

pHrodo[®] Bioparticles[®] for Incucyte[®] Phagocytosis Assays

Product Information

Presentation, Storage and Stability

pHrodo[®] Bioparticles[®] for Incucyte[®] are supplied as lyophilized solid in sufficient quantity capable of performing 100-200 tests (1 test = 1 well of 96-well microtiter plate). The lyophilized solid should be stored at -20°C and once solubilized the suspension should be stored at +4° C. When stored as described, the lyophilized solid will be stable for at least 6 months and the suspension for at least 1 month.

Background and Intended Use

pHrodo[®] Bioparticles[®] for Incucyte[®] are sterile fluorogenic reagents ideally suited to a simple mix-and-read, real-time live-cell quantification of phagocytosis. The unique pHrodo[®]-based system exploits the acidic environment of the phagosome to quantify phagocytosis. As pHrodo[®] Bioparticles[®] for Incucyte[®] residing in the neutral extracellular solution (pH 7.4) are engulfed by phagocytes and enter the acidic phagosome (pH 4.5–5.5), a substantial increase in fluorescence is observed. Application of pHrodo[®] Bioparticles[®] for Incucyte[®] to non-phagocytic cells yields little or no fluorescent signal. With the Incucyte[®] integrated analysis software, background fluorescence is

minimized. These fully sterilized reagents have been validated for use with the Incucyte[®] Live-Cell Analysis System and enable real-time evaluation of phagocytic regulation by pharmacological agents as well as genetic and environmental factors.

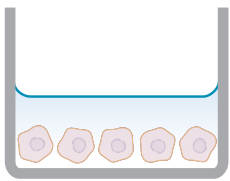
Recommended Use

We recommend that pHrodo[®] Bioparticles[®] for Incucyte[®] are prepared at a stock concentration of 1 mg per mL in full media or PBS. The Bioparticles[®] may then be diluted for direct addition to cells seeded in a 96-well plate to yield 10 µg per well (for *E. coli* and *S. aureus*) or 5 µg per well (for Zymosan). When used in an Incucyte[®] Live-Cell Analysis System, we recommend data collection every 15 minutes.

Product Name	Cat. No.	Amount	Ex. Maxima	Em. Maxima
pHrodo [®] Red <i>E. coli</i> Bioparticles [®] for Incucyte [®]	4615	2 mg	560 nm	585 nm
pHrodo [®] Green <i>E. coli</i> Bioparticles [®] for Incucyte [®]	4616	2 mg	509 nm	533 nm
pHrodo [®] Red Zymosan Bioparticles [®] for Incucyte [®]	4617	1 mg	560 nm	585 nm
pHrodo [®] Green Zymosan Bioparticles [®] for Incucyte [®]	4618	1 mg	509 nm	533 nm
pHrodo [®] Red <i>S. aureus</i> Bioparticles [®] for Incucyte [®]	4619	2 mg	560 nm	585 nm
pHrodo [®] Green <i>S. aureus</i> Bioparticles [®] for Incucyte [®]	4620	2 mg	509 nm	533 nm

Quick Guide

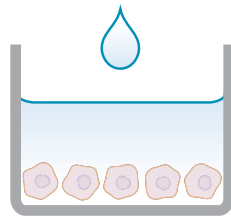
1. Seed target cells



Phagocyte Cell Seeding

Seed phagocytes (50 μ L/well, 1×10^3 to 1×10^4 cells/well) into the 96-well plate and leave to adhere (2–16 h).

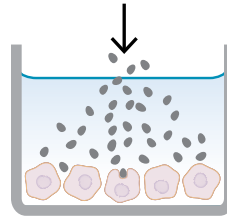
2. Treat cells



Activator|Inhibitor or Molecular Intervention

Add the desired treatments (25 μ L/well) at 4X final assay concentrations.

