

pHrodo™ Red and Green BioParticles® Conjugates

Revision 2.0

MAN0007045

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

BioParticles® Conjugates for Phagocytosis	Ex/Em (nm)	Cat. no.
pHrodo™ Red <i>E. coli</i> BioParticles® Conjugate	560/585	P35361
pHrodo™ Green <i>E. coli</i> BioParticles® Conjugate	509/533	P35366
pHrodo™ Red <i>S. aureus</i> BioParticles® Conjugate	560/585	A10010
pHrodo™ Green <i>S. aureus</i> BioParticles® Conjugate	509/533	P35367
pHrodo™ Red Zymosan A Bioparticles® Conjugate	560/585	P35364
pHrodo™ Green Zymosan A Bioparticles® Conjugate	509/533	P35365

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

1. Resuspend pHrodo™ *E. coli* or *S. aureus* BioParticles® conjugates to 1 mg/mL or Zymosan BioParticles® conjugates to 0.5 mg/mL in Live Cell Imaging Solution or the preferred buffer at pH 7.4. Opsonize, if desired.
2. Vortex well or sonicate to homogeneously disperse the particles.
3. *Optional:* Pretreat cells with effector compound (e.g., 10 μM cytochalasin D to inhibit phagocytosis).
4. Remove the medium from the phagocytic cells and replace it with BioParticles® suspension. Incubate at 37°C for 60–90 minutes for maximal uptake.
5. Analyze the cells using the appropriate Ex/Em maxima. You may store unused BioParticles® suspension for 3–5 days at 4°C.
Do not freeze resuspended BioParticles® conjugates.

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. .

For support visit lifetechnologies.com/support
lifetechnologies.com