



Glowing products for science

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

CF® Dye Dextran 3,500 MW, Anionic and Fixable

CF® Dye labeled dextrans could be used as a fluorescent fluid-phase markers to study cell permeability, endocytosis, or mechanisms of biomolecular delivery. The dextran is 3,500 MW, and contains a formaldehyde-fixable free-amine group.



Product attributes

Probe cellular localization	Fluid phase tracer
Cell permeability	Membrane impermeant
Colors	Blue

Product Description

CF® Dye labeled dextrans could be used as a fluorescent fluid-phase markers to study cell permeability, endocytosis, or mechanisms of biomolecular delivery. The dextran is 3,500 MW, and contains a formaldehyde-fixable free-amine group.

CF® Dye Dextran 3,500 MW, Anionic and Fixable is available in CF®350.

Note: Conjugates of blue-fluorescent dyes like CF®350, CF®405S and CF®405M are not recommended for detecting low abundance targets and may be challenging to use in tissue specimens. Blue dyes have lower fluorescence and photostability, and cells and tissue have high autofluorescence in blue wavelengths, resulting in lower signal to noise compared to other colors.

[Learn more about CF®350 features.](#)

This datasheet was generated on January 23, 2026 at 11:05:10 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cf-dye-dextran-3500-mw-anionic-and-fixable/>