

BactoView™ Live Fluorescent Bacterial Stains

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 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 |

Product Description

BactoView™ 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

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

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](#). For bright and optimized labeling of bacterial endospores, see our [BactoSpore™ Bacterial Stains](#).

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™ Live Green | 500/520 | FITC | 20 uL | 40102-T |
| 100 uL | 40102 | | | |
| BactoView™ Live Red | 572/675 | Cy@3 (microscopy) PE-Cy@5 (flow cytometry) | 20 uL | 40101-T |
| 100 uL | 40101 | | | |

Cy Dye is a registered trademark of Cytiva.

This datasheet was generated on June 8, 2026 at 12:50:44 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/bactoview-live/>