Live/Dead® Fixable Dead Cell Stain Kits

Pub. No. MAN0006891



Revision A.0

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

WARNING! For every chemical, read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from www.lifetechnologies.com/support.

Chemical	CAS	Warning
DMSO	67-68-5	Components of the product may be absorbed into the body through the skin.

- Thaw all reagents. Add 50 µL of Component B to 1 vial of Component A. Mix well. Use as soon as possible.
- 2. Centrifuge 1 \times 10 6 cells. Wash and resuspend the cells in 1 mL PBS.
- Add 1 µL of the dissolved dye (from step 1) to the cells and mix well. Incubate for 30 minutes, protected from light.
- 4. Optional: Wash and fix the cells with 100 μL of 37% formaldehyde for 15 minutes.



For Research Use Only. Not for use in diagnostic procedures.

Excitation Source and Detection Channel

Reactive dye	Excitation	Detection
Blue fluorescent	UV	~450 nm
Violet fluorescent	405 nm	~450 nm
Aqua fluorescent	405 nm	~525 nm
Yellow fluorescent	405 nm	~575 nm
Green fluorescent	488 nm	~530 nm
Red fluorescent	488 nm	~ 585 nm
Far red fluorescent	633/635 nm	~660 nm
Near-IR fluorescent	633/635 nm	~780 nm

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.

DBCLAMER - LIFE TECHNOLOBES COMPORATION AND/OR ITS AFFLIATESI DBCLAMA ALL WARRANTES WITH RESPECT TO THIS DOCUMENT, DEVERSED OR MARLED, INCLUDING UN TO LIMITED TO THOSE OF MERCINARUM FILTERISTS OF A PARTICULAR PURPOSE, OR NON-INTRODUCED. IN THE ORTEN ALL/WED FLAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR TS AFFLIATESI BE LUBLE, WIETHER IN CONTACT, TORY, WARRANTY ON UNCE ARISING FIRM. THIS DOCUMENT, INCLUDIES TECHNOLOGIES AND/OR TS AFFLIATESI BE LUBLE, WIETHER IN CONTACT, TORY, WARRANTY ON UNCE ARISING FIRM. THIS DOCUMENT, INCLUDIES DUT ON LIFE DITE DITE TECHNOLOGIES AND/OR TO A PARTICULAR PURPOSE. INCLUDIES AND/OR TO A PARTICULAR PURPOSE, ON ON-INTRODUCTION UNCER ARISING FIRM. THIS DOCUMENT, INCLUDIES DUT ON LIFE DITE DITE DITE TECHNOLOGIES AND/OR TO A PARTICULAR PURPOSE.

@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

8 October 2014