CF[®] Dyes

Next-Generation Fluorescent Dyes

| Introd | uction to CF® Dyes | |
|--------|---|----|
| | Quick reference table | 2 |
| | Technology overview | 3 |
| | Dye applications at a glance | 4 |
| | CF® Dyes FAQs | 5 |
| Techni | cal Information for CF® Dyes | |
| | Dye profiles by color | 6 |
| | Industry-leading near-infrared dyes | 19 |
| | Dyes for super-resolution imaging | 20 |
| Reacti | ve Dyes and Labeling Kits | |
| | Mix-n-Stain™ antibody labeling kits | 22 |
| | Protein & small ligand labeling kits | 23 |
| | Chemically reactive dyes | 24 |
| CF® Dy | ve Bioconjugates | |
| | Toxins, ligands, tracers, and lectins | 25 |
| | Labeled nucleotides | 25 |
| Antibo | ody Conjugates | |
| | Primary antibodies | 26 |
| | Anti-biotin, anti-GFP, anti-epitope tag Abs | 27 |
| | Secondary antibodies | 27 |
| | Highly cross-adsorbed secondaries | 28 |
| | F(ab') ₂ fragments | 29 |
| | Isotype-specific secondary antibodies | 29 |
| Relate | d Products & Accessories | |
| | Tyramide signal amplification | 30 |
| | Background suppressors & more | 31 |
| | | |



CF® Dyes Quick Reference Table

| | Dye | Ex/Em (nm) | Excitation | Alternative for | Specialized Applications | Brightness* | Photostability* |
|------------------|----------------|------------|------------------|---|---|-------------|-----------------|
| | CF®350 | 347/448 | UV | Alexa Fluor® 350, AMCA, DyLight® 350 | | • | • |
| | CF*405S | 411/431 | 405 nm | Alexa Fluor [®] 405, Cascade Blue™, DyLight [®] 405 | SIM | • | •• |
| | CF®405M | 416/452 | 405 nm | BD Horizon™ V450, eFluor® 450, Pacific Blue™ | SIM, STED, 2-Photon | • | •• |
| | CF®405L | 413/547 | 405 nm | Pacific Orange™, Qdot™ 565, Spark Violet™ 538 | | • | • |
| | CF*430 | 426/498 | 405 nm | Pacific Green™, BD Horizon™ V500, Krome Orange™ | | • | •• |
| E | CF®440 | 440/515 | 405 nm | Alexa Fluor® 430 | | • | •• |
| ectr | CF®450 | 450/538 | 405 nm | Unique dye | | • | • |
| e sp | CF®488A | 490/515 | 488 nm | ATTO 488, Alexa Fluor® 488, Spark Blue™ 515 Cy®2, DyLight® 488, FAM, FITC | SIM, STED, STORM, 2-Photon, TIRF, DNA-PAINT | ••• | ••• |
| Visible spectrum | CF®503R | 503/532 | 488 nm | ATTO 488 | | | |
| S | CF®505 | 505/519 | 488 nm | ATTO 488 | STORM | ••• | ••• |
| | CF*514 | 516/548 | 488 nm | Alexa Fluor [®] 514, Spark Blue [™] 550 | | • • | ••• |
| | CF®532 | 527/558 | 532 nm | Alexa Fluor® 532, ATTO 532, Qdot™ 565 | | | |
| | CF®535ST | 535/568 | 532 nm | Unique dye for STORM | STORM | | |
| | CF®543 | 541/560 | 532 to 546 nm | Alexa Fluor [®] 546, Tetramethylrhodamine (TAMRA) | | • • • | •• |
| | CF®550R | 551/577 | 532 to 568 nm | Unique dye | | ••• | ••• |
| | CF*555 | 555/565 | 532 to 568 nm | Alexa Fluor® 555, ATTO 550, Cy®3, DyLight® 549, TRITC | SIM, STORM | ••• | •• |
| | CF*568 | 562/583 | 532 to 568 nm | Alexa Fluor [®] 568, Spark YG™ 581, ATTO 565, Rhodamine Red | SIM, STED, STORM, TIRF | ••• | ••• |
| | CF*570 | 568/591 | 532 to 568 nm | Alexa Fluor [®] 568, Spark YG [™] 593, ATTO 565, DY-560, Rhodamine Red | | ••• | •• |
| | CF*583 | 583/606 | 532 to 568 nm | Cy®3.5, Texas Red® | | • • | •• |
| | CF®583R | 586/609 | 532 to 568 nm | Cy®3.5, Texas Red® | STORM | ••• | • |
| | CF*594 | 593/614 | 532 to 568 nm | Alexa Fluor® 594, ATTO 594, DyLight® 594, Texas Red® | STED, 2-Photon | | |
| ed | RPE-Astral™616 | 565/617 | 488 nm or 561 nm | PE-Texas Red®, PE/Dazzle™ 594, PE-CF®594 | | | • |
| Far-red | CF®597R | 597/619 | 561 to 568 nm | Alexa Fluor® 594, ATTO 594, DyLight® 594 | STORM | ••• | • |
| | CF*620R | 617/639 | 633 or 635 nm | LightCycler® Red 640 | | • • | ••• |
| | CF*633 | 630/650 | 633 or 635 nm | Alexa Fluor [®] 633, Alexa Fluor [®] 647, Cy [®] 5, DyLight [®] 633 | TIRF, FIONA, gSHRImP, SMT | ••• | ••• |
| | CF®640R | 642/662 | 633 to 640 nm | Alexa Fluor [®] 647, ATTO 647N, Cy [®] 5, DyLight [®] 649 | SIM, STED, TIRF, FLImP, 2-Photon | | ••• |
| | CF®647 | 652/668 | 633 to 640 nm | Alexa Fluor [®] 647, ATTO 647N, Cy [®] 5, DyLight [®] 649 | STORM | ••• | • |
| | CF®647Plus | 652/668 | 633 to 640 nm | Alexa Fluor [®] 647, ATTO 647N, Cy [®] 5, DyLight [®] 649 | | ••• | • |
| | CF*660C | 667/685 | 633 to 640 nm | Alexa Fluor [®] 660, Spark NIR™ 685 | STORM, MINFLUX | • • | •• |
| | CF*660R | 663/682 | 633 to 640 nm | Alexa Fluor® 660 | SMLM, DNA-PAINT | • • | ••• |
| | CF*680 | 681/698 | 680 or 685 nm | Alexa Fluor [®] 680, Cy [®] 5.5, DyLight [®] 680, IRDye [®] 680LT | Near-IR western, STORM, 3D SMLM, MINFLUX | ••• | •• |
| | CF®680R | 680/701 | 680 or 685 nm | Alexa Fluor® 680, Cy®5.5, DyLight® 680, IRDye® 680LT | STED, STORM, SMT, 2-Photon, single molecule spectroscopy | •• | ••• |
| | CF*700 | 695/720 | 680 or 685 nm | Alexa Fluor [®] 700, DyLight [®] 700, BD Horizon [™] Red 718, Spark Red [™] 718 | | • • | •• |
| arec | CF*750 | 755/777 | 680 or 685 nm | Alexa Fluor® 750, Cy®7, DyLight® 750, IRDye® 750 | Photoacoustic imaging, STORM | • • | • |
| infr | CF*770 | 770/797 | 785 nm | DyLight® 800, IRDye® 800CW, ZW800-1 | Near-IR western | •• | • |
| Near-infrared | CF*790 | 784/806 | 785 nm | Alexa Fluor® 790 | | •• | • |
| Z | APC-Astral™813 | 788/813 | 633 to 640 nm | APC/Fire™ 810 | | ••• | • |
| | CF*800 | 797/816 | 785 nm | Spectrally similar to Indocyanine green | | •• | • |
| | CF*820 | 822/835 | 785 nm | DY-820 | | • • | • |
| | CF*850 | 852/570 | 808 nm | Unique dye | | • • | • |
| | CF*870 | 876/896 | 808 nm | Unique dye | | •• | • |

