


pHrodo™ Red, Deep Red, and Green BioParticles™ conjugates for phagocytosis

Catalog Numbers P35361, P35366, P35364, P35365, A10010, P35367, and P35360

Doc. Part No. mp35361 Pub. No. MAN0002447 Rev. A.0

 **WARNING!** 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 thermofisher.com/support.

Contents and storage

Material	Cat. no.	Amount	Ex/Em Maxima™	Storage ^[1]
pHrodo™ Red <i>E. coli</i> BioParticles™ Conjugate	P35361	5 vials each containing 2 mg lyophilized product	560/585 nm	<ul style="list-style-type: none"> • ≤-20°C • Desiccate • Protect from light
pHrodo™ Deep Red <i>E. coli</i> BioParticles™ Conjugate	P35360		640/655 nm	
pHrodo™ Green <i>E. coli</i> BioParticles™ Conjugate	P35366		509/533 nm	
pHrodo™ Red Zymosan A BioParticles™ Conjugate	P35364	5 vials each containing 1 mg lyophilized product	560/585 nm	
pHrodo™ Green Zymosan A BioParticles™ Conjugate	P35365		509/533 nm	
pHrodo™ Red <i>S. aureus</i> BioParticles™ Conjugate	A10010	5 vials each containing 2 mg lyophilized product	560/585 nm	
pHrodo™ Green <i>S. aureus</i> BioParticles™ Conjugate	P35367		509/533 nm	
Number of assays: Sufficient for 100 assays when using described protocol.				

^[1] When stored as directed, the product is stable for at least 6 months.

Product description

pHrodo™ BioParticles™ Conjugates are novel, no-wash fluorogenic reagents developed for quantitative measurements of phagocytosis and its regulation by drugs, genetic, or environmental factors. With an optional no-cell background subtraction method, a large and specific signal is obtained from cells that ingest the particles, providing a specific index of phagocytosis with a variety of pretreatments or conditions. The unique pHrodo™-based system measures phagocytic activity based on acidification of the particles as they are ingested, eliminating the wash and quenching steps that are necessary with nonfluorogenic indicators of bacterial uptake. To achieve this, the particles are conjugated to pHrodo™ dye, a fluorogenic dye that dramatically increases in fluorescence as the pH of its surroundings becomes more acidic (Figure 1). We have included sufficient pHrodo™ BioParticles™ Conjugate for ~100 or 1000 wells (depending on application) in a 96-well format, with step-by-step instructions for performing this assay for both traditional imaging as well as in a fluorescence microplate reader. With proper changes to the protocol, other plate formats can easily be adapted to platforms such as high-content screening (HCS), high-throughput screening (HTS), fluorescent microplate reader, and flow cytometry. With proper settings, these reagents can also be adapted for benchtop instruments, such as EVOS Cell Imaging System, CellInsight CX7 or CX5, and Attune NXT Flow Cytometer.

The methodology for this reagent's use was developed using adherent RAW and MMM (J774A.1) murine macrophage cells, but can be adapted for use with other adherent cells, primary cells, or even cells in suspension. Cells assayed for phagocytic activity with pHrodo™ BioParticles™ conjugates may also be fixed with standard 2–4% paraformaldehyde solutions for later analysis, preserving differences in signal between control and experimental samples with high fidelity for up to 48 hours. pHrodo™ BioParticles™ conjugate preparations are also amenable to opsonization (Cat. Nos. E2870, S2860), which can greatly enhance their uptake and signal strength in the assay.

