

SlowFade® Gold Antifade Reagent

SlowFade® Gold Antifade Reagent with DAPI

Pub. no. MAN0003516

Revision A

Detailed protocol is available online at www.lifetechnologies.com/manuals.

Product	Amount	Cat. no.	Storage*
<i>SlowFade</i> ® Gold Antifade Reagent	1 × 2 mL	S36940	<ul style="list-style-type: none"> • Room temperature (15–30°C) • Protect from light
	5 × 2 mL	S36937	
	1 × 10 mL	S36936	
<i>SlowFade</i> ® Gold Antifade Reagent with DAPI	1 × 2 mL	S36942	
	5 × 2 mL	S36939	
	1 × 10 mL	S36938	

* Product may also be stored at $\leq -20^{\circ}\text{C}$.



1. Add 1 drop of *SlowFade*[®] Gold reagent onto glass slide.



2. Remove excess liquid from coverslip.



3. Place coverslip sample-side down onto *SlowFade*[®] Gold reagent on the glass slide.
Optional: Apply nail polish or hot wax to edges of coverslip to secure it.



4. Cure for 24 hours at room temperature in the dark.



5. Clean coverslip and image cells.

Limited Product Warranty

Life Technologies Corporation and/or its affiliate[s] warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

DISCLAIMER - LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE[S] DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE[S] BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

© 2013 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation and/or its affiliate[s] or their respective owners.

For support visit lifetechnologies.com/support
lifetechnologies.com