



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
<a href="#">CF®488A</a>	490/515 nm	1 mg	<a href="#">80134</a>	<a href="#">CF®488A Features</a>
<a href="#">CF®680</a>	681/698 nm	1 mg	<a href="#">80135</a>	<a href="#">CF®680 Features</a>
<a href="#">CF®750</a>	755/777 nm	1 mg	<a href="#">80136</a>	<a href="#">CF®750 Features</a>
<a href="#">CF®770</a>	770/797 nm	1 mg	<a href="#">80125</a>	<a href="#">CF®770 Features</a>

This datasheet was generated on January 20, 2026 at 05:19:29 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/>