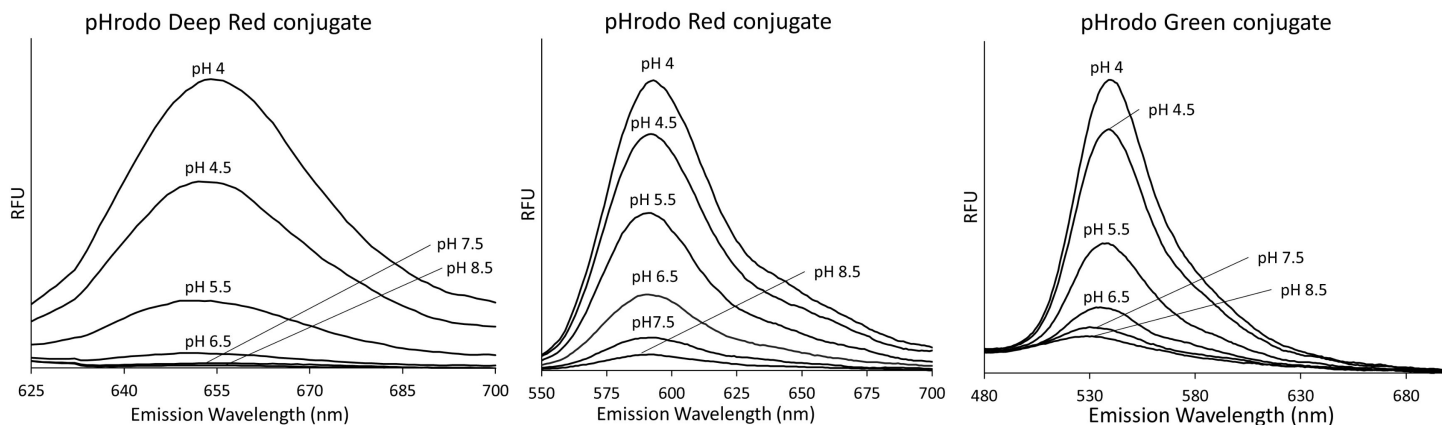


Figure 1 The fluorescence emission spectra of pHrodo™ Deep Red , Red, and Green conjugates

