

LIVE/DEAD™ Cell Imaging Kit

Catalog Number R37601

Pub. no. MAN0006090 Rev. A.0

Contents

Component	Amount
Live Green (Comp. A)	10 × 1 mL (1 µM solution)
Dead Red (Comp. B)	10 × 1 µL (lyophilized stock)



1. Thaw vials.



2. Transfer Live Green (Comp. A) into Dead Red (Comp. B).



3. Mix to create 2X working solution.



4. Add equal volume of 2X working solution to cells within 2 hours.



5. Incubate for 15 minutes at 20–25°C.



6. Image cells.

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

The information in this guide is subject to change without notice.

DISCLAIMER

TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Corporate entity: Life Technologies | Carlsbad, CA 92008 USA | Toll Free in USA 1.800.955.6288

©2016 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 thermofisher.com/support
thermofisher.com

ThermoFisher
 S C I E N T I F I C