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

CF® Dye Biotin

Potential applications of Biotin-CF® Dye include detection of biotin binding sites and the degree of biotinylation of proteins, and for the measurement of avidin and streptavidin in crude biological samples. In addition, biotin-CF® Dye can be used as a polar tracer by microinjection to study the morphology of cells.



Product attributes

Product Description

Potential applications of Biotin-CF® Dye include detection of biotin binding sites and the degree of biotinylation of proteins, and for the measurement of avidin and streptavidin in crude biological samples. In addition, biotin-CF® Dye can be used as a polar tracer by microinjection to study the morphology of cells. CF® Dyes are very photostable which make them suitable for studies where prolonged exposure to light may be necessary. CF® Dye Biotin is available with a selection of 8 CF® Dyes.

Lear more about our bright and photostable <u>CF® Dyes</u>, as well as other available <u>CF® Dye bioconjugates</u>.

Biotin CF® Dye Conjugates

Conjugation	Ex/Em	Size	MW	Catalog No.	Dye Features
CF®488A	490/516 nm	1 mg	~1182	<u>80034</u>	CF®488A Features
<u>CF®532</u>	531/552 nm	1 mg	~954	<u>80038</u>	CF®532 Features
CF®568	562/583 nm	1 mg	~981	80029	CF®568 Features
CF®594	593/615 nm	1 mg	~997	<u>80035</u>	CF®594 Features
CF®633	630/650 nm	1 mg	~1089	<u>80031</u>	CF®633 Features
<u>CF®640R</u>	642/662 nm	1 mg	~1100	<u>80032</u>	CF®640R Features
CF®647	652/668 nm	1 mg	~1254	<u>80036</u>	CF®647 Features
<u>CF®680R</u>	680/701 nm	1 mg	~1180	<u>80037</u>	CF®680R Features

This datasheet was generated on November 12, 2025 at 12:36:36 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/cf-dye-biotin/