



Glowing products for science

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

CF® Dye Dextran 250,000 MW, Anionic and Fixable

This dextran is 250,000 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	Green, 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 250,000 MW, and contains a formaldehyde-fixable free-amine group.

It's available with the following CF® dyes: CF®488A, CF®680, CF®750 and CF®770.

CF® Dye Dextran 250,000 MW

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF@488A	490/515 nm	1 mg	80134	CF@488A Features
CF@680	681/698 nm	1 mg	80135	CF@680 Features
CF@750	755/777 nm	1 mg	80136	CF@750 Features
CF@770	770/797 nm	1 mg	80125	CF@770 Features

This datasheet was generated on June 10, 2026 at 05:21:53 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cf-dye-dextran-250000-mw-anionic-and-fixable/>