

# CellTracker™ Fluorescent Probes

Pub. no. MAN0010620

Revision A.0

Detailed protocol is available online at [www.lifetechnologies.com/manuals](http://www.lifetechnologies.com/manuals).

Product	Amount	Storage
CellTracker™ Blue and CellTracker™ Green BODIPY® dyes	5 mg	<ul style="list-style-type: none"> <li>• ≤ -20°C</li> <li>• Desiccate</li> <li>• Protect from light</li> </ul>
CellTracker™ Violet BMQC dye	5 × 0.1 mg	
CellTracker™ Green CMFDA and CellTracker™ Orange CMTMR dyes	1 mg	
CellTracker™ Green CMFDA, CellTracker™ Orange CMRA, and CellTracker™ Red CMPTX dyes	20 × 50 µg	
CellTracker™ Deep Red dye	20 × 15 µg	

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

## Quick Protocol



1. Remove culture media from cells.



2. Add CellTracker™ (1:1000 dilution) staining solution.



3. Incubate at 37°C for 30 minutes.



4. Remove staining solution.



5. Add imaging media.



6. Image.

### 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](http://www.lifetechnologies.com/termsandconditions). If you have any questions, please contact Life Technologies at [www.lifetechnologies.com/support](http://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.

© 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](http://lifetechnologies.com/support)  
[lifetechnologies.com](http://lifetechnologies.com)