QUICK REFERENCE

Premo[™] Autophagy Sensors



Cat. nos. P36239, P36240, P36241 Pub. no. MAN0007579 Rev. 1.0
Complete detailed protocol is available online at www.lifetechnologies.com.



 Add 2 µL of Premo[™] Autophagy Sensor (Component A) per 10,000 cells in media.



Note: Adjust Premo[™] Autophagy Sensor amount for the desired expression level.



2. Incubate for 16 hours at 37°C.



3. Image cells.



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

Guidelines for Using Premo[™] Autophagy Sensors

- Premo[™] Autophagy Sensors work with most cell types between 10 and 50 particles per cell (PPC).
- For best results, transduce cells at a confluence of no more than 70%.
- For optimal results, you may need to alter the PPC, volume, cell density, temperature, or incubation time.

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 and/or its affiliate(s) or their respective owners.

DISCLAIMER - LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTES 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, PUNTITE, MULTIPE CR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOR.

For support visit lifetechnologies.com/support

lifetechnologies.com

4 December 2012