



Glowing products for science

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

Calcofluor White Stain, 5 mM in Water

A fluorescent blue dye that binds to cellulose and chitin, which can be found in the cell walls of fungi, algae, and plants. In budding yeast Calcofluor White strongly stains bud scars.



Product attributes

Excitation/Emission	360/430 nm
Molecular weight	916.98
Storage Conditions	Store at -10 to -35 °C, Protect from light

Product Description

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. In budding yeast Calcofluor White stains bud scars more strongly, because they have a higher concentration of chitin. Calcofluor White is sometimes used to identify fungal infections in medical samples.

- Fluorescent blue dye
- Binds to cellulose and chitin
- Used to stain cell walls of fungi, algae, and plants
- 5 mM solution in water
- $C_{40}H_{44}N_{12}O_{10}S_2$

Calcofluor White also stains gram-positive bacteria, but not gram-negative bacteria or mammalian cells. See our [Cellular Stains Table](#) for more information on how our dyes stain various organisms.

This datasheet was generated on January 9, 2026 at 06:43:10 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/calcofluor-white-stain-5-mm-water/>