FLIMP: Fluorophore localization imaging with photobleaching; SIM: Structured illumination microscopy; STED: Stimulated emission depletion; STORM: Stochastical optical reconstruction microscopy; TIRF: Total internal reflection fluorescence; FIONA: Fluorescence imaging with one-nanometer accuracy; ExM: Expansion microscopy; SMT: Single-molecule tracking; SMLM: Single-molecule localization microscopy.

The relative brightness and photostability of CF Dyes shown in this table are intended as a general guideline. The values are partially based on extinction coefficients and dye structure, as well as our experience with antibody conjugates in immunofluorescence and flow cytometry experiments. Many factors, such as degree of labeling (DOL), laser power, filters, and gain, influence the performance of fluorescent dyes on a given instrument.

CF® Dyes Technology Overview

Next-Generation Fluorescent Dyes

CF® Dyes are a series of highly water-soluble fluorescent dyes spanning the visible and near-infrared (IR) spectrum for labeling biomolecules, especially proteins and nucleic acids. Developed by scientists at Biotium using new breakthrough chemistries, CF® Dyes rival or exceed the quality of other commercial dyes, such as Alexa Fluor® dyes, due to several novel features.

Novel Rhodamine Chemistry

Rhodamine dyes are known for their excellent photostability and good fluorescence quantum yield; consequently several of the Alexa Fluor® dyes bear the rhodamine core structure. Unfortunately, traditional rhodamine chemistry makes it difficult to extend the fluorescence wavelength into the far-red region and even more challenging to extend into the near-IR region; especially for water-soluble dyes designed for bioconjugation. Recently, Biotium scientists discovered a new way to prepare novel rhodamine dyes of any fluorescence color from green to near-IR. The new chemistry is key to overcoming these challenges and lead to the development of many of our CF Dyes. The new chemistry is a key element in the development of many of our CF® Dyes, which are not only bright and water-soluble but also extremely photostable.

Excellent Labeling Efficiency

Reactive dyes for bioconjugation are generally susceptible to hydrolysis, which can cause problems for shipping, handling and storage, and result in lower labeling efficiency. Heavily sulfonated dyes, such as the Alexa Fluor®, IRDye®, and DyLight® dyes are particularly hygroscopic, worsening the hydrolysis problem. For example, the percent of active Alexa Fluor® 488 succinimidyl ester (SE) could be well below 50% by the time of application (according to the Alexa Fluor® 488 Microscale Labeling Kit product information sheet, provided by Thermo Fisher Scientific). In contrast, all of Biotium's amine-reactive CF® Dyes have a relatively stable form of SE, which is more resistant to hydrolysis than the SE on many of the Alexa Fluor® dyes. Accordingly, CF® Dye SE products generally give consistently higher labeling efficiency, thus providing users better results at a better value.

Mix-n-Stain™ Antibody Labeling Technology

Biotium has developed a breakthrough antibody labeling technology with CF® Dye Mix-n-Stain™ antibody labeling kits. With this technology, you merely need to mix your antibody with the reaction buffer and the CF® Dye provided in the kit. In 30 minutes, you will have an optimally labeled CF® Dye-antibody conjugate ready for immunostaining. The labeling technology provides unprecedented convenience for antibody labeling. Mix-n-Stain™ labeled antibodies can be used for multicolor immunostaining, allowing staining with multiple primary antibodies from the same host species when pre-labeled primary antibodies are not available.

Unrivaled Near-Infrared Dyes

Near-IR dyes are typically much larger in size than dyes in the visible range. The large size often results in serious problems of low dye solubility, dye aggregation/quenching, and poor fluorescence quantum yield. To overcome the problems, many commercial near-IR dyes, such as the near-IR Alexa Fluor®, IRDye®, and DyLight® dyes, are prepared by placing a number of negatively charged sulfonate group on the dyes. While sulfonation improves dye solubility and fluorescence quantum yield to some degree, it creates another even more serious problem: non-specific binding of the bioconjugates prepared from the dyes. For example, conjugation to a highly negatively charged dye can dramatically alter an antibody's isoelectric point, which is essential for maintaining specific antibody-antigen interaction (for examples, see page 19, Figure 3 and page 31, Figure 2).

With this insight, Biotium scientists devised a revolutionary new approach to near-IR dye design using our patented polyethylene glycol dye modification, or pegylated dye chemistry. Dye pegylation offers several key benefits for dye performance:

- · Increases dye solubility without adding charges
- Shields any existing charges on the dye
- Reduces dye aggregation and self-quenching on conjugates for brighter fluorescence
- · Increases both thermal and photostability of the dye
- Perfectly suited for *in vivo* imaging; pegylated dye modification is known to reduce protein immunogenicity and improve biocompatibility

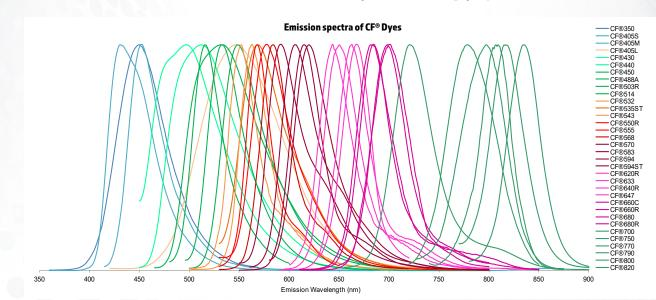
These features, along with a large and growing selection of available wavelengths, make CF® Dyes the industry leaders in near-infrared dye technology. See pages 18-19 to learn more about near-infrared CF® Dyes.

