

## Lipid Coated Microparticles

Product #	Product Name	Lipid	Dye
P-B1PC	PC Lipid Microparticles, NBD	Phosphatidylcholine (PC)	NBD
P-B1PCr	PC Lipid Microparticles, pHrodo™ Red	Phosphatidylcholine (PC)	pHrodo™ Red
P-B1PCnf	PC Lipid Microparticles (no FL)	Phosphatidylcholine (PC)	Unlabeled
P-B1PCPS	PC/PS Lipid Microparticles, NBD	PC to PS ratio of 75:20	NBD
P-B1PE	PE Lipid Microparticles, NBD	Phosphatidylethanolamine (PE)	NBD
P-B1PS	PS Lipid Microparticles, NBD	Phosphatidylserine (PS)	NBD
P-B1PSr	PS Lipid Microparticles, pHrodo™ Red	Phosphatidylserine (PS)	pHrodo™ Red
P-B1PSnf	PS Lipid Microparticles (no FL)	Phosphatidylserine (PS)	Unlabeled

Support: echelon@echelon-inc.com

### Description:

Lipid microparticles are beads with attached phospholipids where the phospholipid headgroups are exposed and available for biological interactions.

### Properties:

Form – Bead suspension

Size – 100  $\mu$ L, 500  $\mu$ L

Storage instructions – Store product at 2-4 °C. Do not freeze.

Storage buffer – Microparticles are in PBS, pH 7.4., 0.02% sodium azide.

Concentration – See certificate of analysis.

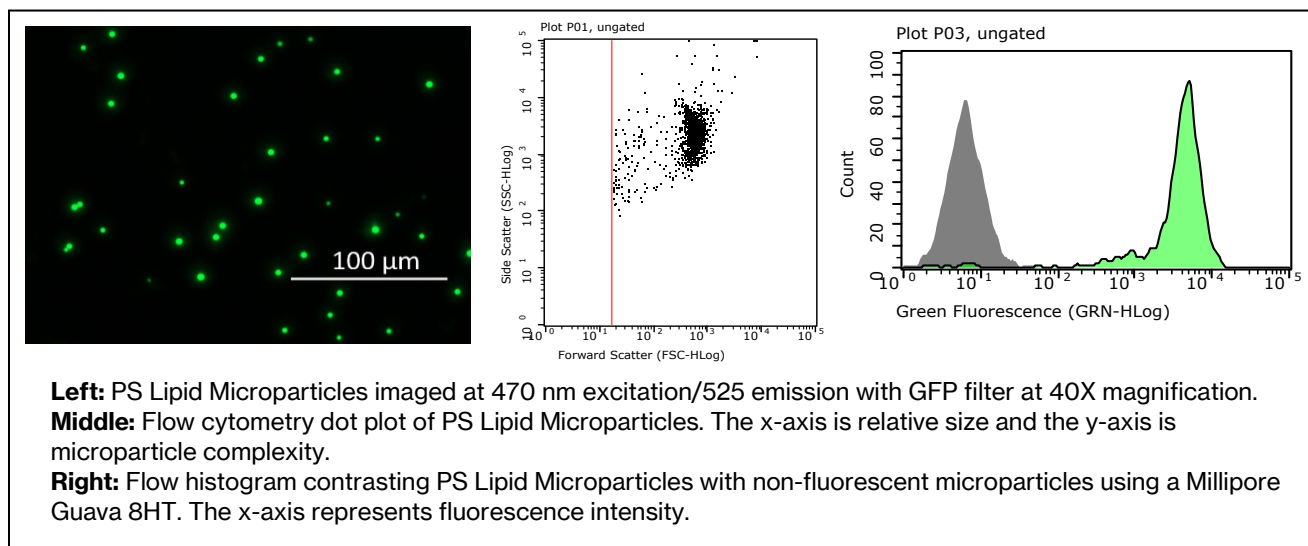
### Applications:

Lipid coated microparticles are intended for use in apoptosis and phagocytosis studies.

