

# CF™660C and CF™660R Dyes

## Superior alternatives to Alexa Fluor® 660

### Technical Summary

#### CF™660C

**Abs/Em Maxima:** 667/685 nm

**Extinction coefficient:** 200,000

**Molecular weight:** 3112

**Flow cytometry laser line:** 633 or 635 nm

**Microscopy laser line:** 633 or 635 nm

**Direct replacement for:** Alexa Fluor® 660, Allophycocyanin (APC)

#### CF™660R

**Abs/Em Maxima:** 663/682 nm

**Extinction coefficient:** 100,000

**Molecular weight:** 888

**Flow cytometry laser line:** 633 or 635 nm

**Microscopy laser line:** 633 or 635 nm

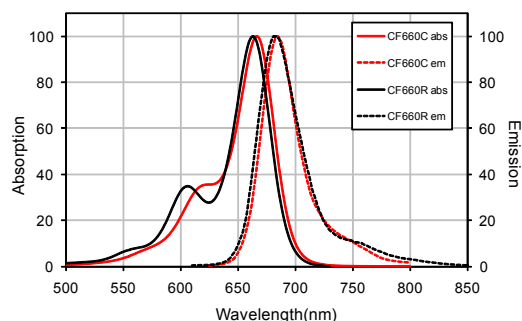
**Direct replacement for:** Alexa Fluor® 660, Allophycocyanin (APC)

### Advantages of CF™660C:

- Much brighter than Alexa Fluor® 660
- More photostable than Alexa Fluor® 660
- Highly water-soluble

### Advantages of CF™660R:

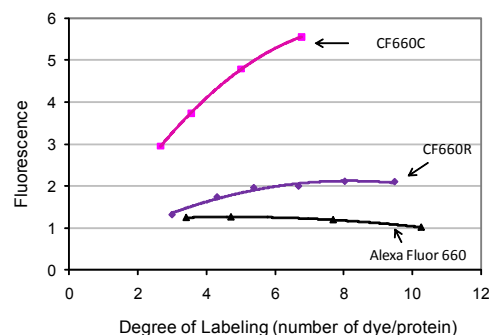
- Brighter than Alexa Fluor® 660
- The most photostable 660 nm dye ideal for confocal microscopy application
- Highly water-soluble



**Figure 1.** Absorption and emission spectra of CF™660C and CF™660R conjugated to goat anti-mouse IgG, respectively, in PBS.

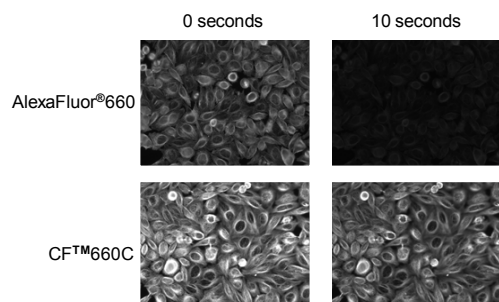
CF™660C and CF™660R are two spectrally similar fluorescent dyes that emit fluorescence at about 685 nm in the borderline spectral region between far-red and near-IR (Figure 1). Although their absorption maxima are at around 660 nm, both dyes can be sufficiently excited by the 633 or 635 nm laser. When combined with other CF™ dyes of shorter wavelengths, CF™660C or CF™660R can serve as a useful long wavelength dye in multi-color detection applications. The two dyes are spectrally similar to Alexa Fluor® 660 but are far superior to the latter in performance. Like Alexa Fluor® 660, CF™660C is a cyanine-based dye. However, when conjugated to protein, CF™660C is several fold brighter and significantly more photostable than Alexa Fluor® 660 (Figures 2 and 3). CF™660R is a rhodamine-based dye. Like rhodamine dyes in general, CF™660R is exceptionally photostable, compared to both Alexa Fluor® 660 and CF™660C (Figure 2). CF™660R is also much brighter than Alexa Fluor® 660, though not as bright as CF™660C (Figure 2). The superior photostability and excellent brightness of CF™660R make the dye an ideal choice for confocal microscopy and other demanding applications.

A list of CF™660C- and CF™660R-based products are shown in Table 1. A full selection of secondary antibodies, antibody labeling kits, and other bioconjugates including phalloidin, annexin V and  $\alpha$ -bungarotoxin are also available for many CF™ dyes. Please visit the Biotium website at [www.biotium.com](http://www.biotium.com) for details.

