BactoView™ Live Fluorescent Bacterial Stains

utotum

Live cell bacteria stains, suitable for flow cytometry or fluorescence microscopy. Stain live and dead, gram-positive and gram-negative bacteria with green or red fluorescence.

Product Description

BactoViewTM Live dyes are fluorogenic live cell bacterial DNA stains. Bright and cell-permeant, they are able to stain both live and dead gram-positive and gram-negative bacteria for microscopy or flow cytometry.

- DNA dyes for bacteria
- Stain live and dead, gram-positive and gram-negative cells
- Choose green or red fluorescence
- Supplied at 500X in DMSO

BactoViewTM Live dyes can stain both live and dead bacteria; dead cells tend to stain more brightly than live cells. BactoViewTM Live Green can stain mammalian cell nuclei, but BactoViewTM Live Red shows weak mitochondrial signal in mammalian cells. In yeast, staining with either BactoViewTM Live dye is weak and not nuclear.

Biotium also offers <u>BactoViewTM Dead Stains</u> and <u>BactoViewTM Viability Kits</u> for highly-selective staining of dead bacteria, available in a selection of 7 colors from green to near-IR. BactoViewTM Dead Stains also can be combined with fluorescent Gram stains like our <u>CF® Dye WGA Conjugates</u>. For bright and optimized labeling of bacterial endospores, see our <u>BactoSporeTM Bacterial Stains</u>.

See our full selection of Microbiology Stains & Kits.

BactoView™ Live Dyes

Product Name	Ex/Em (nm) (with DNA)	Detection Channel	Size (500X in DMSO)	Catalog No.
BactoView TM Live Green	500/520	FITC	20 uL	<u>40102-T</u>
			100 uL	<u>40102</u>
BactoView™ Live Red	572/675	Cy®3 (microscopy) PE-Cy®5 (flow cytometry)	20 uL	<u>40101-T</u>
			100 uL	<u>40101</u>

Cy Dye is a registered trademark of Cytiva.

This datasheet was generated on November 15, 2025 at 02:36:39 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/bactoview-live/

Product attributes

Dye	BactoView™ Live Green, BactoView™ Live Red
Cell permeability	Membrane permeant
Colors	Green, Red
Excitation/Emission	BactoView™ Live Green: 500/520 nm (with DNA), BactoView™ Live Red: 572/675 nm (with DNA)
Concentration	500X in DMSO
Storage Conditions	Store at 2 to 8 °C. Protect from light

Call us: 800-304-5357 Email: btinfo@biotium.com