



Glowing products for science

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

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 [CF® Dyes](#), as well as other available [CF® Dye bioconjugates](#).

Biotin CF® Dye Conjugates

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

This datasheet was generated on January 13, 2026 at 11:37:08 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cf-dye-biotin/>