Glowing products for science

ViaVac™ Red/Green (FUN®1)

A fluorescent vitality dye for yeast, identical in structure to the dye FUN® 1. Living yeast cells show red fluorescent vacuolar tubules.



Product attributes

Excitation/Emission	485/530 nm (cytoplasm), 485/620 nm (vacoule)
Molecular weight	528.84
Storage Conditions	Store at -10 to -35 °C, Protect from light

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

Product Description

ViaVacTM Red/Green (also known as FUN®1) is a fluorescent vitality dye for yeast. It forms red fluorescent vacuolar tubules in metabolically active yeast cells.

- Yeast vitality dye
- Red fluorescent tubules only in live cells
- Diffuse green fluorescence in all cells
- 10 mM solution in DMSO
- Store at -20°C, protected from light
- C₂₄H₁₈ClIN₂S

ViaVacTM dye passively enters cells, giving diffuse, green cytoplasmic staining of both live and dead cells. In metabolically active yeast and fungus cells the dye is actively transported into the vacuole, resulting in red fluorescent staining of tubular intravacuolar structures. ViaVacTM dye is identical in structure to the dye FUN®1.

FUN 1 is a registered trademark of Thermo Fisher Scientific

This datasheet was generated on November 6, 2025 at 12:20:05 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/viavac-redgreen/