

## 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.



### Product Description

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

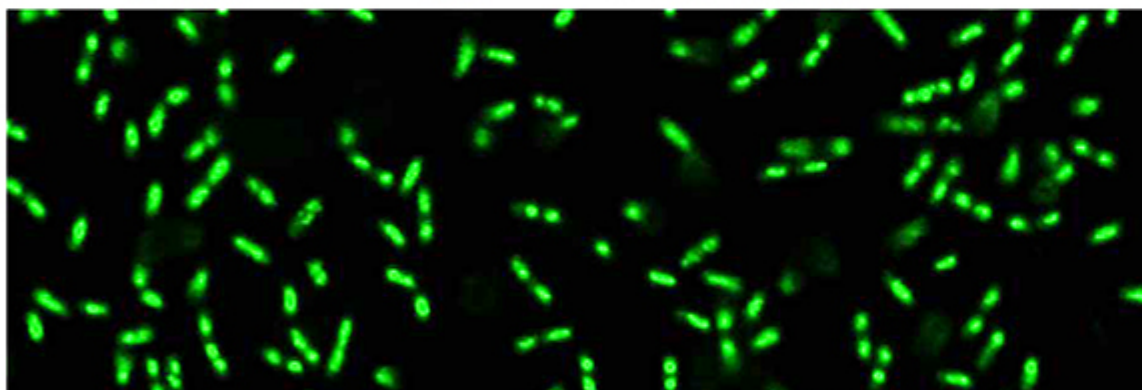
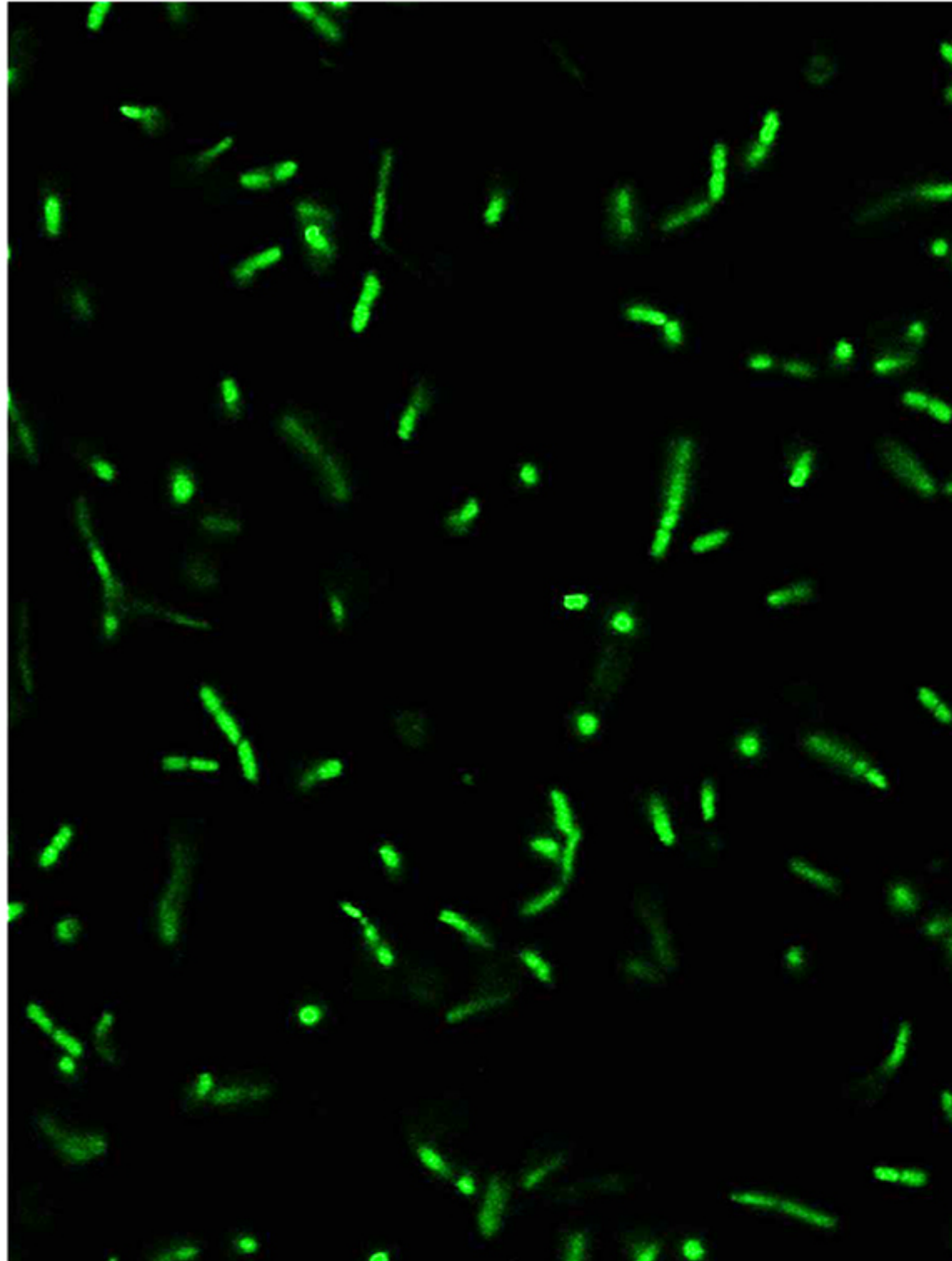
BactoView™ Viability Kits are optimized specifically for live/dead staining of gram-positive strains. The kits include a highly selective BactoView™ Dead Stain for staining dead cells, and a BactoView™ Viability Green Counterstain that labels both live and dead bacteria. The kits are available with your 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. [BactoView™ 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 [BactoView™ 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 [Bacterial Viability and Gram Stain Kit](#) will provide brighter staining of live gram-negative cells than the BactoView Kits.

