

CF™555 Dye

An orange-red dye superior to Cy™3

Technical Summary

Abs/Em Maxima: 555/565 nm

Extinction coefficient: 150,000

Molecular weight: ~900

Flow cytometry laser line: 532 nm

Microscopy laser line: 543, 546, 555 nm

Direct replacement for: Alexa Fluor® 555, Cy™3, Tetramethylrhodamine.

Advantages

- Brighter and more photostable than Cy™3
- Highly water-soluble and pH-insensitive

CF™555 is an orange-red fluorescent dye spectrally similar to Cy™3. Like Cy™3, it is also a cyanine-based dye with excellent water solubility. However, CF™555 is brighter and more photostable than Cy™3 when conjugated to proteins. In general, the performance of CF™555 is comparable to that of Alexa Fluor™ 555, another Cy™3-like dye.

A list of CF™555-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. Visit Biotium website: www.biotium.com for details.

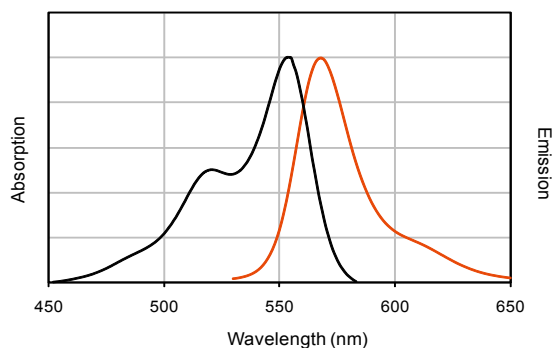


Figure 1. Absorption and emission spectra of CF™555 conjugated to goat anti-mouse IgG in PBS.

Figure 3.

Figure 2.



Biotium, Inc.

3159 Corporate Place, Hayward, CA 94545

www.biotium.com

US orders: 1-800-304-5357

CF™555 fluorescent reagents

Table 1. Partial List of CF™555 Products

| Product Name | Size | Cat No. |
|--|--|--------------|
| CF™555-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 | 20039 |
| 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 | 20037 |
| 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 | 20038 |
| 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 | 20234 |
| 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 | 20034 |
| Goat Anti-Guinea Pig IgG (H+L) whole antibody , 2 mg/mL | 0.5 mL | 20036 |
| Goat Anti-Human IgG (H+L) whole antibody , 2 mg/mL (min X Bovine, Horse, and Mouse) | 0.5 mL | 20230 |
| Goat Anti-Mouse IgG (H+L) whole antibody , 2 mg/mL | 0.5 mL | 20030 |
| Goat Anti-Mouse IgG (H+L) whole antibody , 2 mg/mL (min x Human, Bovine, Horse, Rabbit, and Swine) | 0.5 mL | 20231 |
| Goat Anti-Mouse IgG (H+L), F(ab')₂ fragment , 2 mg/mL | 0.25 mL | 20032 |
| Goat Anti-Rabbit IgG (H+L) whole antibody , 2 mg/mL | 0.5 mL | 20033 |
| Goat Anti-Rabbit IgG (H+L) whole antibody , 2 mg/mL (min X Human, Mouse, and Rat) | 0.5 mL | 20232 |
| Goat Anti-Rabbit IgG (H+L), F(ab')₂ fragment , 2 mg/mL | 0.25 | 20035 |
| Goat Anti-Rat IgG (H+L) whole antibody , 2 mg/mL (min X Human, Bovine, Horse, and Rabbit) | 0.5 mL | 20233 |
| Goat Anti-Swine IgG (H+L) whole antibody , 2 mg/mL | 0.5 mL | 20236 |
| Rabbit Anti-Goat IgG (H+L) whole antibody , 2 mg/mL | 0.5 mL | 20031 |
| Rabbit Anti-Mouse IgG (H+L) whole antibody , 2 mg/mL (min X Human) | 0.5 mL | 20235 |
| Other CF™555-Labeled Products | | |
| Annexin V, 50 µg/mL | 0.5 mL | 29004 |
| α-Bungarotoxin | 0.5 mg | 00018 |
| Phalloidin | 300 U | 00040 |
| Streptavidin | 1 mg | 29038 |
| CF™555 Reactive Dyes and Labeling Kits | | |
| CF™555 hydrazide | 1 mg | 92153 |
| CF™555 maleimide | 1 µmole | 92023 |
| CF™555 succinimidyl ester | 1 µmole | 92130 |
| CF™555 SE protein labeling kit | 3 labelings (for 1 mg protein each) | 92214 |
| Mix-n-Stain™ CF™555 antibody labeling kit | 1x (50-100 µg) labeling | 92234 |
| Mix-n-Stain™ CF™555 antibody labeling kit | 1x (20-50 µg) labeling | 92254 |
| Mix-n-Stain™ CF™555 antibody labeling kit | 1x (5-20 µg) labeling | 92274 |

*Please check Biotium website for updated lists of CF™555 and other CF™ dye products.

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



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