

CellROX® Oxidative Stress Reagents

Cat. Nos. C10422, C10443, C10444, C10448

Pub. No. MAN0007457

Revision 1.0

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.

Spectral Characteristics of CellROX® Oxidative Stress Reagents

CellROX® Reagent	Ex/Em maxima (in nm)
CellROX® Deep Red Reagent	640/665
CellROX® Orange Reagent	545/565
CellROX® Green Reagent	485/520

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

Quick Protocol

1. Treat the cells with the test compound or drug.
2. Add the CellROX® Reagent at a final concentration of 5 µM.
3. Incubate the cells for 30 minutes at 37°C.
4. Remove medium and wash the cells 3 times with PBS.
5. *Optional:* If using CellROX® Deep Red or CellROX® Green, you may fix the cells with 3.7% formaldehyde for 15 minutes.
6. *Optional:* You may stain the cells with NucBlue™ Live Cell Stain, a nuclear counterstain, or another counterstain at this time.
7. *Optional:* If using CellROX® Green, you may permeabilize the cells with 0.5% Triton® X-100 for 10 minutes, if multiplexing with another reagent is desired.
8. 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.

© 2013 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. Triton is a registered trademark of Dow Chemical Co.

For support visit lifetechnologies.com/support
lifetechnologies.com