

CF™633 Dye

Truly the best dye for the 633/635 nm lasers

Technical Summary

Abs/Em Maxima: 630/650 nm

Extinction coefficient: 100,000

Molecular weight: ~820

Flow cytometry laser line: 633 or 635 nm

Microscopy laser line: 633 or 635 nm

Direct replacement for: Alexa Fluor® 633, Alexa Fluor® 647, Cy™5 and DyLight™ 649

Advantages

- Yields the brightest antibody conjugates among spectrally similar dyes when excited by the 633 nm He-Ne laser or the 635 nm red diode laser
- Far more photostable than Alexa Fluor® 647
- Highly water-soluble

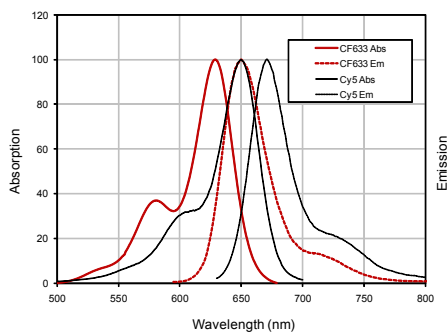


Figure 1. Absorption and emission spectra of goat anti-mouse IgG conjugates labeled with CF™633 and Cy™5, respectively, in PBS.

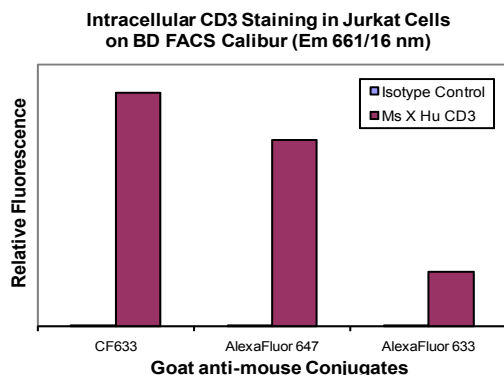


Figure 2. Jurkat cells were stained with intracellular CD3 or isotype control followed by goat anti-mouse IgG conjugates. Fluorescence was analyzed on a BD FACS Calibur in the FL4 channel. The bars represent the relative fluorescence of the geometric means of the population of cells.

Far-red fluorescent dyes offer the advantage of ultra sensitive detection because background signal due to autofluorescence in most biological samples is minimal in this spectral region. For many years, the cyanine dye Cy™5 has been the dye of choice for such detection. More recently, Alexa Fluor® 647 has been developed as a better alternative by having brighter fluorescence and higher photostability. Despite the improvement, Alexa Fluor® 647 still lacks sufficient photostability required for many demanding applications. On the other hand, while Alexa Fluor® 633 is photostable, its fluorescence on proteins is very weak. In fact, it has been a challenging task for dye chemists to develop a far-red fluorescent dye that is both highly fluorescent and photostable for protein and nucleic acid labeling. Using new chemistry, scientists at Biotium have successfully developed a series of far-red and near-IR dyes to overcome these challenges. With its absorption peak at 630 nm, CF™633 can be optimally excited by the 633 nm He-Ne laser or the 635 nm red diode laser with emission maximum at 650 nm. Despite its shorter emission wavelength, which often is not favored by a Cy™5-based detection window on most instruments, CF™633 is still significantly brighter than Alexa Fluor® 647 (Figure 2). However, for detections using non-laser-based excitation and a Cy™5 filter set, the full advantage of CF™633 will not be realized. In this case, we recommend CF™640R, another outstanding far-red CF™ dye that is spectrally closer to Cy™5 or other similar cyanine-based dyes. The most important advantage of CF™633 and other rhodamine-based far-red CF™ dyes are their unmatched photostability (Figure 3). The combination of superior brightness and photostability makes far-red CF™ dyes ideal choices for detections in the long wavelength region on a variety of instruments.

A list of CF™633-based products are shown in Table 1. A full selection of secondary antibodies, antibody labeling kits, and other bioconjugates including phalloidin, annexin V and α -bungarotoxin are also available for many CF™ dyes. Please visit the Biotium website at www.biotium.com for details.

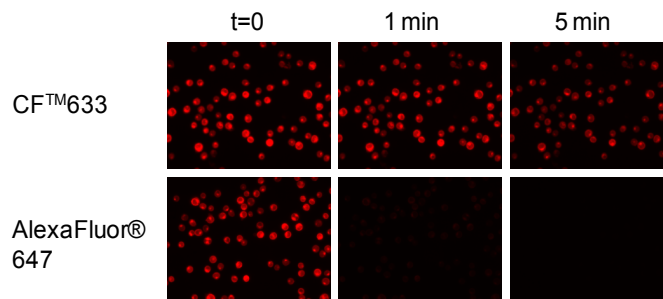


Figure 3. Jurkat cells were fixed, permeabilized and stained with rabbit anti-CD3 (Abcam) followed by CF™633 (Biotium) or AlexaFluor® 647 (Invitrogen) goat anti-rabbit IgG conjugates. Cells were imaged using an Olympus mercury arc lamp microscope equipped with a Cy5 filter set and CCD camera. Images were taken at t=0, 1 min and 5 min.

