

EasyProbes™ Cell Viability Imaging Kit (Blue/Green)

Catalog Number: A064

Table 1. Kit Components and Storage

Material	Amount	Storage	Stability
EasyProbes™ Hoechst 33342 Live Cell Stain	3×2.5 mL	2~8 °C	The product is stable for one year when stored as directed.
EasyProbes™ Green 488 Dead Cell Stain	3×2.5 mL	2~8 °C	

Approximate fluorescence excitation/emission maxima, in nm: Hoechst 33342: 350/461; Green 488 Dead Cell Stain: 500/530, bound to DNA.

Introduction

EasyProbes™ Cell Viability Imaging Kit (Blue/Green) is a ready-to-use assay that can quickly and easily determine the viability of cells. EasyProbes™ Hoechst 33342 reagent stains the nuclei of all the cells and can be detected with a standard DAPI filter. EasyProbes™ Green 488 Dead reagent stains only the nuclei of cells with compromised plasma membrane integrity and is detected using a standard FITC (green) filter set. This kit is appropriate for fluorescence microscopy and flow cytometry.

Quick Start Guide



1. Culture cells in appropriate media and vessel for microscopy



2. Add 2 drops each Hoechst 33342 and Green 488 Dead reagents per mL of media to label cells



3. Incubate 10-30 minutes at RT



4. Image cells