CF® Dyes for Super-Resolution Microscopy

Recent publications comparing synthetic dyes for super-resolution imaging have shown CF® Dyes give the best performance for multiple methods. The superior brightness, photostability, and photochemical switching properties of certain CF® Dyes are ideal for 3-D SIM, 3-D STORM, and other super-resolution and single-molecule imaging techniques. See page 21 for more information.

Multicolor Flexibility

Biotium currently offers more than 30 CF® Dyes, with additional colors in development. The CF® Dye product line includes reactive dyes with a full selection of functional groups (page 24), easy-to-use labeling kits (page 23), CF® Dye-labeled primary and secondary antibodies (pages 26-29), and many other CF® Dye conjugates such as toxins, tracers, ligands, and nucleotides (page 25).

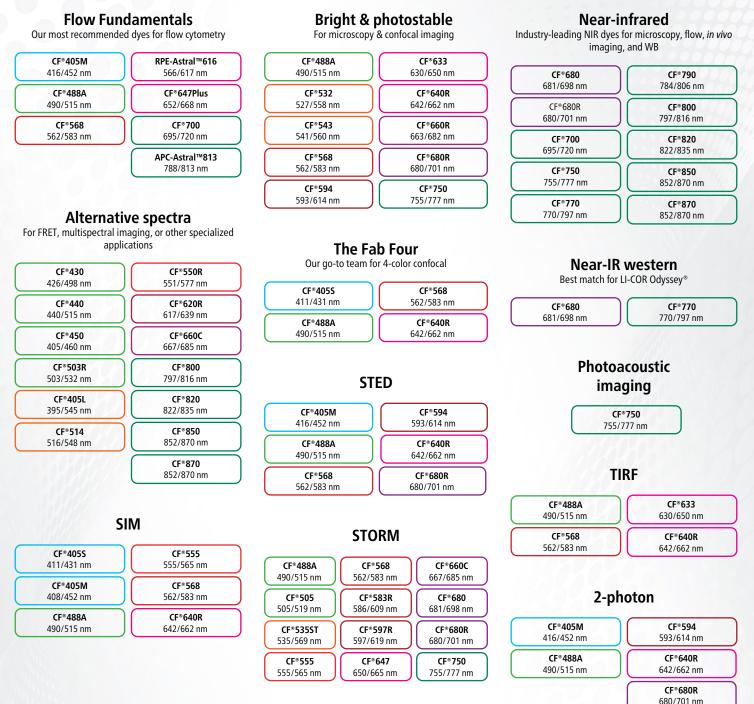


CF® Dyes and conjugates have been cited in hundreds of publications, with new articles published every day. Visit www.biotium.com to download a list of selected references. See page 21 for dyes validated for super-resolution and 2-photon imaging.

Alexa Fluor, Cascade Blue, DyLight, Pacific Blue, Pacific Green, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; ATTO dyes are products of ATTO-TEC GmbH; BD Horizon is a trademark of BD Bioscience; (R) is a registered trademark of GE Healthcare; eFluor is a registered trademark of eBioscience; IRDye and Odyssey are registered trademarks of L-COR Bioscience; Krome Orange is a trademark of Beckman Coulter; LightCycler is a registered trademark of Roche Applied Science.

Dyes At a Glance: Select the Right Dye for Your Application

Use our CF[®] Dye Selection Tool at www.biotium.com to find the best CF[®] Dyes for your application and instrument configuration.



CF[®] Dyes are being tested in new applications all the time, visit **biotium.com** for the most up-to-date information.

Alexa Fluor, Cascade Blue, Pacific Blue, DyLight, and Texas Red are registered trademarks of Thermo Fisher Scientific; Qdot is a trademark of Thermo Fisher Scientific; ATTO dyes are products of ATTO-TEC GmbH; BD Horizon is a trademark of BD Biosciences; Cy is a registered trademark of Cytiva; eFluor is a registered trademark of eBioscience; IRDye and Odyssey are registered trademarks of LI-COR Bioscience; APC/Fire, PE/Dazzle, Spark Blue, Spark Red, Spark YG, and Spark NIR are trademarks of Biolegend; Krome Orange is a trademark of Beckman Coulter; LightCycler is a registered trademark of Roche Applied Science.

See page 20-21 for more information on CF® Dyes for super-resolution imaging and other specialized applications. CF® Dyes are being tested in new applications all the time, visit biotium.com for the most up-to-date information.

Frequently Asked Questions (FAQs)

