

CF™350 Dye

A superior UV-excitable blue fluorescent dye

Technical Summary

Abs/Em Maxima: 347/448 nm

Extinction coefficient: 18,000

Molecular weight: 496

Flow cytometry laser line: UV

Microscopy laser line: UV nm

Direct replacement for: Alexa Fluor® 350, AMCA

Advantages

- Yields the brightest blue fluorescent antibody conjugates when excited at ~350 nm
- Highly water-soluble
- Insensitive to pH

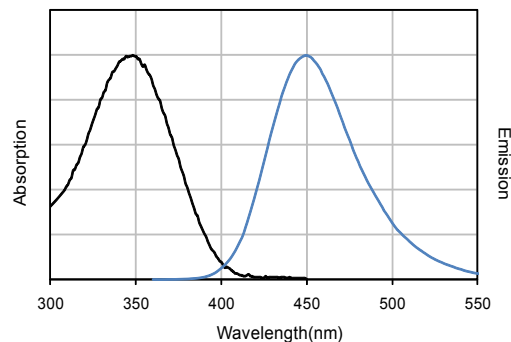


Figure 1. Absorption and emission spectra of CF™350 conjugated to goat anti-mouse IgG in PBS.

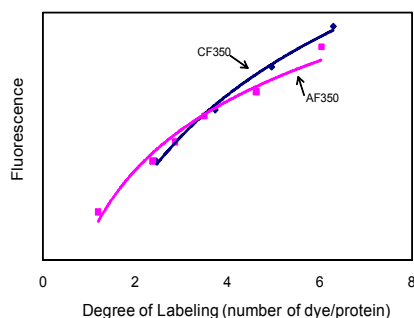


Figure 2. Relative fluorescence of CF™350- and Alexa Fluor® 350-goat anti-mouse conjugates as a function of number of dye per protein (*i.e.*, degree of labeling). Fluorescence was measured at each dye's emission maximum in PBS using 350 nm excitation.

CF™350 is a blue fluorescent dye spectrally similar to the traditional blue fluorescent dye, AMCA. However, compared to AMCA, CF™350 is more water-soluble, more photostable and at least 50% more fluorescent on proteins. CF™350 is at least as bright as Alexa fluor® 350 and thus can be a direct replacement for the latter for all applications (Figure 2).

A list of CF™350-based products are shown in Table 1. A full selection of secondary antibodies, antibody labeling kits, and other bioconjugates including phalloidins, annexin V and α -bungarotoxin are also available for many CF™ dyes. To label your own antibodies, see our revolutionary Mix-n-Stain™ antibody labeling kits, which allow you to prepare CF™ dye-labeled antibodies in 30 minutes without a separation step. Visit Biotium website: www.biotium.com for details.

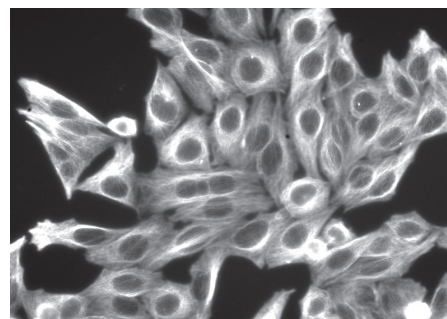


Figure 3. HeLa cells were fixed, permeabilized and stained with mouse anti-alpha-tubulin followed by CF350 goat anti-mouse IgG. Image was taken on an Olympus mercury arc lamp microscope.

If you are looking for an antibody conjugate not listed in our catalog, please let us know. We might be able to add it as a new product, or perform a custom conjugation for you.



Biotium, Inc.

3159 Corporate Place, Hayward, CA 94545

www.biotium.com

US orders: 1-800-304-5357

CF™350 fluorescent reagents

Table 1. CF™350 Product List

Product Name	Size	Cat No.
CF™350-Labeled Secondary Antibody Conjugates		
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	20142
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	20350
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	20351
Donkey Anti-Sheep IgG (H+L) whole antibody , 2mg/mL (min X Chicken, Guinea Pig, Syrian Hamster, Horse, Human, Mouse, Rabbit, and Rat)	0.5 mL	20148
Goat Anti-Mouse IgG (H+L) whole antibody , 2 mg/mL	0.5 mL	20140
Goat Anti-Mouse IgG (H+L) whole antibody , 2 mg/mL (min x Human, Bovine, Horse, Rabbit, and Swine)	0.5 mL	20143
F(ab')₂ fragment of Goat Anti-Mouse IgG (H+L) , 2 mg/mL	0.25 mL	20145
Goat Anti-Rabbit IgG (H+L) whole antibody , 2 mg/mL	0.5 mL	20141
Goat Anti-Rabbit IgG (H+L) whole antibody , 2 mg/mL (min X Human, Mouse, and Rat)	0.5 mL	20144
F(ab')₂ fragment of Goat Anti-Rabbit IgG (H+L) , 2 mg/mL	0.25 mL	20146
Goat Anti-Rat IgG (H+L) whole antibody , 2 mg/mL (min X Human, Bovine, Horse, and Rabbit)	0.5 mL	20147
Rabbit Anti-Mouse IgG (H+L) whole antibody , 2 mg/mL (min X Human)	0.5 mL	20149
Other CF™350-Labeled Products		
Annexin V , 50 µg/mL	0.5 mL	29012
Phalloidin	300 U	00049
Streptavidin	1 mg	29031
CF™350 Reactive Dyes and Labeling Kits		
CF™350 amine	1 mg	92035
CF™350, aminooxy	1 mg	92050
CF™350 hydrazide	1 mg	92151
CF™350 maleimide	1 µmole	92020
CF™350 succinimidyl ester	1 µmole	92109
CF™350 SE protein labeling kit	3 labelings (for 1 mg protein each)	92210
Mix-n-Stain™ CF™350 antibody labeling kit, 1x(50-100 µg) labeling	1 labeling	92230
Mix-n-Stain™ CF™350 antibody labeling kit, 1x(20-50 µg) labeling	1 labeling	92250
Mix-n-Stain™ CF™350 antibody labeling kit, 1x(5-20 µg) labeling	1 labeling	92270

If you are looking for an antibody conjugate not listed in our catalog, please let us know. We might be able to add it as a new product, or perform a custom conjugation for you.