



Glowing products for science

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

CF® Dye Dextran 150,000 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 150,000 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 150,000 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 150,000 MW

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF®488A	490/515 nm	1 mg	80131	CF®488A Features
CF®680	681/698 nm	1 mg	80132	CF®680 Features
CF®750	755/777 nm	1 mg	80133	CF®750 Features
CF®770	770/797 nm	1 mg	80124	CF®770 Features

This datasheet was generated on December 20, 2025 at 06:30:47 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cf-dye-dextran-150000-mw-anionic-and-fixable/>