| Question | Answer |
|---|--|
| What does the CF in CF® Dyes stand for? | CF® was initially an abbreviation for "Cyanine-based Fluorescent dyes". These were the first patented CF® Dyes based on cyanine dye structures. 10 years and more than two dozen dyes later, the CF® Dye portfolio encompasses multiple dye core structures spanning the fluorescence spectrum from UV to near-IR. Today, we believe "CF" more aptly stands for Clear Fluor: dyes that produce superior signal-to-noise. |
| What are pegylated dyes? | Dye pegylation is one of Biotium's patented dye technologies that improves solubility and brightness of near-infrared dyes without introducing excess negative charge, making our near-IR CF® Dyes industry leaders. See page 3 for more information on pegylated dye technology; see pages 18-19 to learn more about near-IR CF® Dyes. |
| What are the chemical structures of CF® Dyes? | The exact chemical structures of CF® Dyes are currently confidential but will be fully disclosed at a later stage when pending patents become granted. In general terms, the structure of a CF® Dye may be divided into two parts: a) dye core structure (i.e. the aromatic ring skeleton that defines the dye's color or absorption/emission wavelengths), and b) core structure-modifying elements. At present, CF® Dyes bear the core structures of coumarin, pyrene, rhodamine, or cyanine dyes. Blue fluorescent CF® Dyes are based on a coumarin or pyrene dye core structure, while green to near-IR CF® Dyes are based on either cyanine or rhodamine dye core structures. Core structure-modifying elements refer to various chemical attachments to the core structure and are a key aspect of the CF® Dye invention that makes CF® Dyes superior to other commercial dyes. |
| What are the quantum yields of CF® Dyes? | The quantum yield of a fluorescent dye can vary widely depending on the dye's micro-environment and if the dye is attached to a protein or other molecule. A good way to compare the relative quantum yields of different dyes is to plot the total fluorescence of the labeled proteins as a function of degree of labeling by the dyes, as we have done with CF® Dyes and other commercial dyes in the dye description pages in this guide. |
| How stable are CF® Dyes? | There are three aspects to dye stability: Chemical stability: CF® Dyes bear the core structures of coumarin, pyrene, rhodamine, or cyanine dyes; all of which are known to have excellent chemical stability. In general, the dyes are far more stable than the antibodies or other biomolecules they label. CF® Dyes are also stable enough for labeled nucleic acids to be used in PCR or nucleic acid hybridization, where high temperature is involved. Reactive group stability: Reactive CF Dyes have a chemically reactive group for bioconjugation. Among the various reactive groups, only amine-reactive succinimidyl ester (SE) and thiol-reactive maleimide groups are susceptible to hydrolysis and therefore are moisture-sensitive. CF® Dye SE products are relatively more stable than other SE dyes. This is because CF® SE Dyes are derived from aliphatic carboxylic groups, which results in a more stable SE form, while other SE dyes usually are derived from aromatic carboxylic acid groups that yield a less stable SE form. Photostability: This refers to the dye's ability to withstand photobleaching. Photobleaching is mainly a concern when dyes are subjected to intense illumination for an extended period of time, such as during confocal microscopy. Among the four types of core structures, rhodamine is the most photostable, followed by cyanine, pyrene, and coumarin cores. The structure-modifying groups and the way they are attached to the dye cores are a key innovative aspect of CF® Dye technologies that contributes to the superior photostability of CF® Dyes over that of other dyes. In general, rhodamine-based CF® Dyes, whose wavelengths range from green to the near-IR region, offer the best photostability, making these dyes ideal for microscopy applications. |
| Are CF® Dyes sensitive to pH? | CF® Dyes are chemically stable within the range of at least pH 2 to pH 11. The fluorescence of most CF® Dyes is relatively insensitive to pH, except for that of CF®405M, CF®568, CF®620R, and CF®633. The fluorescence of these four CF® Dyes becomes weaker when pH drops below 4.5. |
| Are CF® Dyes fixable? | CF® Dyes can tolerate formaldehyde fixation. However, whether a CF® Dye-labeled probe is fixable will depend on the fixability of the probe itself. Proteins with free amine groups that bind other proteins generally are formaldehyde-fixable. |
| What is the difference between CF®405S, CF®405M, and CF®405L? | All three of these dyes can be excited by the 405 nm laser (or UV mercury lamp). They differ in their emission wavelengths. CF®405S has the shortest blue fluorescence emission at 431 nm, while CF®405M has a longer wavelength blue fluorescence emission at 452 nm. CF®405L has orange fluorescence emission at 545 nm. We recommend choosing the dye that best fits your instrument's detection settings (see pages 6-7 for more information). |
| For several CF® Dye colors, there is an R form and a C form, both having similar absorption and emission spectra. In such a case, which of the two CF® Dyes should I choose? | Rhodamine-based CF® Dyes (designated R) generally have better photostability but weaker fluorescence than their cyanine-based equivalents (designated C). Therefore, rhodamine-based near-IR CF® Dyes are a better choice for microscopy, while cyanine-based CF® Dyes are more ideal for flow cytometry, western blotting, and other applications where photobleaching is less of a concern. Another factor to consider is the size of the dyes. Some of the cyanine-based near-IR CF® Dyes are much larger than the rhodamine-based equivalents. For antibody labeling, either version of the CF® Dyes is suitable. However, for applications where the dye size may cause a steric problem, the smaller dye may be a better choice. |
| How soluble are CF® Dyes? | CF® Dyes are highly water-soluble (>100 mg/mL). They are also very soluble in other polar solvents, such as DMSO, DMF, methanol, and ethanol. However, CF® Dyes are poorly soluble or insoluble in non-polar solvents. |
| What are the charges on CF® Dyes? | Most CF® Dyes carry 1-2 negative charges while some cyanine-based near-IR CF® Dyes carry 3-4 negative charges. However, the more negatively charged CF® Dyes have unique structural features that shield the biomolecules from the negative charges; such that the biomolecules (such as antibodies) do not lose specificity due to excess negative charge. |
| Can CF® Dyes be used for STORM, STED, SIM, or TIRF? | Many of our CF® Dyes have been validated in multiple super-resolution techniques. Biotium also offers dyes specifically designed for STORM imaging. See pages 20-21 for more information. |
| Can CF® Dyes be used for 2-photon microscopy? | Some CF® Dyes have been validated for 2-photon excitation. See pages 20-21 for more information. |
| What are the major applications of CF® Dyes? | CF® Dyes are ideal for protein labeling because of their high water solubility, which reduces fluorescence quenching. They are also useful for labeling oligonucleotides that require multiple copies of a dye for maximal fluorescence, such as the preparation of FISH probes, where water-soluble dyes can minimize fluorescence quenching. Finally, CF® Dyes make excellent polar tracers that can be used for visualizing the morphology or long-term tracing of neurons. Several CF® Dyes have been validated in specialized applications, including spectral flow cytometry, SIM, TIRF, STORM, and other super-resolution imaging techniques, as well as photoacoustic imaging and 2-photon microscopy. See pages 20-21 for more information about CF® Dyes in super-resolution and other specialized imaging applications. |
| | |

CF® 350 A bright UV-excitable blue fluorescent dye

Technical Summary

Abs/Em maxima: 347/448 nm

- Extinction coefficient: 18,000
- Molecular weight: ~ 496
- Excitation source: UV

Alternative for: Alexa Fluor® 350, AMCA, DyLight® 350

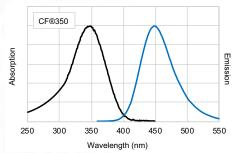


Figure 1. Absorption and emission spectra of CF \circledast 350 goat anti-mouse conjugate in PBS.

CF®405S and CF®405M

Improved brightness and photostability for the 405 nm laser line

Technical Summary

CF®4055

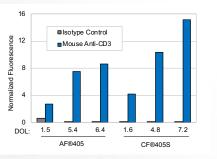
F[®]405S & CF[®]405M

Abs/Em maxima: 404/431 nm Extinction coefficient: 33,000 Molecular weight: ~ 1,169 Excitation laser line: 405 nm Alternative for: Alexa Fluor® 405, Cascade Blue™, DyLight® 405

CF®405M

Abs/Em maxima: 408/452 nm Extinction coefficient: 41,000 Molecular weight: ~ 503 Excitation laser line: 405 nm Alternative for: Pacific Blue™, BD Horizon™ V450

Figure 2. Intracellular staining of Jurkat cells was performed with mouse anti-CD3 or isotype control followed by goat anti-mouse IgG conjugated to Alexa Fluor® 405 (AF405) or CF®405S. Fluorescence was analyzed on a BD LSRII flow cytometer with 405 nm excitation and 450/50 nm emission filter. Bars represent geometric mean fluorescence.



