



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

## Viability/Cytotoxicity Assay Kit for Bacteria Live & Dead Cells



This kit provides two-color fluorescence staining of both live bacteria (green) and dead bacteria (red) using two DNA dyes, DMAO and EthD-III.

### Product Description

This kit provides two-color fluorescence staining of both live bacteria (green) and dead bacteria (red) using two DNA probes, DMAO and EthD-III.

DMAO is a green-fluorescent nucleic acid dye which stains both live and dead bacteria. Ethidium Homodimer III (EthD-III) is a red-fluorescent nucleic acid dye that selectively stains dead bacteria with damaged cell membranes. When stained with a mixture of DMAO and EthD-III, bacteria with intact cell membranes show green fluorescence, whereas bacteria with damaged cell membranes show red fluorescence. The kit is suitable for use with fluorescence microscopes and flow cytometers. The assay principles are general and applicable to most bacteria types. For 100 flow cytometry assays or 1000 microscopy assays.

Biotium also offers [BactoView™ Dead Stains](#) and [BactoView™ Viability Kits](#) for highly-selective staining of dead bacteria, available in a selection of 7 colors from green to near-IR. BactoView™ Dead Stains also can be combined with fluorescent Gram stains like our [CF® Dye WGA Conjugates](#).

[BactoView™ Live Green](#) and [BactoView™ Live Red](#) are designed for staining both live and dead 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.

See our full selection of [Microbiology Stains & Kits](#).

This datasheet was generated on June 17, 2026 at 12:20:28 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/viabilitycytotoxicity-assay-kit-for-bacteria-live-and-dead-cells/>

### Product attributes

Apoptosis/viability marker	Live/dead assay
For live or fixed cells	For live/intact cells
Detection method/readout	Fluorescence microscopy
Assay type/options	Endpoint assay, Homogeneous assay, No-wash staining, Real-time imaging
Colors	Green/Red
Storage Conditions	Store at 2 to 8 °C, Protect from light