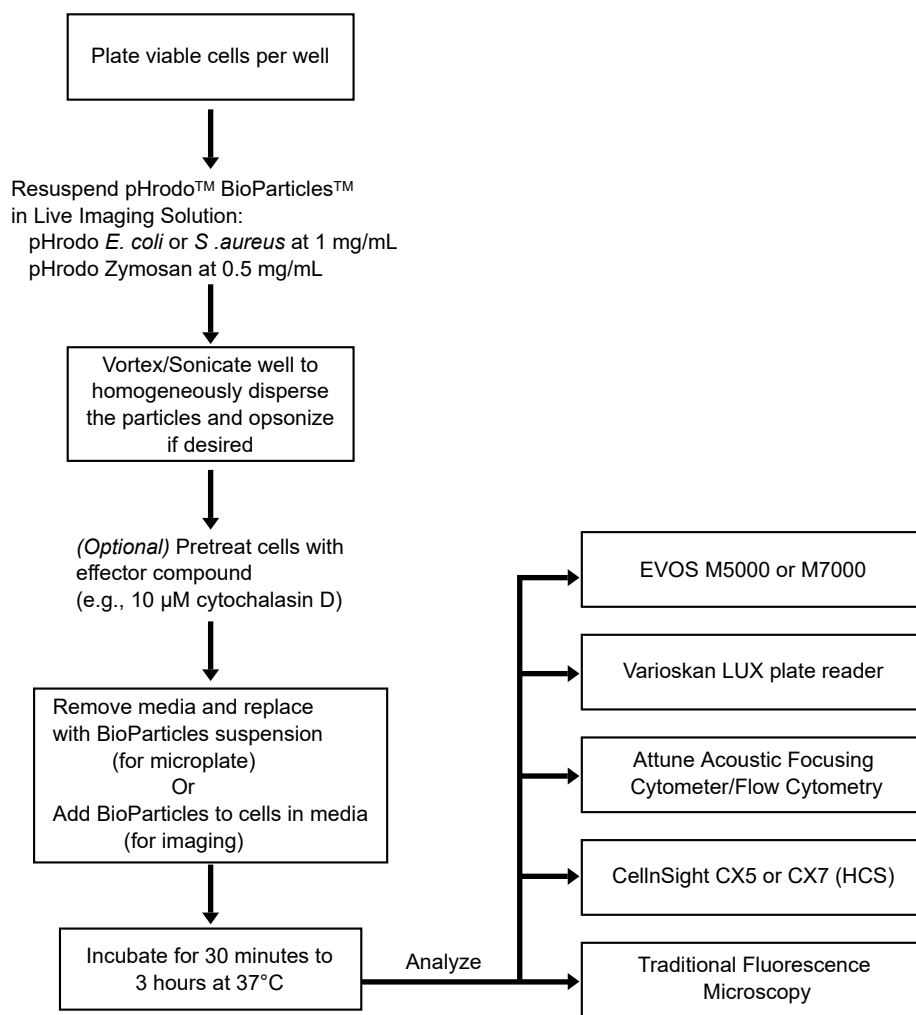


Figure 2 Workflow for pHrodo™ BioParticles™ conjugates

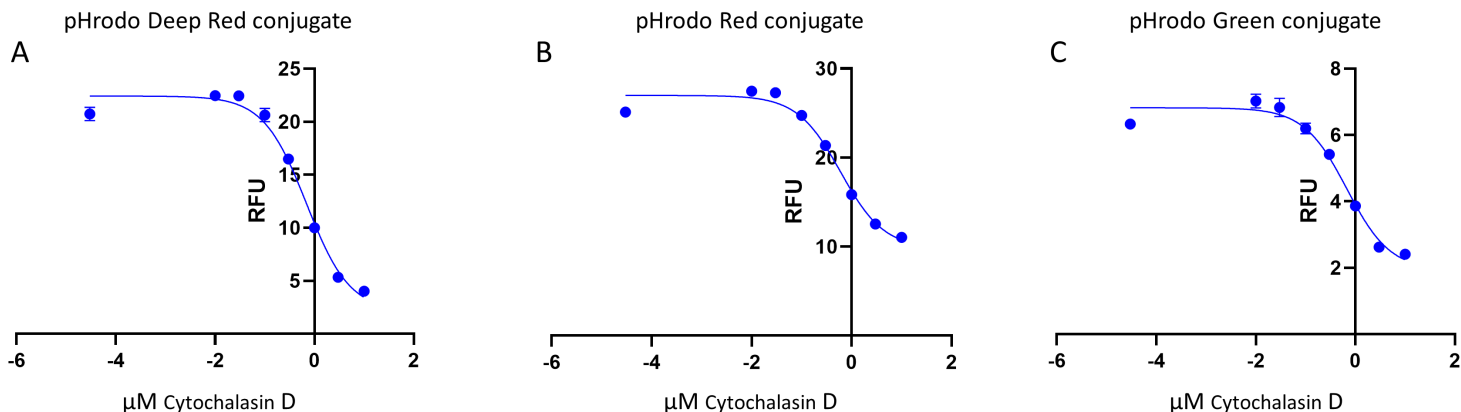


Figure 3 Effect of different concentrations of Cytochalasin D in RAW cells detected with (A) pHrodo™ Deep Red *E. coli* BioParticles™, (B) pHrodo™ Red *E. coli* BioParticles™, (C) pHrodo™ Green *E. coli* BioParticles™ conjugates in RAW cells while using a fluorescent microplate reader (shown here without the background subtraction).

Required materials not provided

- Macrophage RAW and MMM cells cultured in cell growth medium. Other cell lines may also be used, if preferred.
- Uptake Buffer; we recommend using Live Cell Imaging Solution (Cat. No. A14291DJ) for best results. Alternatively, you may use any other appropriate buffer at pH 7.4.
- 96-well microplate or any other plate, with proper installment capable of detecting fluorescence according to emission/excitation maxima given in Table 1.
- Bath Sonicator
- (Optional) Opti-MEM™ culture medium (Cat. No. 31985-062).
- (Optional) Stock solutions of experimental effector compounds for testing their influence on phagocytosis (for example, cytochalasin D inhibits phagocytosis by inhibiting actin cytoskeletal rearrangements).

Experimental protocol for imaging

The following protocol describes an experimental test of phagocytic function with appropriate controls in a 96-well plate; however, other plate formats can be adopted with proper adjustments to the protocol. This protocol describes the use of one vial of particles, prepared at 1 mg/mL in the buffered saline solution of your choice then added to cells in complete media at 1:10 for a final concentration of 0.1 mg/mg in the imaging assay. To minimize background fluorescence from non-ingested pHrodo™ BioParticles™ Conjugate, we strongly recommend controlling the extracellular pH by adding Live Cell Imaging Solution (Cat. No. A14291DJ) for best results. Alternatively, you may use any other appropriate buffer at pH 7.4

Procedural guidelines

- Work with cells in aseptic conditions to prevent microbial contamination
 - Warm buffer and pHrodo bioparticles™ to 20–25°C
- Note:** The kit will use 2 mL of buffer per vial to generate a 10X solution of particles that is enough for 1,000 tests at 100 $\mu\text{g}/\text{mL}$ working concentration.

- (Optional) Remove cell culture medium, wash once and replace with pH neutral saline solution warmed to 37°C. Ensure it contains 5–10 mM glucose to support metabolism during the assay.

