

FxCycle™ PI/RNase Staining Solution

Pub. no. MAN0007554

Rev. Date 26 November 2012

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



1. Begin with 1×10^6 cells that have been ethanol fixed. Wash cells once and remove the buffer from the cell pellet.



2. Add 0.5 mL of FxCycle™ PI/RNase Solution to the cell pellet.



3. Vortex gently to resuspend the cell pellet.



4. Incubate for 15–30 minutes at room temperature in the dark.



5. Analyze using 488-nm, 532-nm, or similar excitation, and collect using 585/42 bandpass filter or equivalent.

For research use only. Not for use in diagnostic procedures.

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.

© 2012 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.

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.

For support visit www.lifetechnologies.com/support
www.lifetechnologies.com