



Glowing products for science

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

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 (vacuole)
---------------------	--

Molecular weight of antigen	528.84
-----------------------------	--------

Storage Conditions	Store at -10 to -35 °C, Protect from light
--------------------	--

Product Description

ViaVac™ 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
- $C_{24}H_{18}ClIN_2S$

ViaVac™ 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. ViaVac™ 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 July 20, 2026 at 07:58:26 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/viavac-redgreen/>