Note: This step will remove the cell culture medium, which can be a source of autofluorescence in the assay. If the assay is done in saline solution, be sure that the 37°C incubation step happens in a benchtop incubator, and not in the 5% CO₂ atmospheric conditions of a cell culture incubator required for carrying out the experiment in cell culture media.

Preparing the cells

1. Subculture the RAW or MMM macrophage cells (or preferred cell type) in complete medium for 3–4 days in advance of performing the assay.
2. On the day of the assay, harvest the cells from the culture plates and centrifuge the suspension. Resuspend the pellet in FluoroBrite™ DMEM (Cat. No. A1896701) medium or preferred culture medium at 10⁶ cells/mL. Scale your culture to aim for $\sim 2 \times 10^6$ cells per vial of pHrodo™ BioParticles™ Conjugate. Alternatively, cells can be plated into the 96-well plate a day or more in advance, with the aim of having 10,000 viable cells per well on the day of the assay.
3. Plate the cells into a 96-well plate at 10,000 cells/well using 100 μL per well. We recommend plating your positive control and experimental wells in triplicate or greater. Be sure to leave one well empty of cells for every positive control well, so that a no-cell control background subtraction may be performed.

If you are using cells other than RAW or MMM, you may need to determine optimal cell culture conditions and densities for your specific cell type. In general, better signals in the plate reader are obtained with maximal cell densities.

For example, plate four columns of four wells, leaving the fifth column of four empty for the no-cell control. Note that higher background fluorescence levels may be seen with acidic poly-D-lysine coated microplates.

4. Add 100 μL of FluoroBrite™ DMEM (Cat. No. A1896701) medium or complete culture medium to the wells left aside for the no-cell background determination.
5. Cover the loaded microplate and allow the cells to settle and adhere to the microplate for at least one hour in a humidified incubator with 5% CO_2 at 37°C.
6. Prepare the experimental wells by adding the experimental phagocytosis effector (e.g., Cytochalasin D) at the desired concentrations, taking care to add vehicle controls to untreated wells. Note that the time and concentration of experimental effector pretreatment may vary greatly with the agent or treatment under study.

Preparing the BioParticles™ conjugate

1. Resuspend the pHrodo BioParticles™ in pH neutral saline solution by adding 2 mL of buffered saline solution to a vial to generate a 1 mg/mL stock suspension or 0.5 mg/mL for pHrodo™ Zymosan A BioParticles™.

2. Triturate the particles well to ensure complete hydration.

Note: For best results, place the vial of resuspended particles in a bath sonicator for 10 minutes.

Aggregation of the particles immediately after resuspension is normal and monodispersion of particles will aid in phagocytic uptake and subsequent microscopic examination. If particle aggregates are still evident after 10 minutes, repeat the 10 minute sonication step.

3. (Optional) Opsonization of the bioparticles may be desired. If so, pretreat the particles with opsonizing reagent after sonication as detailed in Cat No. E2870.

Adding the Fluorescent Particles

1. Add the suspension of particles to the cell culture microplate in a 1:10 dilution, or 10 μL of particles added to 100 μL of cell culture medium for a particle concentration of 100 $\mu\text{g}/\text{mL}$.
2. Mix well to ensure complete dispersion of the particles into the preparation.

Typical imaging and high content analysis protocols will vary widely based on cell type and density. Most will achieve the best signal with particles present between 50 and 200 $\mu\text{g}/\text{mL}$.

3. Place the cells at 37°C for 30 minutes to 3 hours.

This will vary with the cell type in use and the cellular activity, but fluorescence from internalized particles should become visible after 30 to 60 minutes, when the pH of the phagosome begins to drop below 6. The experiment can be left to run as long as necessary to achieve signal maximum. Typical phagocytosis experiments are left to run for two hours.

Fluorescence Measurements and Results

- Image all experimental, control, and no-cell control wells using the appropriate imaging conditions according to the excitation and emission maxima according to the table in “Contents and storage” on page 1.

- To visualize signal onset in a time lapse experiment, it is best to generate a set of data at the endpoint to determine the optimal exposure time and illumination intensity that will be used in a follow up study to monitor internalization and signal generation from the particles. Typical exposure conditions should keep the signal between 25% and 30% of maximum RFUs for the system to avoid overexposure.

