

pHrodo™ Red and Green Dextran, 10,000 MW

Revision Date 27 June 2012

Catalog Nos. P10361, P35368

MAN0007046

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

BioParticles®	Ex/Em (nm)	Cat. no.
pHrodo™ Red Dextran, 10,000 MW	560/585	P10361
pHrodo™ Green Dextran, 10,000 MW	509/533	P35368

Quick Protocol

1. Resuspend pHrodo™ Dextran in 0.5 mL deionized water for a 1 mg/mL stock solution. You may store unused pHrodo™ Dextran stock solutions at -20°C for up to six weeks.

For research use only. Not for human or animal therapeutic or diagnostic use.

2. Dilute the stock solution in Live Cell Imaging Solution (LCIS) or the preferred buffer at pH 7.4 for a working solution at a final concentration of 50 to 100 $\mu\text{g}/\text{mL}$.
3. *Optional:* Pretreat cells with effector compound (e.g., 100 μM Dynasore to inhibit endocytosis).
4. Remove growth medium from the cells and wash 1X with LCIS. Replace with working pHrodo™ Dextran solution and incubate at 37°C for 10–30 minutes.
5. Analyze the cells using the appropriate Ex/Em maxima.

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 or their respective owners. 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