Features

- Brighter and more photostable than AMCA
- Direct replacement for Alexa Fluor® 350
- Highly water-soluble and pH-insensitive

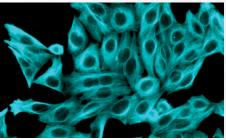


Figure 2. HeLa cells stained with mouse anti-tubulin antibody and CF®350 goat anti-mouse IgG (cyan).

Features

- CF®405S: Brighter than Alexa Fluor® 405 (Fig. 2)
- CF®405M: More photostable than Pacific Blue™, with less spillover in the green channel
- Validated for super-resolution imaging by SIM (see p. 21)

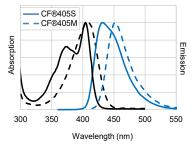


Figure 1. Absorption and emission spectra of CF®405S and CF®405M goat anti-mouse conjugates in PBS.

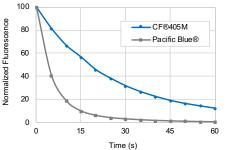


Figure 3. Relative photostability of CF®405M and Pacific Blue™. CF®405M and Pacific Blue™ dye solutions were continuously exposed to mercury arc lamp microscope excitation with a DAPI filter set. Images were captured every 5 seconds for one minute. Fluorescence intensity was normalized to time 0.

5 • www.biotium.com

CF®405L

A 405 nm-excitable dye with orange fluorescence emission

Technical Summary

Abs/Em maxima: 395/545 nm Extinction coefficient: 24,000 Molecular weight: ~ 1573 Excitation laser line: 405 nm Alternative for: Pacific Orange™

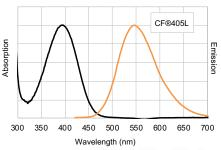


Figure 1. Absorption and emission spectra of CF®405L goat anti-mouse conjugate in PBS.

CF®430 and CF®440

Photostable 405 nm-excitable dyes with green fluorescence

Technical Summary

CF®430

Abs/Em maxima: 426/498 nm Extinction coefficient: 40,000 Molecular weight: ~ 429 Excitation laser line: 405 nm Alternative for: Pacific Green™, BD Horizon™ V500, Krome Orange™

CF®440

Abs/Em maxima: 440/515 nm Extinction coefficient: 40,000 Molecular weight: ~ 716 Excitation laser line: 405 nm Alternative for: Alexa Fluor® 430

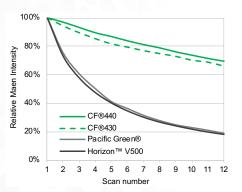


Figure 2. Relative photostability of CF®430 and CF®440 compared to spectrally-similar dyes. Cells were stained with biotinylated primary antibodies followed by streptavidin conjugates of CF®430, CF®440, Pacific Green™, or BD Horizon™ V500. Fluorescence was imaged on a Zeiss LSM700 confocal microscope in the FITC channel using 405 nm excitation. Images were acquired every 5 seconds for 12 consecutive scans of the same field of view using the same imaging settings for each dye. The mean fluorescence intensity of each image was normalized to the first scan for each dye.

Features

- Photostable dyes suitable for microscopy
- CF®430 is a perfect match for the CFP filter set
- Suitable for flow cytometry in the AmCyan channel
- Highly water-soluble and pH-insensitive

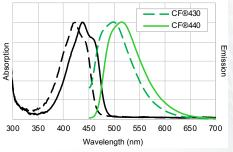


Figure 1. Absorption and emission spectra of CF®430 and CF®440 goat anti-mouse conjugates in PBS.

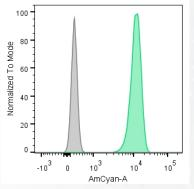


Figure 3. Flow cytometry analysis of Jurkat cells stained with isotype control (gray peak) or mouse anti-CD3 (green peak) followed by CF®430 goat anti-mouse IgG, analyzed in the AmCyan channel of a BD LSRII flow cytometer. F[®]488A

CF®450 405 nm-excitable green dye with unique spectral properties

Technical Summary

Abs/Em maxima: 448/533 nm Extinction coefficient: 40,000 Molecular weight: ~ 689 Excitation laser line: 405 nm

CF®488A

A superior green fluorescent dye

Technical Summary

Abs/Em maxima: 490/515 nm

Extinction coefficient: 70,000

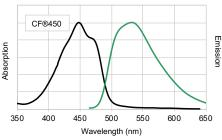
Molecular weight: ~ 914

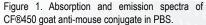
Excitation laser line: 488 nm

Alternative for: Alexa Fluor® 488, DyLight® 488, fluorescein (FITC, FAM), Cy®2

Features

- Minimally charged, for less non-specific binding than Alexa Fluor® 488
- Narrower emission spectrum for less bleed into the red channel
- Very photostable
- Compatible with STED, TIRF, and 2-photon microscopy (p. 21)





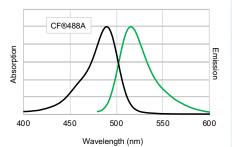


Figure 1. Absorption and emission spectra of CF®488A goat anti-mouse conjugate in PBS.

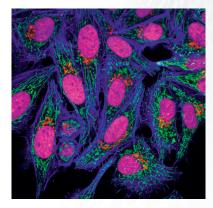


Figure 2. HeLa cells stained with rabbit anti-COXIV and CF®488A goat anti-rabbit IgG (mitochondria, green), mouse anti-Golgin 97 and CF®555 goat anti-mouse IgG (Golgi, red), CF®405M phalloidin (actin filaments, blue), and RedDot™2 (nuclei, magenta). See p. 31 for more information on RedDot™2.

CF®503R & CF®514

Alternative green fluorescent dyes for spectral imaging

Technical Summary CF®503R

Abs/Em maxima: 503/532 nm Extinction coefficient: 90,000 Molecular weight: ~ 1100 Excitation laser line: 488 nm

Technical Summary CF®514

Abs/Em maxima: 516/548 nm Extinction coefficient: 105,000 Molecular weight: ~ 1216 Excitation laser line: 488 nm Alternative for: Alexa Fluor® 514

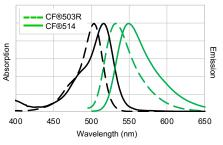


Figure 1. Absorption and emission spectra of CF®503R or CF®514 goat anti-mouse conjugates in PBS.

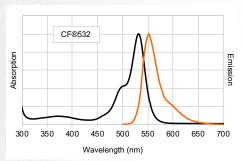
www.biotium.com • 9

CF®535S1

Technical Summary

CF®532

Abs/Em maxima: 527/558 nm Extinction coefficient: 96,000 Molecular weight: ~ 685 Excitation laser line: 532 nm Alternative for: Alexa Fluor® 532, Atto 532



A bright orange fluorescent dye for the 532 nm laser

Figure 1. Absorption and emission spectra of CF®532 goat anti-mouse IgG conjugate in PBS.

