

BactoView™ Dead Stains

Highly-selective dead cell stains for bacteria cultures. The stains are suitable for both gram-positive and gram-negative strains.



Product Description

BactoViewTM Dead Stains are novel DNA binding dyes for live/dead discrimination in bacteria. These bright fluorogenic DNA binding dyes are cell membrane-impermeant, so they selectively stain dead bacteria with compromised cell membranes without the need to wash after staining.

- Highly-selective dead cell stains for bacteria cultures
- Superior live/dead differentiation for gram-positive strains than traditional dyes
- Also provide excellent results in gram-negative strains
- Quick 30-minute staining, no wash required
- Available in 7 colors from green to near-IR

Note: As of 6/16/2025, the formulation of BactoView™ Dead Stains (Cat. No. 40107-40113) was updated from 500X in water to 1000X in DMSO.

Highly-Selective Dead Cell Staining for Gram-Positive and Gram-Negative Strains

Traditional vital nucleic acid dyes like propidium iodide or ethidium homodimer are efficiently excluded from live gram-negative bacteria but are taken up by live gram-positive bacteria, resulting in high background in live cells and poor live/dead discrimination. BactoViewTM Dead Stains have novel chemical structures that are efficiently excluded from both gram-positive and gram-negative strains, for highly selective live/dead discrimination. In addition to staining dead bacteria, BactoViewTM Dead Stains also stain Bacillus subtilis endospores, but with dimmer fluorescence. BactoViewTM Dead Stains have low fluorescence until they bind DNA, allowing bright, no-wash staining. The stains are available with a wide selection of emission wavelengths ranging from green to near-infrared, for microscopy or flow cytometry analysis.

Call us: 800-304-5357

Product attributes

Apoptosis/viability marker	Dead cell stain
For live or fixed cells	For live/intact cells
Detection method/readout	Fluorescence microscopy, Flow cytometry
Assay type/options	Endpoint assay
Colors	Green, Red, Far-red, Near-infrared
Storage Conditions	Store at -10 to -35 °C, Protect from light

BactoView™ Dead Stains & Viability Kits

Product Name	Ex/Em (nm)	Detection Channel	Size (1000X in DMSO)	Catalog No.
BactoView TM Dead 500/515	497/515	FITC	20 uL	<u>40107-T</u>
100 uL	<u>40107</u>			
BactoView™ Dead 560/570	559/570	Rhodamine, PI, PE	20 uL	<u>40108-T</u>
100 uL	<u>40108</u>			
BactoView™ Dead 570/585	572/583	Rhodamine, PI, PE	20 uL	<u>40109-T</u>
100 uL	<u>40109</u>			
BactoView™ Dead 600/615	603/613	Texas Red® or PE-Texas Red®	20 uL	<u>40110-T</u>
100 uL	<u>40110</u>			
BactoView™ Dead 655/670	653/671	Cy®5, APC	20 uL	<u>40111-T</u>
100 uL	<u>40111</u>			
BactoView™ Dead 690/710	683/707	Cy®5.5	20 uL	<u>40112-T</u>
100 uL	<u>40112</u>			
BactoView™ Dead 760/780	759/780	Cy®7, APC-Cy®7	20 uL	40113-T
100 uL	<u>40113</u>			
BactoView™ Viability Kit (Green/Red)	Green (498/522)/Red (572/583)	FITC (Green)/Rhodamine, PI, PE (Red)	1 kit	<u>32019</u>
BactoView™ Viability Kit (Green/Far-Red)	Green (498/522)/Far-Red (653/671)	FITC (Green)/Cy®5, APC (Far-Red)	1 kit	32020

Texas Red is a registered trademark of Thermo Fisher Scientific; Cy Dye is a registered trademark of Cytiva.

Staining of Gram-Negative E. Coli

Staining of Gram-Positive B. Subtilis

BactoView™ Dead Also Stains B. Subtilis Endospores

Biotium also offers BactoView™ Viability Kits, which include a choice of red or far-red BactoView™ Dead Stain for dead bacteria and BactoView™ Viability Green Counterstain to stain all bacteria. BactoView™ Dead Stains also can be combined with fluorescent Gram stains like our CF® Dye WGA Conjugates.

We also offer <u>BactoViewTM Live Green</u> and <u>BactoViewTM Live Red</u> for staining live bacteria. Note that BactoViewTM Stains cannot be used to distinguish bacteria from eukaryotic cells, because they will stain other cell types as well. For bright and optimized labeling of bacterial endospores, see our <u>BactoSporeTM Bacterial Stains</u>.

For staining mammalian cells, see our NucSpot® Nuclear Stains for live/dead discrimination or nuclear counterstaining of fixed mammalian cells. For live nuclear staining of mammalian cells, see our NucSpot® Live Stains. Also, view our Cellular Stains Table for more information on how our dyes stain various organisms.

This datasheet was generated on November 20, 2025 at 08:03:34 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/bactoview-dead-stains/