

SlowFade® Antifade Kit

Quick Facts

Storage upon receipt:

- Room temperature

Introduction

When exposed to excitation light, all fluorescent dyes fade (photobleach). The photobleaching is dependent on: 1) the intensity; and 2) the duration of illumination. The photon output of a dye represents the average number of cycles of excitation followed by fluorescence emission that the dye undergoes before it is irreversibly photobleached. The average photon output is defined by the ratio of the probability that the dye will fluoresce (fluorescence quantum efficiency or Q_f) and the probability that it will photoreact irreversibly to become a nonfluorescent species (photobleaching quantum efficiency or Q_b). For example, fluorescein, which is very photolabile, has a Q_f/Q_b of about 30,000 in alkaline solution. Both Q_f and Q_b are properties of the dye that may be affected significantly by the dye's environment. The primary environmental influence on Q_b is the presence of singlet oxygen and free radical species. The main purpose of any antifade reagent is to sustain dye fluorescence. This is usually accomplished by inhibiting the generation and diffusion of reactive oxygen species, thereby reducing Q_b (preferably without any accompanying decrease in Q_f , so that fluorescence will persist).

The active ingredient in Molecular Probes' SlowFade® Antifade Kit is 1,4-diazabicyclo[2.2.2]octane (DABCO), which appears to act as a free radical scavenger that extends useful fluorescence emission. Our original SlowFade® formulation (S2828) was designed to reduce the fading rate of fluorescein to almost zero (Figure 1); it does, however, decrease fluorescein's fluorescence intensity. Because it provides a nearly constant emission intensity, the original SlowFade® is especially useful for quantitative measurements, as well as applications that employ extreme, prolonged, or repeated excitation intensities such as confocal microscopy. This SlowFade® formulation can extend the useful life of fluorescein's fluorescence by more than 50-fold, and can preserve the signal in cell and tissue mounts for up to two years (Dr. Robert Bacallao, Northwestern University, personal

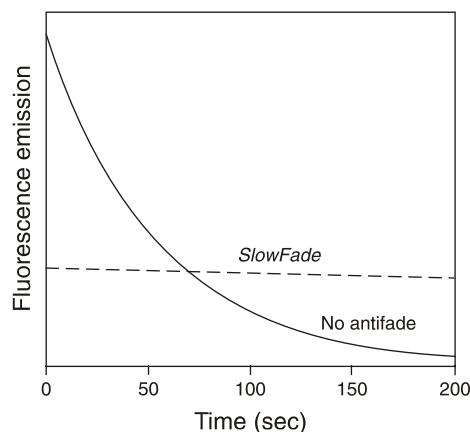


Figure 1. The fluorescence intensity of fluorescein as a function of illumination time, under the following conditions: in the presence of the SlowFade® antifade reagent ("SlowFade") and in the absence of an antifade reagent ("No antifade"). In these experiments, we added free fluorescein directly to a solution, and then examined the mixture in a capillary tube using a 20H / 0.4 lens. In real samples such as cells and tissues, we find that the local environment influences the bleaching rates, yielding results that are sometimes qualitatively different from those shown here.

communication). The practical limits will depend significantly on the particular dye and its surrounding environment. We must caution that the original SlowFade® formulation appears to excessively quench the fluorescence of 7-amino-4-methylcoumarin-3-acetic acid (AMCA) and Cascade Blue® fluorophores. We have also found that it is not generally effective as an antifade reagent for phycobiliproteins.

Our experiments show that SlowFade® provides better protection against fading when used at an appropriate pH. To maximize the performance of this product, we include a buffered wash solution with each kit.

For maximum resistance to photobleaching for a wide range of fluorescent dyes, we now offer SlowFade® Gold antifade reagent (S36936, S36937). SlowFade® Gold reagent is provided as a premixed and ready-to-use solution. Because SlowFade® Gold reagent does not cure over time, samples can be viewed immediately—simply tack the corners of the slide with hot wax or nail polish, then image. SlowFade® Gold reagent is intended for short-term use (3–4 weeks) only; samples mounted using SlowFade® Gold reagent may degrade over time. To save samples for months after mounting, we offer ProLong® Gold antifade reagent (P36930, P36934), which cures within 24 hours. ProLong® Gold and SlowFade® Gold antifade reagents are also available

with DAPI (P36931, P36935, S36938, S36939). Detailed protocols for *SlowFade*[®] Gold and ProLong[®] Gold reagents are available on our website at probes.invitrogen.com.

