

Product Information

EasyProbes™ Antifade Reagent with DAPI

Catalog Number	Unit Size
C068	10 mL
C069	5 × 2 mL

Storage upon receipt:

- -20~4°C
- Protect from light

The product is stable for at least 6 months.

Quick Start Guide



1. Add 1 drop of EasyProbes™ Antifade Reagent onto glass slide.



2. Remove excess liquid from coverslip.



3. Place coverslip sample-side down onto antifade reagent onto the glass slide.



4. Apply nail polish or hot wax to edges of coverslip to secure it.
5. Clean coverslip and image cells.