BactoView™ Viability Green Cou

Untreated  
*B. subtilis*



## Product attributes

<b>Apoptosis/viability marker</b>	Dead cell stain, All cell stain
<b>For live or fixed cells</b>	For live/intact cells
<b>Detection method/readout</b>	Fluorescence microscopy, Flow cytometry
<b>Assay type/options</b>	Endpoint assay
<b>Colors</b>	Green, Red, Far-red
<b>Storage Conditions</b>	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.
<a href="#">BactoView™ Dead 500/515</a> 100 uL	497/515 <a href="#">40107</a>	FITC	20 uL	<a href="#">40107-T</a>
<a href="#">BactoView™ Dead 560/570</a> 100 uL	559/570 <a href="#">40108</a>	Rhodamine, PI, PE	20 uL	<a href="#">40108-T</a>
<a href="#">BactoView™ Dead 570/585</a> 100 uL	572/583 <a href="#">40109</a>	Rhodamine, PI, PE	20 uL	<a href="#">40109-T</a>
<a href="#">BactoView™ Dead 600/615</a> 100 uL	603/613 <a href="#">40110</a>	Texas Red® or PE-Texas Red®	20 uL	<a href="#">40110-T</a>
<a href="#">BactoView™ Dead 655/670</a> 100 uL	653/671 <a href="#">40111</a>	Cy®5, APC	20 uL	<a href="#">40111-T</a>
<a href="#">BactoView™ Dead 690/710</a> 100 uL	683/707 <a href="#">40112</a>	Cy®5.5	20 uL	<a href="#">40112-T</a>
<a href="#">BactoView™ Dead 760/780</a> 100 uL	759/780 <a href="#">40113</a>	Cy®7, APC-Cy®7	20 uL	<a href="#">40113-T</a>
<a href="#">BactoView™ Viability Kit (Green/Red)</a>	Green (498/522)/Red (572/583)	FITC (Green)/Rhodamine, PI, PE (Red)	1 kit	<a href="#">32019</a>
<a href="#">BactoView™ Viability Kit (Green/Far-Red)</a>	Green (498/522)/Far-Red (653/671)	FITC (Green)/Cy®5, APC (Far-Red)	1 kit	<a href="#">32020</a>

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 June 8, 2026 at 12:15:50 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/bactoview-viability-kits/>