Materials

Kit Contents

- *SlowFade*[®] antifade reagent (Component A), 10 mL in 50% glycerol (v/v) ready-to-use and sufficient for at least 200 coverslip-size experiments
- **2X concentrated *SlowFade*[®] antifade reagent** (Component B), 2.5 mL provided for those applications in which glycerol may be incompatible
- **Equilibration buffer** (Component C), 60 mL

All components contain 2 mM sodium azide as a preservative.

Storage and Handling

Store the *SlowFade*[®] kit at room temperature. The components should be stable for at least one year.

Application

The *SlowFade*[®] Antifade Kit is for nonliving specimens only, including fixed cells, tissues, and cell-free preparations. A representative specimen should be tested for compatibility with the antifade reagent prior to research applications. Careful fixation is important for maintaining the sample and staining integrity. The effects of *SlowFade*[®] antifade reagent on binding affinities of dyes and ligands is currently undetermined.

This product is provided in two different solutions to allow its use in applications in which glycerol may be incompatible, such as mounting specimens containing lipophilic plasma membrane stains like DiI. The glycerol-containing solution is ready-to-use and recommended for most applications. The 2X concentrated solution can be diluted 1:1 in filtered distilled water to obtain a 1X glycerol-free solution. If glycerol is acceptable for the application, simply dilute the 2X concentrated solution with an equal volume of glycerol to generate additional glycerol-containing solution.

We have found that the *SlowFade*[®] formulation provides much better protection against fading when used at basic pH. Prior to adding *SlowFade*[®] antifade reagent, pre-equilibrate the specimen for five or ten minutes in the *SlowFade*[®] equilibration buffer provided. To speed the equilibration process, the researcher may wish to rinse the specimen three or four times in the equilibration buffer. However, please use the equilibration buffer sparingly; we have supplied only enough for a few rinses. Remove as much equilibration buffer as possible before adding the *SlowFade*[®] reagent to the sample so as not to dilute the antifade reagent. If you choose not to employ the equilibration buffer, the sample should be almost dry before applying the antifade reagent. After applying one drop of antifade reagent, mount and seal the slide. Sealed samples may turn yellow within about a month of preparation, even when stored in the refrigerator. This color does not appear to affect the fluorescence or morphology of the specimen.

To further reduce photobleaching, minimize the exposure of fluorescently labeled specimens to light with neutral density filters and expose samples only when observing or recording a signal. Maximize collection of fluorescence by using a minimum of optics, high-numerical aperture objectives, relatively low magnification, high-quality optical filters, and high-speed film or high-efficiency detectors.

General References

- “Guiding principles of specimen preservation for confocal fluorescence microscopy,” R. Bacallao *et al.*, in *Handbook of Biological Confocal Microscopy*, J. Pawley, Ed., pp 197–205, Plenum Press, New York (1990).
- “Photometric analysis of antifading reagents for immunofluorescence with laser and conventional illumination sources,” G. Bock *et al.*, *J Histochem Cytochem* 33, 699 (1985).
- “1,4-Diazobizyklo[2.2.2]oktan (DABCO) verzogest das Ausbleichen von immunofluoreszenzpreparaten,” G. Langanger, J. De Mey and H. Adam, *Mikroskopie* 40, 237 (1983).

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

Cat #	Product Name	Unit Size
P36930	ProLong [®] Gold antifade reagent.....	10 mL
P36934	ProLong [®] Gold antifade reagent *special packaging*	5 x 2 mL
P36931	ProLong [®] Gold antifade reagent with DAPI	10 mL
P36935	ProLong [®] Gold antifade reagent with DAPI *special packaging*	5 x 2 mL
S2828	<i>SlowFade</i> [®] Antifade Kit	1 kit
S36936	<i>SlowFade</i> [®] Gold antifade reagent	10 mL
S36937	<i>SlowFade</i> [®] Gold antifade reagent *special packaging*	5 X 2 mL
S36938	<i>SlowFade</i> [®] Gold antifade reagent with DAPI	10 mL
S36939	<i>SlowFade</i> [®] Gold antifade reagent with DAPI *special packaging*	5 X 2 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 Web site — 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.

Several Molecular Probes products and product applications are covered by U.S. and foreign patents and patents pending. Our products are not available for resale or other commercial uses without a specific agreement from Molecular Probes, Inc. We welcome inquiries about licensing the use of our dyes, trademarks or technologies. Please submit inquiries by e-mail to probesbusdev@invitrogen.com. 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.