



Glowing products for science

## 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

BactoView™ 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. 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. In addition to staining dead bacteria, BactoView™ Dead Stains also stain *Bacillus subtilis* endospores, but with dimmer fluorescence. BactoView™ 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](tel:800-304-5357)

### Product attributes

<b>Apoptosis/viability marker</b>	Dead 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, Near-infrared
<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>	497/515	FITC	20 uL	<a href="#">40107-T</a>
100 uL	<a href="#">40107</a>			
<a href="#">BactoView™ Dead 560/570</a>	559/570	Rhodamine, PI, PE	20 uL	<a href="#">40108-T</a>
100 uL	<a href="#">40108</a>			
<a href="#">BactoView™ Dead 570/585</a>	572/583	Rhodamine, PI, PE	20 uL	<a href="#">40109-T</a>
100 uL	<a href="#">40109</a>			
<a href="#">BactoView™ Dead 600/615</a>	603/613	Texas Red® or PE-Texas Red®	20 uL	<a href="#">40110-T</a>
100 uL	<a href="#">40110</a>			
<a href="#">BactoView™ Dead 655/670</a>	653/671	Cy®5, APC	20 uL	<a href="#">40111-T</a>
100 uL	<a href="#">40111</a>			
<a href="#">BactoView™ Dead 690/710</a>	683/707	Cy®5.5	20 uL	<a href="#">40112-T</a>
100 uL	<a href="#">40112</a>			
<a href="#">BactoView™ Dead 760/780</a>	759/780	Cy®7, APC-Cy®7	20 uL	<a href="#">40113-T</a>
100 uL	<a href="#">40113</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.

## 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 [BactoView™ Live Green](#) and [BactoView™ Live Red](#) for staining live bacteria. 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 January 8, 2026 at 12:34:45 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/bactoview-dead-stains/>