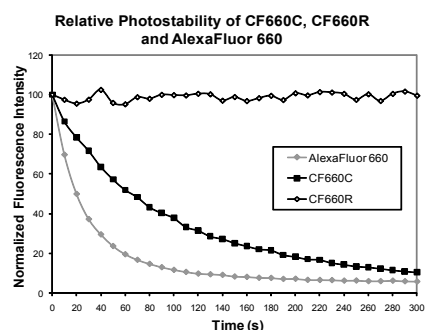


**Figure 2.** Relative fluorescence of CF™660C-, CF™660R and Alexa Fluor® 660-goat anti-mouse conjugates as a function of number of dye per protein (*i.e.*, degree of labeling). Fluorescence was measured at each dye's emission maximum in PBS using 633 nm excitation.

## CF™660C and CF™660R fluorescent reagents



**Figure 3.** HeLa cells were fixed, permeabilized and stained with mouse  $\alpha$ -tubulin followed by CF™660C or AlexaFluor® 660 goat anti-mouse IgG conjugates. Cells were imaged using an Olympus mercury arc lamp microscope equipped with a Cy5 filter set and CCD camera. Images were taken at t=0 and 10 seconds.



**Figure 4.** HeLa cells were fixed, permeabilized and stained with mouse  $\alpha$ -tubulin followed by CF™660C, CF™660R or AlexaFluor® 660 goat anti-mouse IgG conjugates. Cells were imaged using an Olympus mercury arc lamp microscope equipped with a Cy5 filter set and CCD camera. The graph illustrates the relative fluorescence intensities of sequential images taken every 10 seconds for 5 minutes.

Table 1. CF™660C and CF™660R Product List

Product Name	Size	Cat No.
<b>CF™660C- and CF™660R-Labeled Secondary Antibody Conjugates</b>		
CF™660C Donkey Anti-Goat IgG (H+L) whole antibody, 2mg/mL (min X Chicken, Guinea Pig, Syrian Hamster, Horse, Human, Mouse, Rabbit, and Rat)	0.5 mL	20051
CF™660C Goat Anti-Mouse IgG (H+L) whole antibody, 2 mg/mL	0.5 mL	20050
CF™660C Goat Anti-Mouse IgG (H+L) whole antibody, 2 mg/mL (min x Human, Bovine, Horse, Rabbit, and Swine)	0.5 mL	20052
CF™660C Goat Anti-Rabbit IgG (H+L) whole antibody, 2 mg/mL	0.5 mL	20053
CF™660R Goat Anti-Mouse IgG (H+L) whole antibody, 2 mg/mL	0.5 mL	20054
CF™660R Goat Anti-Rabbit IgG (H+L) whole antibody, 2 mg/mL	0.5 mL	20055
<b>Other CF™660R-Labeled Products</b>		
CF™660R Phalloidin	300 U	00047
CF™660R Streptavidin	1 mg	29040
<b>CF™660C and CF™660R Reactive Dyes and Labeling Kits</b>		
CF™660C maleimide	1 $\mu$ mole	92028
CF™660C succinimidyl ester	1 $\mu$ mole	92137
CF™660C SE protein labeling kit	3 labelings (for 1 mg protein each)	92219
Mix-n-Stain™ CF™660C antibody labeling kit, 1x(50-100 $\mu$ g) labeling	1 labeling	92239
Mix-n-Stain™ CF™660C antibody labeling kit, 1x(20-50 $\mu$ g) labeling	1 labeling	92260
Mix-n-Stain™ CF™660C antibody labeling kit, 1x(5-20 $\mu$ g) labeling	1 labeling	92280
CF™660R, aminoxy	1 mg	92059
CF™660R maleimide	1 $\mu$ mole	92031
CF™660R succinimidyl ester	1 $\mu$ mole	92134
CF™660R SE protein labeling kit	3 labelings (for 1 mg protein each)	92223
Mix-n-Stain™ CF™660R antibody labeling kit, 1x(50-100 $\mu$ g) labeling	1 labeling	92243
Mix-n-Stain™ CF™660R antibody labeling kit, 1x(20-50 $\mu$ g) labeling	1 labeling	92261
Mix-n-Stain™ CF™660R antibody labeling kit, 1x(5-20 $\mu$ g) labeling	1 labeling	92281

Listed products are for research use only. Not for use in diagnostic or therapeutic procedures. CF is a trademark of Biotium; CF dye technologies are covered by pending US and international patents. Alexa Fluor is registered trademark of Invitrogen.



**Biotium, Inc.**  
3159 Corporate Place, Hayward, CA 94545  
www.biotium.com  
US orders: 1-800-304-5357