

ReadiDrop™ Cell Viability Assays

Bio-Rad's ReadiDrop Cell Viability Assays are designed for ease of use in cell sorting and cell analysis applications. Excluding dead cells in a flow cytometry experiment is critical for accurate result interpretation. When the cell membrane collapses, impermeant dyes, such as propidium iodide (PI) and 7-aminoactinomycin D (7-AAD), can enter and bind to double-stranded DNA. Simply add one or two drops to assess the health of your cells.



- Phosphate buffered saline (PBS) solution is gentle on cells
- Room temperature stability allows easy storage
- No weighing, pipetting, and diluting needed

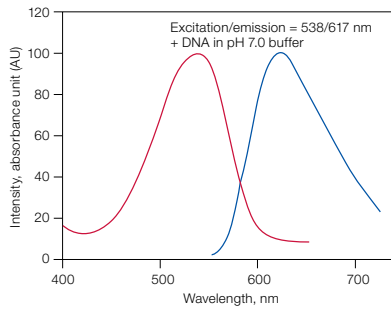
Visit bio-rad.com/web/ReadiDrop1 for more information.



Ideal for Cell Sorting and Flow Cytometry

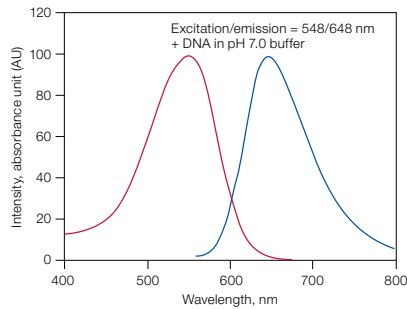
ReadiDrop Propidium Iodide

Cell impermeant nucleic acid-binding dye that binds to double-stranded DNA/RNA.

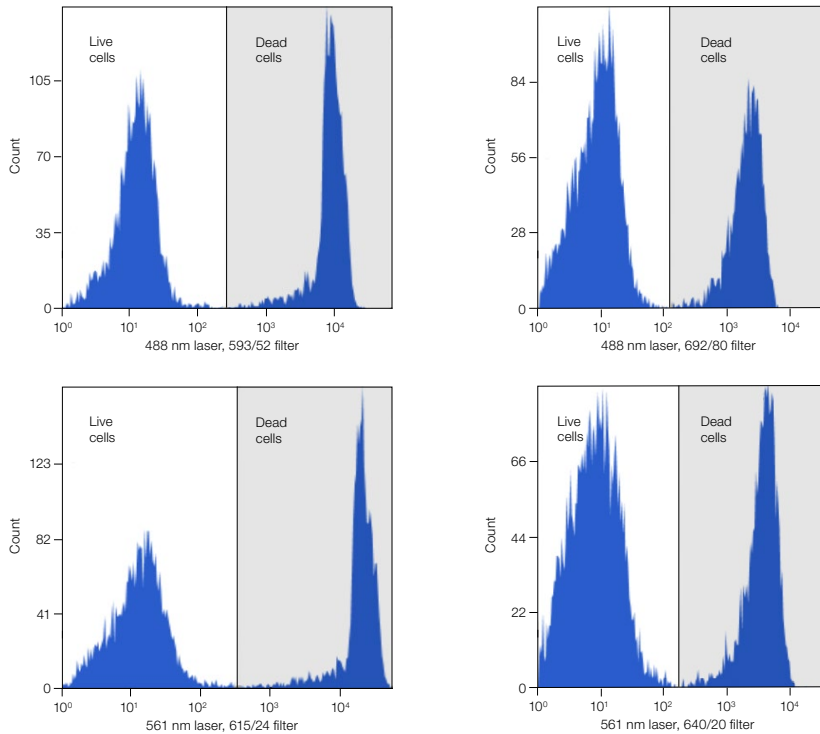


ReadiDrop 7-AAD

Cell impermeant DNA-binding dye that has a high affinity for GC-rich regions of double-stranded DNA.



Excitation and emission curves for propidium iodide and 7-AAD.



Assessment of cell viability using ReadiDrop reagents and the ZE5™ Cell Analyzer. Jurkat cells were heat shocked at 65°C for 5 min. Cells were washed and resuspended at 2×10^6 cells/ml in PBS. Twenty μ l of ReadiDrop Propidium Iodide or 7-AAD were added to 0.5 ml cells and incubated for 15 min at room temperature in the dark. Cells were washed once in PBS prior to acquisition on the ZE5 Cell Analyzer.

Ordering Information

Catalog #	Description	Maximum Excitation, nm	Maximum Emission, nm	Optimal Excitation Laser, nm	S3e™ Cell Sorter	ZE5 Cell Analyzer	ZOE™ Fluorescent Cell Imager
1351101	ReadiDrop Propidium Iodide	538	617	561 or 488	•	•	
1351102	ReadiDrop 7-AAD	548	648	561 or 488	•	•	

Each assay includes 3 x 3 ml bottles of ready-to-use reagent suspended in PBS for dead cell exclusion in cell sorting and flow cytometry applications.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 01 877 89019 **Belgium** 32 03 710 53 00 **Brazil** 55 11 3065 7550
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 36 01 459 6192 **Denmark** 45 04 452 10 00 **Finland** 35 08 980 422 00 **France** 33 01 479 593 00
Germany 49 089 3188 4393 **Hong Kong** 852 2789 3300 **Hungary** 36 01 459 6190 **India** 91 124 4029300 **Israel** 972 03 963 6050 **Italy** 39 02 49486600
Japan 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 0 318 540 666 **New Zealand** 64 9 415 2280 **Norway** 47 0 233 841 30
Poland 36 01 459 6191 **Portugal** 351 21 4727717 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 36 01 459 6193 **Spain** 34 091 49 06 580
Sweden 46 08 555 127 00 **Switzerland** 41 0617 17 9555 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300
United Kingdom 44 01923 47 1301

