



Glowing products for science

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

## Live Bacterial Gram Stain Kit

A fluorescent staining kit for live cell gram staining in bacteria.  
Suitable for fluorescence microscopy and flow cytometry.



### Product attributes

## Product Description

The Live Bacterial Gram Stain Kit is a fluorescent staining kit for gram staining in bacteria. The assay is suitable for detection by fluorescence microscopy or flow cytometry.

WGA binds specifically to N-acetylglucosamine in the peptidoglycan layer of gram-positive bacteria. The Live Bacterial Gram Stain Kit contains WGA conjugated to CF@594 to stain the surface of gram-positive cells with red fluorescence. CF@594 is spectrally similar to Texas Red®, but with brighter fluorescence. The kit also includes DAPI to stain the DNA of all bacteria with blue fluorescence. Gram-positive bacteria stained using this kit will show a fluorescent blue interior and fluorescent red cell surface, while gram-negative bacteria will be stained with blue fluorescence only.

This kit is suitable for detection by fluorescence microscopy and flow cytometry.

We also offer the [Bacterial Viability and Gram Stain Kit \(32001\)](#) that also includes a dead-cell dye to simultaneously measure cell viability.

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](#).

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

Texas Red is a registered trademark of Thermo Fisher Scientific.

## References

Science Journal of Chemistry (2018) 6;5 pp77-82 [doi:10.11648/j.sjc.20180605.11](https://doi.org/10.11648/j.sjc.20180605.11)

This datasheet was generated on January 11, 2026 at 08:50:13 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/live-bacterial-gram-stain-kit-800-assays/>