
Product Information

EasyProbes™ Green 488 Dead Cell Stain

Catalog Number	Unit Size
C063	6 × 2.5 mL

Storage upon receipt:

- 2-8°C
- Protect from light

Ex/Em: 500/530 nm, bound to DNA

Quick Start Guide



1. Add 2 drops per mL of media



2. Incubate 5-15 minutes at RT



3. Image cells