

CellTrace™ Yellow Cell Proliferation Kit

Cat. No. C34567, C34573

Pub. No. MAN0014056

Revision D

Detailed protocol is available online at www.thermofisher.com.

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.thermofisher.com/support.

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



1. Remove supernatant from the cell pellet.



2. Add CellTrace™ Yellow (1:500 dilution) staining solution and gently re-suspend the cells.



3. Incubate at 37°C for 20 minutes.

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

ThermoFisher
SCIENTIFIC



4. Add complete culture medium and mix.



5. Incubate at 37°C for 5 minutes.



6. Pellet the cells and remove the supernatant.



7. Resuspend the cell pellet in fresh, pre-warmed complete culture medium.



8. Analyze by flow cytometry or proceed with stimulation or incubation.

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.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses.

Manufacturer: Life Technologies Corporation | 29851 Willow Creek Road | Eugene, OR 97402

© 2017 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
SCIENTIFIC