Features

- Designed for the 532 nm laser
- Brighter than Alexa Fluor® 532 (Fig. 2)
- Highly water-soluble and pH-insensitive

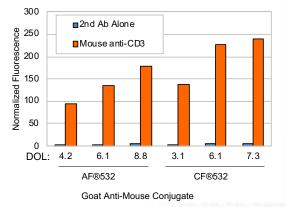


Figure 2. Flow cytometry analysis of Jurkat cells stained with Alexa Fluor® 532 (AF532) antibody or CF®532 secondary antibody conjugates. Intracellular staining was performed with mouse anti-CD3 antibody followed by goat anti-mouse IgG conjugates. Background was determined by staining with secondary antibody (2nd Ab) alone. Fluorescence was analyzed on a BD FACSCalibur flow cytometer in the FL2 channel. The bars represent the relative fluorescence of the geometric means of the cell populations.

CF[®]535ST

An orange fluorescent dye designed for STORM super-resolution imaging

Technical Summary

Abs/Em maxima: 535/568 nm Extinction coefficient: 95,000 Molecular weight: ~ 728 Excitation laser line: 532 nm

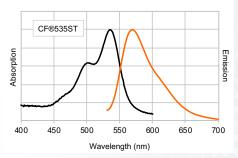


Figure 1. Absorption and emission spectra of CF®535ST goat anti-mouse IgG conjugate in PBS.

See page 21 for more information about CF® Dyes for super-resolution imaging.

CF®543 Bright orange dye ideal for the 543 nm laser

Technical Summary

Abs/Em maxima: 541/560 nm Extinction coefficient: 100,000 Molecular weight: ~ 870 Excitation laser line: 532 to 546 nm Alternative for: Alexa Fluor® 546, TAMRA

Features

- Optimized for the 543 nm laser
- Brightest conjugates among similar dyes
- Highly water-soluble and pH-insensitive

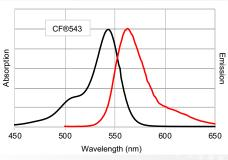


Figure 1. Absorption and emission spectra of CF\$543 goat anti-mouse conjugate in PBS.

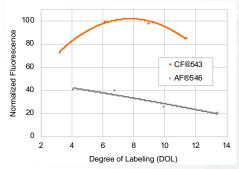


Figure 2. Relative fluorescence of CF®543 and Alexa Fluor® 546 (AF546) goat anti-mouse conjugates as a function of the number of dye molecules per protein (degree of labeling).

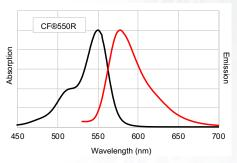
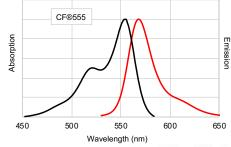


Figure 1. Absorption and emission spectra of CF\$550R goat anti-mouse conjugate in PBS.



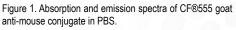


Figure 2. Rat intestine stained with CF®555 Mix-n-Stain™ labeled mouse anti-ZO1 (tight junctions, red) and NucSpot® 470 (nuclei, green). See p. 22 for more information on Mix-n-Stain™ kits.

CF®550R

Alternative orange/red dye for spectral imaging

Technical Summary

Abs/Em maxima: 551/577 nm Extinction coefficient: 100,000 Molecular weight: ~ 686 Excitation laser line: 532 nm or 568 nm

CF®555

A bright and photostable orange-red dye

Technical Summary

Abs/Em maxima: 555/565 nm

Extinction coefficient: 150,000

Molecular weight: ~ 959

Excitation laser line: 532 nm or 568 nm

Alternative for: Alexa Fluor® 555, ATTO 550, Cy®3, DyLight® 549, Rhodamine

Features

- Brighter than Cy®3
- Highly water-soluble
- Validated in multicolor STORM (see p. 21)

CF[®]555

CF®568 Outshines Alexa Fluor®568

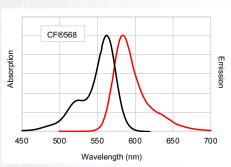
Technical Summary

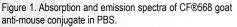
Abs/Em maxima: 562/583 nm

Extinction coefficient: 100,000

Molecular weight: ~ 714

- Excitation laser line: 532 nm or 568 nm
- Alternative for: Alexa Fluor® 568, ATTO 565, Rhodamine Red





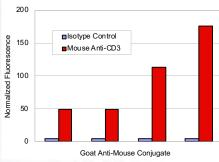


Figure 2. Intracellular staining of Jurkat cells was performed using mouse anti-CD3 or isotype control followed by goat anti-mouse IgG conjugates. Fluorescence was analyzed on a BD FACSCalibur flow cytometer in the FL2 channel. Bars represent the relative fluorescence of the geometric means of the cell populations.

Features

- Much brighter antibody conjugates than Alexa Fluor® 568
- Extremely photostable
- Excellent choice for multiplexing with CF®488A and CF®640R
- Compatible with TIRF and multicolor STORM (see p. 21)

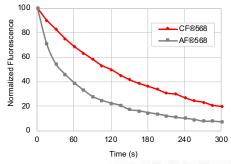


Figure 3. Photostability of CF®568 and Alexa Fluor® 568 (AF568) streptavidin conjugates. Intracellular staining of Jurkat cells was performed using anti-CD3-biotin followed by streptavidin conjugated to CF®568 or AF568. Cells were continuously exposed to mercury arc lamp microscope excitation with a Cy®3 filter set. Images were captured every 15 seconds for 5 minutes and fluorescence intensity was normalized to time 0.

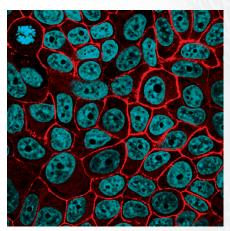


Figure 4. MCF-7 cells stained with CF®568 monoclonal anti-Ep-CAM (clone EGP40/826) at 5 ug/mL (red). Nuclei are counterstained with Hoechst 33342 (blue). See p. 26 for more information on primary antibody conjugates.

CF®570 Red fluorescent dye with superior brightness

Technical Summary

Abs/Em maxima: 568/591 nm

Extinction coefficient: 150,000

Molecular weight: ~ 2998

Excitation laser line: 561 to 568 nm

Alternative for: Alexa Fluor® 568, ATTO 565, DY-560, Rhodamine Red

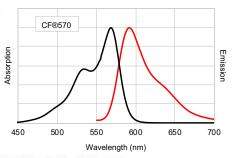


Figure 1. Absorption and emission spectra of CF®570 goat anti-mouse conjugate in PBS.

