## CF® Dye Dextran 40,000 MW, Anionic and Fixable

This dextran is 40,000 MW, and contains a formaldehyde-fixable free-amine group; available with a selection of CF® dyes.



## **Product attributes**

Conjugation	CF®488A, CF®680, CF®750, CF®770
Probe cellular localization	Fluid phase tracer
Cell permeability	Membrane impermeant
Colors	Green Near-infrared

Call us: 800-304-5357 Email: btinfo@biotium.com

## **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 40,000 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 40,000 MW

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF®488A	490/515 nm	1 mg	<u>80126</u>	CF®488A Features
CF®680	681/698 nm	1 mg	<u>80127</u>	CF®680 Features
CF®750	755/777 nm	1 mg	<u>80128</u>	CF®750 Features
<u>CF®770</u>	770/797 nm	1 mg	<u>80122</u>	CF®770 Features

This datasheet was generated on November 7, 2025 at 07:11:07 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/cf-dye-dextran-40000-mw-anionic-and-fixable/">https://biotium.com/product/cf-dye-dextran-40000-mw-anionic-and-fixable/</a>