Experimental Protocol for Microplates

Phagocytosis assay protocol

The following protocol describes an experimental test of phagocytic function with appropriate controls in a 96-well plate; however, other plate formats can be adopted with proper adjustments to the protocol. Cellular auto fluorescence background is determined with cells plated free of pHrodo™ BioParticles™ Conjugates (but otherwise under control and experimental conditions), and reagent background fluorescence is determined using wells that contain the pHrodo™ BioParticles™ Conjugates but no cells. This protocol describes the use of one vial of particles, prepared at 1 mg/mL in the buffered saline solution of your choice. To minimize background fluorescence from non-ingested pHrodo™ BioParticles™ Conjugate, we strongly recommend controlling the extracellular pH by adding Live Cell Imaging Solution (Cat. no. A14291DJ) for best results. Alternatively, you may use any other appropriate buffer at pH 7.4.

Assay controls

To minimize experimental errors, we recommend making measurements from a minimum of three replicates of positive control, experimental, and no-cell control samples, though the numbers of experimental and control wells can be adjusted as required to meet the needs of the particular study.

Amount of BioParticles™

A single vial of pHrodo™ BioParticles™ conjugate dilutes to 2 mL of volume that will be used in the assay, which is distributed across 20 wells. 100 μL of this suspension is used per sample well, including no-cell background controls. The average fluorescence value of these no-cell background control wells is subtracted from all cell-containing wells at the end of the assay to yield a cell-specific, net phagocytosis signal. Note that a large specific signal can readily be obtained with or without background subtraction, as shown in Figure 3.

Preparing the cells

1. Subculture the RAW or MMM macrophage cells (or preferred cell type) in complete medium for 3–4 days in advance of performing the assay.
2. On the day of the assay, harvest the cells from the culture plates and centrifuge the suspension. Resuspend the pellet in Opti-MEM™ medium or preferred culture medium at 10^6 cells/mL. Scale your culture to aim for $\sim 2 \times 10^6$ cells per vial of pHrodo™ BioParticles™ Conjugate. Alternatively, cells can be plated into the 96-well plate a day or more in advance, with the aim of having 100,000 viable cells per well on the day of the assay.

If you are using cells other than RAW or MMM, you may need to determine optimal cell culture conditions and densities for your specific cell type. In general, better signals in the plate reader are obtained with maximal cell densities.

3. Plate the cells into a 96-well plate at 100,000 cells/well using 100 μ L per well. We recommend plating your positive control and experimental wells in triplicate or greater. Be sure to leave one well empty of cells for every positive control well, so that a no-cell control background subtraction may be performed.

For example, plate four columns of four wells, leaving the fifth column of four empty for the no-cell control. Note that higher background fluorescence levels may be seen with acidic poly-D-lysine coated microplates.

4. Add 100 μ L of Opti-MEM™ medium or complete culture medium to the wells left aside for the no-cell background determination.
5. Cover the loaded microplate and allow the cells to settle and adhere to the microplate for at least one hour in a humidified incubator with 5% CO₂ at 37°C.
6. Prepare the experimental wells by adding the experimental phagocytosis effector (e.g., Cytochalasin D) at the desired concentrations, taking care to add vehicle controls to untreated wells. Note that the time and concentration of experimental effector pretreatment may vary greatly with the agent or treatment under study.

Preparing the BioParticles™ conjugate

1. Thaw one vial each of the pHrodo™ BioParticles™ fluorescent particles for every 20 wells to be tested. This number includes the no-cell control wells that will receive fluorescent particles, but no cells. Pipette 2 mL Uptake Buffer such as the Live Cell Imaging Solution (Cat. No. A14291DJ) into the vial containing lyophilized product and briefly vortex the solution to completely resuspend the particles so that you have 1 mg/mL for pHrodo™ *E. coli* or *S. aureus* BioParticles™, or 0.5 mg/mL for pHrodo™ Zymosan A BioParticles™.
2. Transfer the suspension into a clean glass tube and sonicate for 10 minutes, until all the fluorescent particles are homogeneously dispersed.