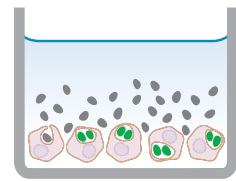
3. Add pHrodo[®] Bioparticles[®] for Incucyte[®]



pHrodo[®] Bioparticles[®] Addition

Add your choice of Bioparticle[®] (e.g., *E. coli*, *S. aureus*, Zymosan) to the 96-well plate (approximately 10 μ g per well depending on Bioparticle[®] 25 μ L/well at 4X final assay concentrations).

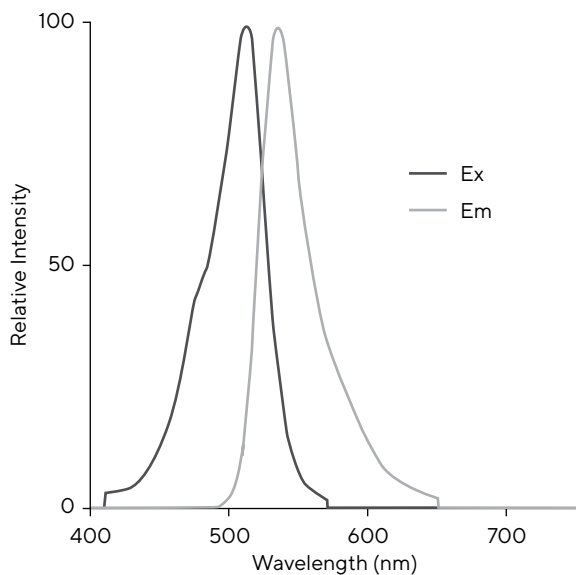
4. Live-cell fluorescent imaging



Automated Imaging and Quantitative Analysis

Capture images every 10–30 minutes (20X or 10X) in Incucyte[®] Live-Cell Analysis System for 2–48 hours. Analyze using integrated software.

A. pHrodo[®] Green



B. pHrodo[®] Red

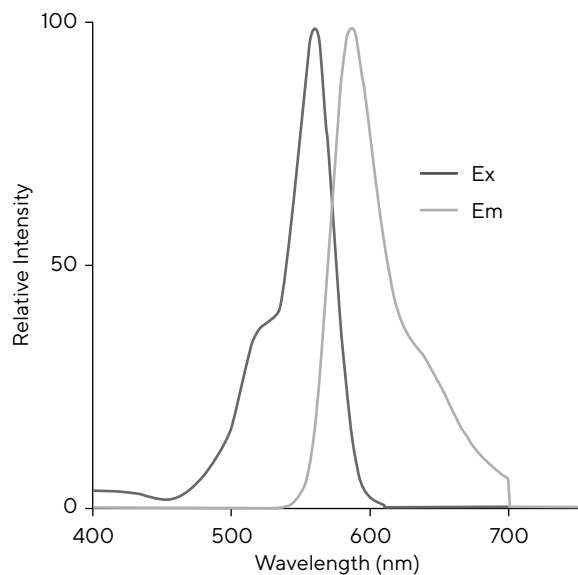


Figure 1

Excitation and Emission Spectra for the (A) pHrodo[®] Green and (B) pHrodo[®] Red Fluorophores, Determined in pH 4.0 Buffer

