



Glowing products for science

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

Yeast Viability Staining Kit

A rapid and selective fluorescent assay to determine yeast viability in microscopy or flow cytometry applications. Contains two fluorescent components, CF® Dye-labeled Con A and Live-or-Dye™.



Product Description

The Yeast Viability Staining Kit is a fluorescent assay to determine yeast viability. Each kit contains two fluorescent components: CF® Dye-labeled Concanavalin A (Con A) and Live-or-Dye™ viability dye. We offer 3 different color combinations to choose from.

The kit has been validated in the following species: *Saccharomyces cerevisiae*, *Candida albicans*, *Aspergillus niger*, and *Zygosaccharomyces bailii*.

If there is a different color combination that you would prefer, email us at techsupport@biotium.com.

How It Works

When both dyes are used to stain yeast, live cells display bright fluorescent cell wall staining (Con A), while dead cells with compromised membrane activity exhibit both cell wall staining and cytoplasmic staining (Live-or-Dye™). The dye labeling is extremely stable, allowing the cells to be fixed and permeabilized without loss of fluorescence or dye transfer between cells.

Live-or-Dye™ viability dyes are cell membrane impermeable amine-reactive dyes designed for discrimination between live and dead cells during flow cytometry or microscopy. They are able to enter into dead cells that have compromised membrane integrity and covalently label free amines on intracellular proteins.

Concanavalin A (Con A) is a widely used lectin that selectively binds to α -mannopyranosyl and α -glucopyranosyl residues, found in the cell wall of yeast and fungi. It is labeled with one of our superior bright and photostable CF® Dyes.

Product attributes

Colors	Green/Red, Red/Blue, Red/Far-red
Cell permeability	Membrane impermeant
For live or fixed cells	For live/intact cells
Probe cellular localization	Cytoplasmic, Membrane/cell surface
Detection method/readout	Fluorescence microscopy, Flow cytometry

This datasheet was generated on January 18, 2026 at 08:01:49 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/yeast-viability-staining-kit/>