Yeast Vitality Staining Kit

A fluorogenic assay that allows you to determine whether your yeast cells are metabolically active. Contains two fluorescent components, Calcofluor White and ViaVacTM Red/Green.



Product attributes

Call us: 800-304-5357

Email: btinfo@biotium.com

Product Description

The Yeast Vitality Staining Kit is a fluorogenic assay that allows you to determine whether your yeast cells are metabolically active. The kit contains two fluorescent components: Calcofluor White and ViaVacTM Red/Green.

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

How It Works

ViaVacTM Red/Green is a fluorescent viability dye for yeast, identical in structure to the dye FUN® 1. ViaVac passively diffuses into cells, showing green cytoplasmic staining. In many species of yeast and fungi, in metabolically active cells the dye is actively transported into the vacuole, resulting in red fluorescent staining of tubular intravacuolar structures.

Calcofluor White is a fluorescent blue dye that binds to cellulose and chitin, which can be found in the cell walls of fungi, algae, and plants.

After staining yeast with this kit, both living and dead cells will be outlined in blue, while only living cells will show red fluorescent, vacuolar tubules.

FUN 1 is a registered trademark of Thermo Fisher Scientific

This datasheet was generated on November 5, 2025 at 07:08:33 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/yeast-vitality-staining-kit/