. Visualization of myelin is useful for understanding myelin distribution, identifying brain structures and relative location of other labels, and mapping a particular section within the brain. Traditional methods require the use of antibodies, such as anti-myelin basic protein, or chromogenic (transmitted-light) methods, such as the Loyez method or Schmued's gold chloride technique, all of which are time consuming, requiring multiple steps over one to three days. The FluoroMyelin stains, in contrast, require only a single 20-minute labeling step plus washes. These stains can be used in combination with antibodies and other dyes, and with standard histochemical methods for cryosectioned material. They are believed to work via lipophilic affiliation, taking advantage of the high lipid content of myelin in axon sheaths, but they are more selective than traditional lipophilic dyes such as DiI. The lipophilic nature of FluoroMyelin stains does, however, lead to faint labeling of all cell membranes. In addition to being used on their own, FluoroMyelin stains can be combined with other dyes in a single labeling step as in our BrainStain™ Imaging Kit, which includes NeuroTrace® 530/615 red fluorescent Nissl stain and DAPI nucleic acid stain for three-color visualization.

Materials

Contents

The FluoroMyelin Green and FluoroMyelin Red stains are provided in a volume of 1 ml. Each is a 300X solution in water.

Storage

Upon receipt, the stains should be stored at 4°C, upright and protected from light. The stock solutions are stable for at least 6 months if stored properly.

Spectral Characteristics

A standard FITC filter set is suitable for imaging FluoroMyelin Green stain (excitation/emission maxima ~479/598 nm). There should be little or no bleedthrough into standard DAPI, TRITC, or far-red filter sets.

A standard TRITC or Texas Red® filter set is suitable for imaging FluoroMyelin Red stain (excitation ~558/654 nm). There should be little or no bleedthrough into standard DAPI, FITC, or far-red filter sets.

Protocol for Staining Mouse Brain Cryosections

This protocol has been optimized for staining 12 µm mouse brain cryosections on Superfrost Plus slides (Erie Scientific Co.), using 200 µl of staining solution within a PAP pen well at room temperature. Optimal staining conditions may vary slightly depending on sample conditions and should be determined by the user.

1. Rehydrate and permeabilize specimens. Bring the tissue sections on slides to room temperature, then rehydrate them in either PBS¹ or PBT (PBS + 0.2% TRITON X-100) for at least 20 minutes. Permeabilization with TRITON X-100 is not necessary for staining with FluoroMyelin stains, but is likely to be necessary for other counterstains or antibodies to be used.

2. Prepare staining solution. Prepare the staining solution by diluting the stock solution 300-fold into PBS.

3. Stain specimens. Flood the sections with staining solution and stain them for 20 minutes at room temperature.

4. Wash. When staining is complete, remove the solution, rinse in PBS, and wash 3 times for 10 minutes each with PBS.

5. Counterstain and mount. At this point the sections can be counterstained as necessary, or mounted with an aqueous antifade mounting medium such as ProLong® or ProLong Gold antifade reagent.

Note

1. PBS (phosphate-buffered saline) = 137 mM NaCl, 2.7 mM KCl, 10 mM phosphate ($\text{Na}_2\text{HPO}_4/\text{KH}_2\text{PO}_4$), pH 7.4

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

Cat #	Product Name	Unit Size
F34651	FluoroMyelin™ Green fluorescent myelin stain *solution in water*	1 ml
F34652	FluoroMyelin™ Red fluorescent myelin stain *solution in water*	1 ml

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 — 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

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

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.