

Product Information

EasyProbes™ Propidium Iodide Dead Cell Stain

Catalog Number	Unit Size
C059	6 × 2.5 mL

Storage upon receipt:

- 2-25°C
- Protect from light

Ex/Em: 535/617 nm, bound to DNA

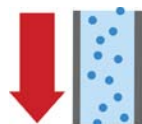
Quick Start Guide



1. Add 2 drops per 10^6 cells in 1 mL



2. Incubate 15-30 minutes at RT



3. Proceed with flow cytometry