
Product Information

EasyProbes™ Hoechst 33342 Live Cell Stain

Catalog Number	Unit Size
C057	6 × 2.5 mL

Storage upon receipt:

- 2-25°C
- Protect from light

Ex/Em: 350/461 nm, bound to DNA

Quick Start Guide



1. Add 2 drops per mL of media



2. Incubate 15-30 minutes at RT



3. Image cells