

SlowFade*[®] Diamond Antifade Mountant**SlowFade*[®] Diamond Antifade Mountant with DAPI**

Pub. no. MAN0010262

Revision B.0

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

Product	Amount	Cat. no.	Storage*
<i>SlowFade</i> [®] Diamond Antifade Mountant	5 × 2 mL	S36963	<ul style="list-style-type: none"> • 2–8°C • Protect from light
	10 mL	S36972	
<i>SlowFade</i> [®] Diamond Antifade Mountant with DAPI	5 × 2 mL	S36964	
	10 mL	S36973	

Refractive index of mountant: 1.42

* Product may also be stored at ≤−20°C. When stored as directed, the product is stable for at least 6 months.



1. Add 1 drop of *SlowFade*[®] Diamond Antifade Mountant onto glass slide.



2. Remove excess liquid from coverslip.



3. Place coverslip sample-side down onto *SlowFade*[®] Diamond mountant on the slide.

Optional: Apply nail polish or hot wax to edges of coverslip to secure it.



4. 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.

©2014 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. .

For support visit lifetechnologies.com/support
lifetechnologies.com