CF®583 & CF®583R

Brighter than Cy®3.5

CF®583 Technical Summary

Abs/Em maxima: 583/606 nm Extinction coefficient: 150,000 Molecular weight: ~ 3127 Excitation laser line: 561 to 568 nm Alternative for: Cy®3.5

CF®583R Technical Summary

Abs/Em maxima: 586/609 nm Extinction coefficient: 100,000 Molecular weight: ~ 773 Excitation laser line: 561 to 568 nm Alternative for: Cy®3.5, Texas Red®

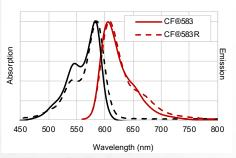


Figure 1. Absorption and emission spectra of CF®583 & CF®583R

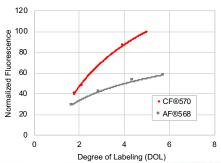


Figure 2. Relative fluorescence of CF®570 and Alexa Fluor® 568 (AF568) goat anti-mouse conjugates as a function of the number of dye molecules per protein (degree of labeling).

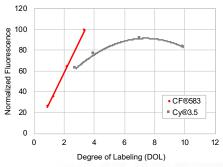


Figure 2. Relative fluorescence of CF®583 and Cy®3.5 goat anti-mouse conjugates as a function of the number of dye molecules per protein (degree of labeling).

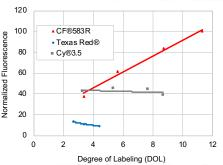


Figure 3. CF®583R produces brighter conjugates at a lower degree of labeling than Cy®3.5 and Texas Red®. Relative fluorescence of goat anti-mouse conjugates of the indicated dyes at varying degrees of labeling (DOL, or dye molecules per antibody).

CF[®]594 & CF[®]594ST

Truly the brightest deep red dye, with STORM-compatible option

Technical Summary

Abs/Em maxima: 593/614 nm Extinction coefficient: 115,000 Molecular weight: ~ 729 Excitation laser line: 532 to 594 nm Alternative for: Alexa Fluor® 594, DyLight® 594, Texas Red®

Features

- Brightest antibody conjugates among spectrally similar dyes, with excellent photostability
- Compatible with 2-photon microscopy (see p. 21)
- CF®594ST matches CF®594 spectrally, but is compatible with STORM imaging (see p. 21)

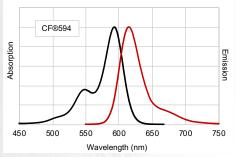


Figure 1. Absorption and emission spectra of CF@594 goat anti-mouse conjugate in PBS.

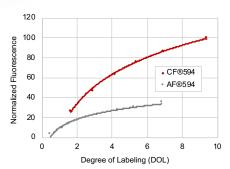


Figure 2. Relative fluorescence of CF®594 and Alexa Fluor® 594 (AF594) goat anti-mouse conjugates as a function of the number of dye molecules per protein (degree of labeling).

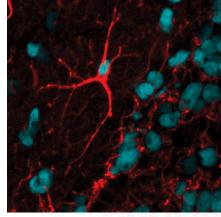


Figure 3. Glial cells in frozen section of rat brain stained with rabbit anti-GFAP antibody and CF®594 goat antirabbit IgG (red). Nuclei are stained with RedDot™2 (pseudocolored cyan). Mounted with Everbrite™ Mounting Medium. See p. 31 for more information on RedDot™2 and EverBrite™ Mounting Medium.

CF®620R A bright and photostable far-red dye

Technical Summary Abs/Em maxima: 617/639 nm Extinction coefficient: 115,000 Molecular weight: ~ 738 Excitation laser line: 633 nm or 635 nm Spectrally similar to: LightCycler® Red 640

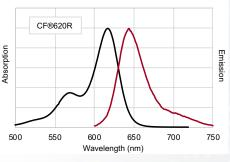


Figure 1. Absorption and emission spectra of CF®620R free acid in PBS.

- Features
- Highly fluorescent and extremely photostable
- Absorption/emission at 617/639 nm for use in FRET or other specialized applications

CF®620F

CF®633 The best dye for 633/635 laser lines

Technical Summary

Abs/Em maxima: 630/650 nm

Extinction coefficient: 100,000

Molecular weight: ~ 821

Excitation laser line: 633 nm or 635 nm

Alternative for: Alexa Fluor® 633, Alexa Fluor® 647, Cy®5, DyLight® 633, DyLight® 649

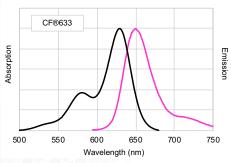


Figure 1. Absorption and emission spectra of CF@633 goat anti-mouse conjugate in PBS.

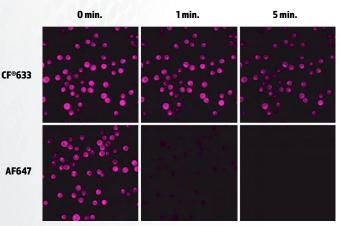


Figure 3. Relative photostability of CF®633 and Alexa Fluor® 647 (AF647) goat anti-mouse conjugates. Jurkat cells were fixed, permeabilized, and stained with rabbit anti-CD3 followed by CF®633 or Alexa Fluor® 647 goat anti-rabbit IgG conjugates. Cells were imaged using a mercury arc lamp microscope equipped with a Cy®5 filter set and CCD camera. Sequential images were captured at 0, 1, and 5 minutes.

Features

- Yields the brightest antibody conjugates among spectrally similar dyes
- Far more photostable than Alexa Fluor® 647
- Compatible with TIRF, FIONA, and gSHRImP super-resolution imaging methods (see p. 21)

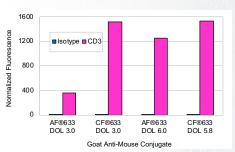


Figure 2. CF®633 yields the brightest far-red conjugates. Jurkat cells were stained with mouse anti-CD3 or isotype control antibody, followed by goat anti-mouse conjugates with varying degrees of labeling (DOL, or dye molecules per antibody). Fluorescence was measured in the APC channel of a BD FACSCalibur™ flow cytometer; bars represent geometric mean fluorescence.

CF®640R

A highly photostable far-red dye for the 640 nm laser

Technical Summary

Abs/Em maxima: 642/662 nm Extinction coefficient: 105,000

- Molecular weight: ~ 832
- Excitation laser line: 633 to 640 nm
- Alternative for: Alexa Fluor® 647, ATTO 647N, Cy®5, DyLight® 649

Figure 1. Absorption and emission spectra of CF®640R goat anti-mouse conjugate in PBS.

