

## Product Information

### Pluronic F-127®

Catalog Number	Unit Size
C020	2 g
C021	1 mL

#### Storage upon receipt:

- Room temperature

is normally kept at or below 0.1%. More weakly fluorescent indicators may require more concentrated loading solutions and correspondingly greater amounts of Pluronic® F-127. In general it is desirable to use the minimum amount of AM ester needed to achieve adequate fluorescence signal to noise levels. Loading may be done at any temperature that is tolerable for the cells. Note that the incubation temperature generally affects the extent of intracellular dye compartmentalization. After labeling, the cells are washed with **fresh medium** before beginning the experiment.

### Product Description

Pluronic F-127 is a nonionic, surfactant polyol (molecular weight approximately 12,500 daltons) that has been found to facilitate the solubilization of water-insoluble dyes and other materials in physiological media. Pluronic F-127 is commonly used to help disperse the acetoxymethyl (AM) esters of ion indicators as well as the lipophilic tracers.

### Guidelines for Use

#### Handling Solid Pluronic® F-127 and Stock Solutions

Dissolve 2 g of Pluronic® F-127 in 10 mL of anhydrous dimethyl sulfoxide (DMSO) to give a 20% (w/v) stock solution. This may require heating at ~40°C for about 20 minutes.

Store solution at room temperature. Do not refrigerate or freeze the solutions because Pluronic® F-127 may come out of solution. If the product does crystallize, it can be reheated until it goes back into solution. If precipitation is observed in purchased solutions of Pluronic® F-127, the product may be resolubilized by heating at ~40°C and vortexing before use.

#### Using Pluronic® F-127

The experimental conditions for loading cells with AM esters varies with cell type due to differences both in probe uptake and in the intracellular esterase activity required for hydrolysis of the AM esters. Solutions of the AM esters in DMSO must be kept anhydrous since the solvent will readily take up moisture, leading to loss of cell-loading efficacy. Pluronic® F-127 should be added only to working solutions. Typically, a small volume of the AM ester, dissolved at 1–5 mM in DMSO, is mixed with the 20% (w/v) Pluronic® F-127 stock solution in DMSO at a ratio of 1:1 immediately before use. The solution of AM ester and Pluronic® F-127 is then diluted into the cell-loading buffer to achieve a final AM ester concentration of between 1 µM and 10 µM and the cells are incubated for between 10 minutes and 1 hour. The final concentration of Pluronic® F-127