BactoView™ Viability Kits

A two-color viability staining kit optimized specifically for gram-positive bacteria strains. Offers a simple 30-minute staining protocol, no wash required.



Call us: 800-304-5357

Product Description

BactoViewTM Viability Kits are optimized specifically for viability staining of gram-positive bacteria. The kits include Biotium's novel BactoViewTM Stains for two-color staining of live and dead cells. BactoViewTM Dead Stains are novel membrane-impermeant DNA binding dyes that selectively stain dead bacteria with compromised cell membranes. BactoViewTM Viability Green Counterstain labels both live and dead bacteria with green fluorescence and is formulated for two-color staining with BactoViewTM Dead.

Note: As of 6/16/2025, the formulation of kit components BactoView[™] Dead 570/585 (Cat. No. 40109) and BactoView[™] Dead 655/670 (Cat. No. 40111) was updated from 500X in water to 1000X in DMSO.

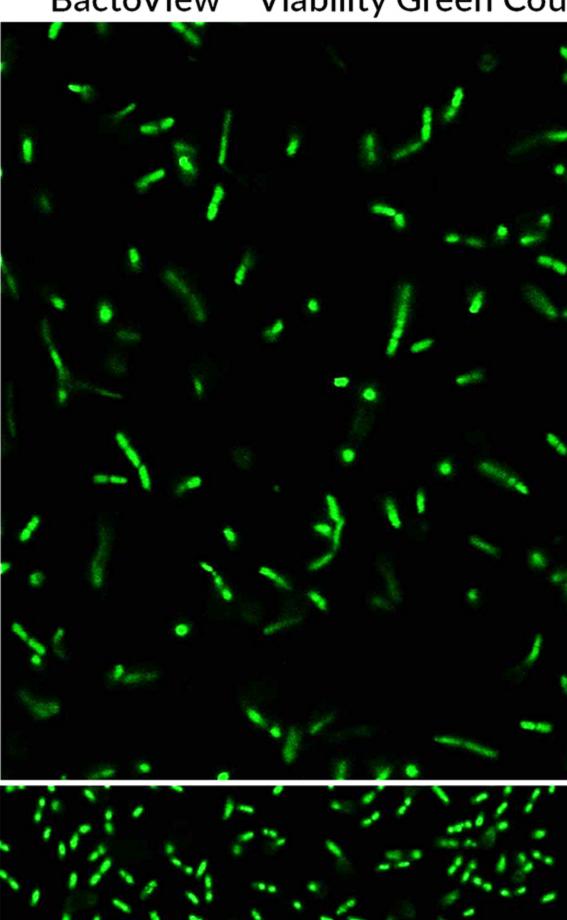
Robust Viability Assessment for Gram-Positive Bacteria

BactoViewTM Viability Kits are optimizes specifically for live/dead staining of gram-positive strains. The kits include a highly selective BactoViewTM Dead Stain for staining dead cells, and a BactoViewTM Viability Green Counterstain that labels both live and dead bacteria. The kits are available with your choice of dye combination for green/visible red or green/far-red fluorescence detection. The stains are fluorogenic for no-wash staining and can be used to stain cells in culture medium or buffer.

Traditional vital nucleic acid dyes like propidium iodide or ethidium homodimer are efficiently excluded from live gram-negative bacteria, which have an outer membrane protecting the cell wall. But these dyes often are taken up by live gram-positive bacteria that lack an outer membrane, 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. See our full selection of BactoViewTM Dead Stains in colors ranging from green to near-infrared fluorescence in the table below.

Note: While the BactoView™ Viability Kits can be used with gram-negative strains or mixed cultures of gram-positive and gram-negative bacteria, our <u>Bacterial Viability and Gram Stain Kit</u> will provide brighter staining of live gram-negative cells than the BactoView Kits.

BactoView™ Viability Green Cou



Untreated B. subtilis

Product attributes

Apoptosis/viability marker	Dead cell stain, All 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		
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™ 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	<u>40113-T</u>
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.

We also offer BactoView™ Live stains with green or red fluorescence. Note that BactoView™ 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 BactoSpore™ Bacterial Stains. 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 08:03:37 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/bactoview-viability-kits/