CF™633 fluorescent reagents

Table 1. CF™633 Product List

Product Name	Size	Cat No.
CF™633-Labeled Secondary Antibody Conjugates		
Donkey Anti-Goat IgG (H+L) whole antibody, 2 mg/mL (min X Chicken, Guinea Pig, Syrian Hamster, Horse, Human, Mouse, Rabbit, and Rat)	0.5 mL	20127
Donkey Anti-Mouse IgG (H+L) whole antibody, 2 mg/mL (min X Bovine, Chicken, Goat, Guinea Pig, Syrian Hamster, Horse, Human, Rabbit, and Sheep)	0.5 mL	20124
Donkey Anti-Rabbit IgG (H+L) whole antibody, 2 mg/mL (min X Bovine, Chicken, Goat, Guinea Pig, Syrian Hamster, Horse, Human, Mouse, Rat, and Sheep)	0.5 mL	20125
Donkey Anti-Rat IgG (H+L) whole antibody, 2 mg/mL (min X Bovine, Chicken, Goat, Guinea Pig, Syrian Hamster, Horse, Human, Rabbit, and Sheep)	0.5 mL	20137
Donkey Anti-Sheep IgG (H+L) whole antibody, 2 mg/mL (min X Chicken, Guinea Pig, Syrian Hamster, Horse, Human, Mouse, Rabbit, and Rat)	0.5 mL	20134
Goat Anti-Chicken IgY (IgG) (H+L) whole antibody, 2 mg/mL (min X Bovine, Goat, Guinea Pig, Syrian Hamster, Horse, Human, Mouse, Rabbit, Rat, and Sheep)	0.5 mL	20126
Goat Anti-Guinea Pig IgG (H+L) whole antibody, 2 mg/mL	0.5 mL	20129
Goat Anti-Human IgG (H+L) whole antibody, 2 mg/mL (min X Bovine, Horse, and Mouse)	0.5 mL	20132
Goat Anti-Mouse IgG (H+L) whole antibody, 2 mg/mL	0.5 mL	20120
Goat Anti-Mouse IgG (H+L) whole antibody, 2 mg/mL (min x Human, Bovine, Horse, Rabbit, and Swine)	0.5 mL	20121
Goat Anti-Mouse IgG (H+L), F(ab')₂ fragment, 2 mg/mL	0.25 mL	20130
Goat Anti-Rabbit IgG (H+L) whole antibody, 2 mg/mL	0.5 mL	20122
Goat Anti-Rabbit IgG (H+L) whole antibody, 2 mg/mL (min X Human, Mouse, and Rat)	0.5 mL	20123
Goat Anti-Rabbit IgG (H+L), F(ab')₂ fragment, 2 mg/mL	0.25 mL	20131
Goat Anti-Rat IgG (H+L) whole antibody, 2 mg/mL (min X Human, Bovine, Horse, and Rabbit)	0.5 mL	20133
Goat Anti-Swine IgG (H+L) whole antibody, 2 mg/mL	0.5 mL	20138
Rabbit Anti-Goat IgG (H+L) whole antibody, 2 mg/mL	0.5 mL	20128
Rabbit Anti-Mouse IgG (H+L) whole antibody, 2 mg/mL (min X Human)	0.5 mL	20136
Rabbit Anti-Rat IgG (H+L) whole antibody, 2 mg/mL (min X Human)	0.5 mL	20135
Other CF™633-Labeled Products		
Annexin V, 50 µg/mL	0.5 mL	29008
α-Bungarotoxin	0.5 mg	00009
Phalloidin	300 U	00046
Streptavidin	1 mg	29037
Wheat germ agglutinin (WGA)	5x1 mg	29024
CF™633 Reactive Dyes and Labeling Kits		
CF™633, aminoxy	1 mg	92053
CF™633 hydrazide	1 mg	92156
CF™633 maleimide	1 µmole	92026
CF™633 succinimidyl ester	1 µmole	92133
CF™633 SE protein labeling kit	3 labelings (for 1 mg protein each)	92217
Mix-n-Stain™ CF™633 antibody labeling kit, 1x(50-100 µg) labeling	1 labeling	92237
Mix-n-Stain™ CF™633 antibody labeling kit, 1x(20-50 µg) labeling	1 labeling	92257
Mix-n-Stain™ CF™633 antibody labeling kit, 1x(5-20 µg) labeling	1 labeling	92277

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. Cy is a trademark of GE Healthcare. DyLight is a trademark of ThermoFisher.



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