### Background:

Three lipids and two fluorescent dyes are available. Phosphatidylserine (PS) is an anionic, intracellular phospholipid component of the cell membrane and is involved in cell signaling. Phosphatidylcholine (PC) is primarily an extracellular phospholipid component of cell membranes and is a key structural lipid. Phosphatidylethanolamine (PE) is primarily found in the inner leaflet of the plasma membrane and plays a role in cell signaling, membrane structure, and fluidity.

### Data:



#### Technical Notes:

1. Total lipid concentration is approximately 1  $\mu\text{mol}$  per 1 mg microparticles.
2. Fluorescent containing microparticles contain trace amounts of a fluorescent lipid.
  - a. NBD has maximal excitation/emission at approximately 460/540 nm.
  - b. pHrodo™ Red has maximal excitation/emission at approximately 560/590 nm.
3. The microparticle is comprised of a silica core and is 3  $\mu\text{m}$ .
4. Lipid coated microparticles contain 0.02% sodium azide. If your experiments are sensitive to sodium azide, remove it before use. eg. 1 x PBS buffer, pH 7.4, followed by centrifugation @10,000 xg for 10 minutes.
5. Centrifuge the microparticles at 10,000 x g or lower. Higher centrifugation may damage microparticles.
6. AnnexinV demonstrates binding to the PS containing microparticles. AnnexinV does not bind the PC containing microparticles.
7. pHrodo is the trademark of ThermoFisher Scientific.

#### Preparation Notes:

We provide the following section as a starting point and strongly encourage researchers to consult the scientific literature and to conduct optimization experiments to establish the most favorable conditions for their experiments. Other buffers and conditions may show improved results.

1. Microparticles can aggregate with storage. Before each use, sonicate for 5-10 minutes in a water bath sonicator followed by mixing with pipette to ensure uniform suspension.
2. Pre-treating the microparticles with a blocking buffer reduced non-specific binding of AnnexinV to PC coated microparticles and appears to reduce non-specific uptake of the microparticles in RAW 264.7 macrophages. 10 mM HEPES, 140 mM NaCl, 2.5 mM  $\text{CaCl}_2$ , 5% BSA and 0.25% Tween20, followed by centrifugation at 10,000 X g for 10 minutes.
3. Dilute the microparticles with your preferred assay buffer to roughly  $2.5 - 1 \times 10^6$  microparticles / mL. This microparticle suspension should be further diluted 1:4 - 1:10 in cell media before addition to the cells. This is to dilute any item contained in the assay

buffer, such as detergent, that may negatively impact the cells. We added microparticles to the RAW 264.7 macrophages at a 1:4 ratio in cell media. This was roughly  $0.25 \times 10^6$  microparticles in 1 mL cell media.

4. After incubation with microparticles, we suggest removing non-ingested microparticles with washing. This can be followed by a quenching step.
  - a. For NBD containing microparticles, NBD can be quenched with 0.4% trypan blue.

#### References:

1. Neil Paterson, Tim Lämmermann (2022) Macrophage network dynamics depend on haptokinesis for optimal local surveillance eLife 11:e75354
2. Nabil Rabhi, Kathleen Desevin (2022) Obesity-induced senescent macrophages activate a fibrotic transcriptional program in adipocyte progenitors Life Science Alliance DOI: 10.26508/lsa.202101286
3. Brando Cieniewicz, Ankit Bhatta (2023) Chimeric TIM-4 receptor-modified T cells targeting phosphatidylserine mediates both cytotoxic anti-tumor responses and phagocytic uptake of tumor-associated antigen for T cell cross-presentation Molecular Therapy Vol. 31 No 7
4. Rachel Grazda, Allison N. Seyfried (2023) Impaired inflammation resolution in murine bone marrow failure is rescued by Resolvin E1 treatment bioRxiv preprint <https://doi.org/10.1101/2023.02>