Adding the fluorescent particles

1. After the cells have adhered and the phagocytosis effectors have been added, remove the culture medium from each of the microplate wells by vacuum aspiration.
2. Quickly replace the culture medium with 100 μ L of the prepared pHrodo™ BioParticles™ suspension from step 2.2 on page 5, adding it to the positive control, experimental, and no-cell background subtraction wells. Experimental effector solutions may be prepared ahead of time with separate vials of pHrodo™ BioParticles™ suspension to keep them present throughout the assay.
3. Cover and transfer the microplate to an incubator warmed to 37°C for 1–2 hours to allow phagocytosis and acidification to reach its maximum. Although aseptic techniques have been used to produce pHrodo™ BioParticles™ conjugates, these products are not sterile and incubation of more than 4 hours is not recommended. Do not use an elevated CO₂ cell culture incubator unless the Uptake Buffer in use has a bicarbonate buffering system, because elevated CO₂ levels will artificially acidify the buffer and elevate the background fluorescence.

Fluorescence measurements and results

1. Scan or image all experimental, control, and no-cell control wells of the microplate in the fluorescence plate reader or any other appropriate fluorescent instrument using the appropriate settings according to the excitation and emission maxima given in Table 1.
2. If using a fluorescence plate reader, calculate the net phagocytosis and the response to the phagocytosis effector agent. Net phagocytosis is calculated by subtracting the average fluorescence intensity of the no-cell negative-control wells from all positive control and experimental wells. The mean and standard deviation of the net positive control and net experimental wells should then be calculated. The phagocytosis response to the experimental effector (% Effect) can then be calculated as a fraction of the net positive control phagocytosis as follows:

$$\% \text{ Effect} = \frac{\text{Net experimental phagocytosis} \times 100\%}{\text{Net positive control phagocytosis}}$$

Product list

Current prices may be obtained from our website or from our Customer Service Department.

Cat #	Product Name	Unit Size
A10010	pHrodo™ Red <i>S. aureus</i> BioParticles™ conjugate for phagocytosis	5 × 2 mg
P35361	pHrodo™ Red <i>E. coli</i> BioParticles™ conjugate for phagocytosis	5 × 2 mg
P35364	pHrodo™ Red Zymosan A BioParticles™ conjugate *for phagocytosis*	5 × 1 mg
P35365	pHrodo™ Green Zymosan A BioParticles™ conjugate *for phagocytosis*	5 × 1 mg
P35366	pHrodo™ Green <i>E. coli</i> BioParticles™ conjugate *for phagocytosis*	5 × 2 mg
P35367	pHrodo™ Green <i>S. aureus</i> BioParticles™ conjugate *for phagocytosis*	5 × 2 mg
P35360	pHrodo™ Deep Red <i>E. coli</i> BioParticles™ conjugate for phagocytosis	5 × 2 mg
Related Products		
P10361	pHrodo™ Red dextran, 10,000 MW *for endocytosis*	0.5 mg
P35368	pHrodo™ Green dextran, 10,000 MW *for endocytosis*	0.5 mg
P36600	pHrodo™ Red, succinimidyl ester (pHrodo™ Red, SE)	1 mg
P35370	pHrodo™ Green C ₂ -maleimide	1 mg
P35371	pHrodo™ Red C ₂ -maleimide	1 mg
A10025	pHrodo™ Red <i>E. coli</i> BioParticles™ Phagocytosis Kit *for flow cytometry* *100 tests*	1 kit
A10026	pHrodo™ Red Phagocytosis Particle Labeling Kit *for flow cytometry* *100 tests*	1 kit
P35362	pHrodo™ Red Avidin *Fluorogenic pH sensor*	1 mg
P35363	pHrodo™ Red Microscale Labeling Kit *Fluorogenic pH sensor* *3 labelings*	1 kit
A14291DJ	Live Cell Imaging Solution	500 mL
R37602	Image-iT™ Fixation/Permeabilization Kit	1 kit
R37603	BackDrop™ Background Suppressor *for live cells*	1 kit
R37605	NucBlue™ Live Cell Stain *Hoechst 33342 special formulation*	1 kit
R37606	NucBlue™ Fixed Cell Stain *DAPI special formulation*	1 kit
E2870	<i>Escherichia coli</i> BioParticles™ opsonizing reagent	1 unit
S2860	<i>Staphylococcus aureus</i> BioParticles™ opsonizing reagent	1 unit

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 at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Life Technologies Corporation | 29851 Willow Creek | Eugene, OR 97402
For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Revision history: Pub. No. MAN0002447

Revision	Date	Description
5.0	29 October 2020	Rebranding and addition of imaging protocol
4.0	30 May 2013	Baseline for revision

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

