

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.

Learn 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 June 17, 2026 at 03:50:39 AM. Visit product page to check for updated information before use.

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