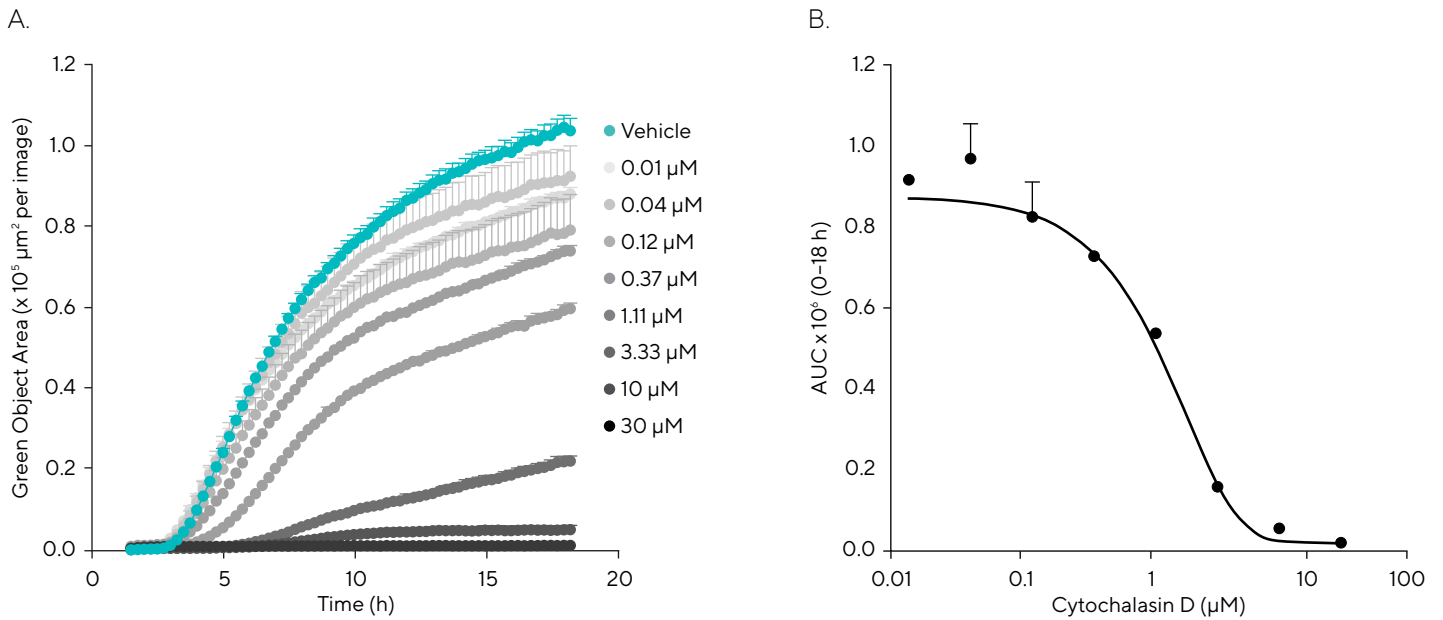


Figure 2
Concentration-Dependent Attenuation of pHrodo® Green *E. Coli* Bioparticles® Phagocytosis by the Actin Polymerization Inhibitor Cytochalasin D in J774A.1 Murine Macrophages

Note. (A) Time-course of phagocytosis in the absence (open symbols) and increasing concentrations of cytochalasin D (progressively darker gray symbols). Phagocytosis has been quantified as the fluorescence area for each time-point. (B) Concentration response curve to cytochalasin D. Area under the curve (AUC) values have been determined from the time-course shown in panel A (0–18 hours) and are presented as the mean ± SEM, n=3 wells.

Protocols and Procedures

Required Materials

- pHrodo® Red *E. coli* Bioparticles® for Incucyte® (Sartorius Cat. No. 4615)
or
- pHrodo® Green *E. coli* Bioparticles® for Incucyte® (Sartorius Cat. No. 4616)
or
- pHrodo® Red Zymosan Bioparticles® for Incucyte® (Sartorius Cat. No. 4617)
or
- pHrodo® Green Zymosan Bioparticles® for Incucyte® (Sartorius Cat. No. 4618)
or
- pHrodo® Red *S. aureus* Bioparticles® for Incucyte® (Sartorius Cat. No. 4619)
or
- pHrodo® Green *S. aureus* Bioparticles® for Incucyte® (Sartorius Cat. No. 4620)

General Guidelines

- We recommend medium with low levels of riboflavin to reduce the green fluorescence background. EBM, F12-K, and Eagles MEM have low riboflavin (< 0.2 mg/L). DMEM and RPMI have high riboflavin (> 0.2 mg/L).
- Following cell seeding, place plates at ambient temperature (15 minutes for adherent cell lines and 45 minutes for non-adherent cell lines) to ensure homogenous cell settling.
- When preparing the pHrodo® Bioparticles®, we recommend using a glass vial to vortex and sonicate the solution for 10–15 min (longer sonification may be required for Zymosan Bioparticles®).
- Remove bubbles from all wells by gently squeezing a wash bottle (containing 70–100% ethanol with the inner straw removed) to blow vapor over the surface of each well.
- After placing the plate in the Incucyte® Live-Cell Analysis System, allow the plate to warm to 37° C for 30 minutes prior to scanning.

