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

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

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 <u>Bacterial Viability and Gram Stain Kit (32001)</u> that also includes a dead-cell dye to simultaneously measure cell viability.

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>.

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

This datasheet was generated on November 11, 2025 at 06:53:25 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/live-bacterial-gram-stain-kit-800-assays/