

## EasyProbes™ Cell Viability Imaging Kit (Blue/Red)

Catalog Number: A065

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™ Propidium Iodide Dead Cell Stain	3×2.5 mL	2~8 °C	

**Approximate fluorescence excitation/emission maxima, in nm:** Hoechst 33342: 350/461; Propidium Iodide: 535/617, bound to DNA.

### Introduction

EasyProbes™ Cell Viability Imaging Kit (Blue/Red) 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™ Propidium Iodide reagent stains only the nuclei of cells with compromised plasma membrane integrity and is detected using a standard TRITC/RFP (orange) 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 Propidium Iodide reagents per mL of media to label cells



3. Incubate 10-30 minutes at RT



4. Image cells