

ProLong[®] Live Antifade Reagent

Pub. no. MAN0011089

Revision A.0

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

Product	Amount	Cat. no.	Storage*
ProLong [®] Live Antifade Reagent	5 × 1 mL	P36974	• ≤ -20°C or 2–8°C
	1 mL	P36975	• Protect from light

* When stored at ≤ -20°C, the product is stable for at least 6 months with up to 4 freeze-thaw cycles.

When stored at 2–8°C, the product should be used within 30 days (a precipitate may form at this temperature).

1. Prepare imaging working solution by diluting the ProLong® Live Antifade Reagent 1:100 with Live Cell Imaging Solution (Cat. no. A14291DJ) or with FluoroBrite™ DMEM Media (Cat. no. A1896701).

Note: You may also use other suitable imaging media.

2. Stain the cells with the desired fluorescent dye and/or use cells expressing a desired fluorescent protein.
3. Wash the cells 1× with PBS.
4. Add imaging working solution (from step 1) to the cells.
5. Incubate cells in the dark for 15 minutes to 2 hours.
Note: For better results, 2 hour incubation is recommended.
6. Image or analyze the 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.

©2015 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