

CF® Dye Free Acids

Non-reactive, free acid forms of CF® Dyes that can be used as controls for the corresponding CF® Dye conjugates.



Product attributes

Functional group	Free acid
Storage Conditions	Store at -10 to -35 °C, Protect from light
Shelf life	Guaranteed for at least 12 months from date of receipt when stored as recommended
Reconstitution	Reconstitute in DMSO. The stock solution is stable for at least 12 months when stored at -10 to -35 °C

Product Description

CF® Dyes are a series of highly water-soluble fluorescent dyes spanning the visible and near-infrared spectrum. They offer superior brightness, photostability, and signal-to-noise when compared to other widely used fluorescent dyes for life science research. These free acid forms of CF® Dyes are non-reactive molecules that can be used as reference standards or controls for corresponding CF® Dye conjugates such as antibodies or other proteins for cell-based or *in vivo* assays.

- Non-reactive, free acid forms of the CF® Dyes
- Used as controls for corresponding CF® Dye conjugates

Learn more about Biotium's superior [CF® Dyes](#).

CF® Dye Free Acids

Dye	Ex/Em	MW	ϵ^1	Size	Catalog No.	Dye Features
CF@405M	416/452 nm	~503	41,000	1 mg	92116	CF@405M Features
CF@450	448/533 nm	~689	40,000	1 mg	98004	CF@450 Features
CF@488A	490/516 nm	~914	70,000	1 mg	92067	CF@488A Features
CF@532	531/552 nm	~685	96,000	1 mg	92072	CF@532 Features
CF@594	593/615 nm	~729	115,000	1 mg	96013	CF@594 Features
CF@640R	642/663 nm	~832	105,000	1 mg	92063	CF@640R Features
CF@647	652/668 nm	~985	240,000	1 mg	92062	CF@647 Features
CF@680	681/698 nm	~3153	210,000	1 mg	92070	CF@680 Features
CF@750	755/779 nm	~2921	250,000	1 mg	92071	CF@750 Features
CF@770	770/797 nm	~3091	220,000	1 mg	92066	CF@770 Features
CF@790	783/808 nm	~3179	210,000	0.25 umol	92159	CF@790 Features
CF@820	822/835 nm	~2711	253,000	1 mg	98023	CF@820 Features

1. Extinction Coefficient (ϵ).

References

Download a list of [CF® dye references](#).

This datasheet was generated on December 17, 2025 at 07:13:49 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/near-ir-cf-dye-free-acids/>