

## ReadiDrop™ Cell Viability Assays



Bio-Rad's ReadiDrop Cell Viability Assays are designed for ease of use in cell sorting and cell analysis applications. Simply add one or two drops of the cell viability assay to assess the health of your cells. ReadiDrop Assays are available with the two most commonly used dyes, propidium iodide and 7-aminoactinomycin D (7-AAD), for dead cell exclusion.

### Key Benefits of ReadiDrop Cell Viability Assays

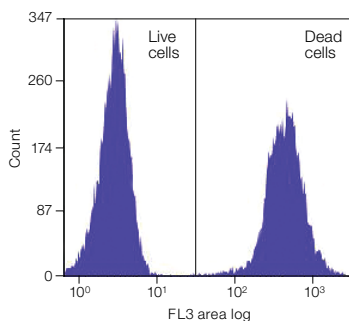
- **Ideal for cell sorting** — fluorescent molecules are suspended in phosphate buffered saline (PBS) without preservatives
- **Stable at room temperature** — keep a bottle of any ReadiDrop Cell Viability Assay by your S3™ Cell Sorter or your cell analyzer for easy access
- **No more weighing, pipetting, and diluting** — ready-to-use solution is provided at a final concentration of 1 µg/ml

Visit [www.bio-rad.com/web/ReadiDrop1](http://www.bio-rad.com/web/ReadiDrop1) for more information.

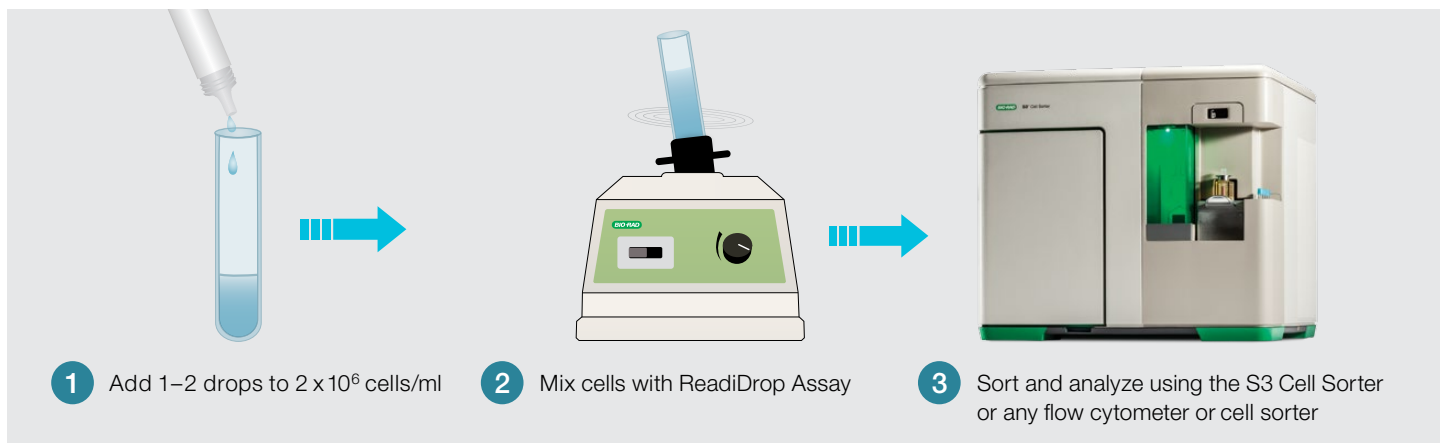
**BIO-RAD**

## Simplifying Cell Viability

Excluding dead cells in a flow cytometry experiment is critical for accurate result interpretation (Figure 1). One identifier of a dead cell is the collapse of the cell membrane, which allows impermeable dyes, such as propidium iodide and 7-AAD, to enter. Using ReadiDrop Cell Viability Assays, easily evaluate the health of your cells in three easy steps (Figure 2).



**Fig. 1. Easily identify dead cells in your experiment with ReadiDrop Cell Viability Assays.** A mixture of live and heat-killed Jurkat cells were stained with ReadiDrop Propidium iodide for 1 min prior to sorting.



**Fig. 2. Evaluate the health of your cells in three steps.**

## Related Products

### Fixable Cell Viability Dyes

Easily incorporate a VivaFix™ Cell Viability Assay into your flow cytometry experiment without losing fluorescence intensity.

Catalog Number	Labeling Dye Description	Optimal Excitation Laser, nm	S3 Cell Sorter (488/561 nm) Filter Channel
135-1111	VivaFix 353/442	355	—
135-1112	VivaFix 410/450	405	—
135-1113	VivaFix 408/512	405	—
135-1114	VivaFix 398/550	405	—
135-1115	VivaFix 498/521	488	FL1 (525/30)
135-1116	VivaFix 547/573	561	FL2 (586/25)
135-1117	VivaFix 583/603	561	FL3 (615/25)
135-1118	VivaFix 649/660	640	—

## Ordering Information

Catalog Number	ReadiDrop Cell Viability Assay	Excitation Maximum, nm	Emission Maximum, nm	S3 Cell Sorter (488/561 nm) Optimal Filter Set
135-1101	ReadiDrop Propidium Iodide	538	617	FL3 (585/25)
135-1102	ReadiDrop 7-AAD	548	648	FL4 (655LP)



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [www.bio-rad.com](http://www.bio-rad.com) **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 11 3065 7550  
**Canada** 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00  
**France** 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300  
**Israel** 03 963 6050 **Italy** 39 02 216091 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666  
**New Zealand** 64 9 415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04  
**Singapore** 65 6415 3188 **South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 026 674 55 05  
**Taiwan** 886 2 2578 7189 **Thailand** 1800 88 22 88 **United Kingdom** 020 8328 2000