Note: The mouse macrophage cell line J774A.1 was used to optimize the described conditions; however, the methodology can be adapted to accommodate any phagocyte.

1 Seed Target Cells

- 1.1 Seed phagocytic cells (50 μ L per well) at an appropriate density into a 96-well flat bottom plate (Corning Cat. No. 3595) such that by Day 1, the cell confluence is approximately 10%–20%. The seeding density will need to be optimized for cell type used; however, we have found that 1×10^3 to 1×10^4 cells per well are reasonable starting points.

Note: Phagocyte cell growth can be monitored by recording phase images using the Incucyte® Live-Cell Analysis System and confluence algorithm.

2 Treat Cells

- 2.1 Once the target cells have reached appropriate confluence, remove the cell plate from the incubator and add desired treatments. The volumes | dilutions may be varied; however, we recommended 25 μ L, prepared at 4X final assay concentration.
- 2.2 Incubate the treatments for the desired duration.

3 Prepare pHrodo® Bioparticles® and Add to Cells.

- 3.1 Prepare pHrodo® Bioparticles® for Incucyte® by resuspending to 1 mg/mL in PBS or complete media of choice. Transfer this solution to a glass vial, vortex and sonicate for a minimum of 10–15 minutes (longer sonication may be required for Zymosan).

Note: The formation of a homogeneous suspension may be improved by initial reconstitution in PBS, followed by subsequent dilution in assay media (PBS final assay concentration of 5%).

- 3.2 After incubation with the treatments, add the pHrodo® Bioparticles® for Incucyte® of your choice to the plate; we recommend 10 μ g per well for *E. coli*/*S. aureus* or 5 μ g for Zymosan.

Note: Remove bubbles at the liquid surface by gently squeezing a wash bottle (containing 100% ethanol with the inner straw removed) to blow vapor over the surface of each well.

4 Live-Cell Imaging

- 4.1 In the Incucyte® integrated software, schedule 24-hour repeat scanning for every 15 minutes, 2 images per well, for 2–48 hours (until the fluorescence area and intensity plateaus).
 - a. Scan on schedule, standard.
 - b. Channel selection: select “phase” and “red” or “green” (depending on Bioparticle® used)
 - c. Objective: 10X or 20X

For Research Use Only. Not for Therapeutic or Diagnostic Use.

Product Label License

This product is provided under an intellectual property license from Thermo Fisher Scientific. The transfer of this product is conditioned on the buyer using the purchased product solely in research conducted by the buyer, excluding contract research or any fee for service research, and the buyer must not sell or otherwise transfer this product or its components for (a) diagnostic, therapeutic or prophylactic purposes; (b) testing, analysis or screening services, or information in return for compensation on a per-test basis; (c) manufacturing or quality assurance or quality control, or (d) resale, whether or not resold for use in research. For information on purchasing a license to this product for purposes other than as described above, contact Thermo Fisher Scientific customer service.

North America

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
USA
Phone +1 734 769 1600

Europe

Sartorius UK Ltd.
Longmead Business Centre
Blenheim Road
Epsom
Surrey, KT19 9QQ
United Kingdom
Phone +44 1763 227400

Asia Pacific

Sartorius Japan K.K.
4th Floor, Daiwa Shinagawa North Bldg.
1-8-11, Kita-Shinagawa 1-chome
Shinagawa-Ku
Tokyo 140-0001
Japan
Phone +81 3 6478 5202