



Glowing products for science

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

CF® Dye Dextran 70,000 MW, Anionic and Fixable

This dextran is 70,000 MW, and contains a formaldehyde-fixable free-amine group; available with a selection of CF® Dyes.



Product attributes

Probe cellular localization	Fluid phase tracer
Cell permeability	Membrane impermeant
Colors	Blue, Green, Far-red, Near-infrared

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 70,000 MW, and contains a formaldehyde-fixable free-amine group.

It's available with the following CF® Dyes: CF@350, CF@488A, CF@633, CF@680, CF@750 and CF@770.

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.

CF® Dye Dextran 70,000 MW

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF@350	347/448 nm	1 mg	80138	CF@350 Features
CF@488A	490/515 nm	1 mg	80117	CF@488A Features
CF@633	630/650 nm	1 mg	80141	CF@633 Features
CF@680	681/698 nm	1 mg	80129	CF@680 Features
CF@750	755/777 nm	1 mg	80130	CF@750 Features
CF@770	770/797 nm	1 mg	80123	CF@770 Features

This datasheet was generated on June 22, 2026 at 06:46:58 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cf-dye-dextran-70000-mw-anionic-and-fixable/>