

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

1. Extinction Coefficient ( $\epsilon$ ).

## References

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

This datasheet was generated on August 2, 2025 at 04:43:08 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/near-ir-cf-dye-free-acids/>