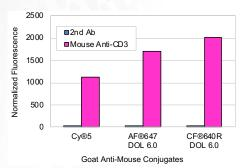


Figure 2. Intracellular staining of Jurkat cells was performed using mouse anti-CD3 or isotype control followed by goat anti-mouse IgG conjugates. Fluorescence was analyzed on a BD FACSCalibur flow cytometer in the FL4 channel. Bars represent the relative fluorescence of the geometric means of the cell populations.

Features

- Best photostability among Cy®5-like dyes
- Yields highly fluorescent protein conjugates
- Compatible with TIRF and FLImP super-resolution microscopy (see p. 21)

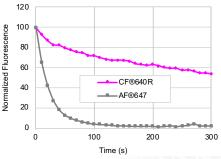


Figure 3. Relative photostability of CF®640R and Alexa Fluor® 647 (AF647). HeLa cells were stained with anti-tubulin antibody conjugates of each dye. Cells were continuously illuminated by a mercury arc lamp microscope and sequential images were captured at 0, 1, and 3 minutes. Mean fluorescence was normalized to time 0.

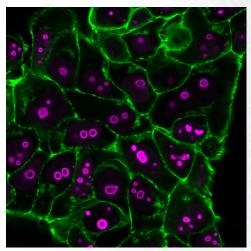
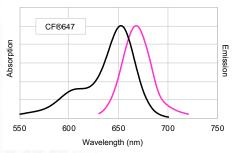


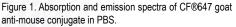
Figure 4. MCF-7 cells stained with CF®640R monoclonal anti-Cyclin B1 (clone CCNB1/1098) at 5 ug/mL (magenta). Actin filaments are stained with CF®488A phalloidin (green). See p. 26 for more information on primary antibody conjugates.

CF®647 A highly fluorescent far-red dye

Technical Summary

Abs/Em maxima: 650/665 nm Extinction coefficient: 240,000 Molecular weight: ~ 1058 Excitation laser line: 633 to 640 nm Alternative for: Cy®5, Alexa Fluor® 647, DyLight® 649





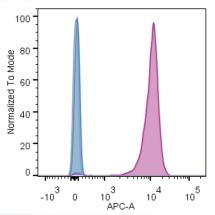


Figure 2. Intracellular staining of Jurkat cells with CF®647 monoclonal anti-nucleolin (clone 365-2) (pink) or CF®647 IgG1 isotype control (blue) at 1 ug/tube, compared to unstained cells (yellow). Cells were analyzed in the APC channel of a BD LSRII flow cytometer. See p. 26 for more information on primary antibody conjugates.

Features

- Brighter than Cy®5
- Highly water-soluble and pH-insensitive
- Validated in multi-color STORM imaging (see p. 21)

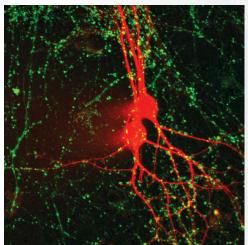


Figure 3. Cultured rat hippocampal neurons microinjected with CF®647 hydrazide (red) and stained with SynaptoGreen™ C4 (FM®1-43) (green, synaptic vesicles). Image courtesy of Professor Guosong Liu, Tsinghua University, Beijing, China.

CF®660C & CF®660R

CF®660C and CF®660R

Superior alternatives to Alexa Fluor® 660

Technical Summary

CF®660C

Abs/Em maxima: 667/685 nm Extinction coefficient: 200,000 Molecular weight: ~ 3112 Excitation laser line: 633 to 640 nm Alternative for: Alexa Fluor® 660, APC

CF®660R

Abs/Em maxima: 663/682 nm Extinction coefficient: 100,000 Molecular weight: ~ 888 Excitation laser line: 633 to 640 nm Alternative for: Alexa Fluor® 660, APC

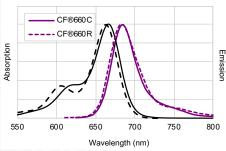


Figure 1. Absorption and emission spectra of CF@660C and CF@660R goat anti-mouse conjugates in PBS.

CF®660C Features

- Much brighter and more photostable than Alexa Fluor® 660
- Compatible with multicolor super-resolution imaging by STORM (see p. 21)

CF®660R Features

- Brighter than Alexa Fluor® 660
- Unrivaled photostability among spectrally similar dyes

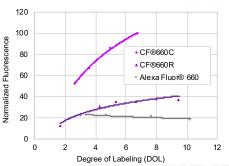


Figure 2. Relative fluorescence of CF®660, CF®660R, and Alexa Fluor® 660 (AF660) goat anti-mouse conjugates as a function of the number of dye molecules per protein (degree of labeling).

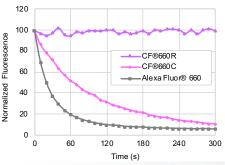


Figure 3. Relative photostability of CF®660C, CF®660R, and Alexa Fluor® 660 (AF660) conjugates. HeLa cells were stained with mouse anti-tubulin followed by CF®660C, CF®660R or AF660 goat anti-mouse IgG conjugates. Cells were continuously exposed to mercury arc lamp microscope excitation using a Cy®5 filter set. Images were captured every 10 seconds for five minutes and fluorescence intensity was normalized to time 0.

CF®680 and CF®680R

Two outstanding 680 nm-excitable dyes

Technical Summary

CF®680

Abs/Em maxima: 681/698 nm Extinction coefficient: 210,000 Molecular weight: ~ 3241 Excitation laser line: 680 nm or 685 nm Alternative for: Alexa Fluor® 680, Cy®5.5, IRDye® 680

CF®680R

Abs/Em maxima: 680/701 nm Extinction coefficient: 140,000 Molecular weight: ~ 912 Excitation laser line: 680 nm or 685 nm Alternative for: Alexa Fluor® 680, Cy®5.5, IRDye® 680

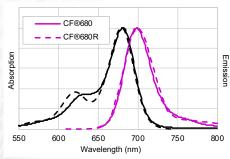


Figure 1. Absorption and emission spectra of CF®680 and CF®680R goat anti-mouse conjugates in PBS.

CF®680 Features

- The brightest among spectrally similar dyes
- Validated in multicolor STORM and 3D super-resolution microscopy (see p. 21)
- Compatible with LI-COR® Odyssey®

CF®680R Features

- Unrivaled photostability among spectrally similar dyes
- Compatible with STED, STORM, single molecule spectroscopy, and 2-photon microscopy (see p. 21)
- Molecular weight compatible with nucleic acid labeling
- Compatible with LI-COR® Odyssey®

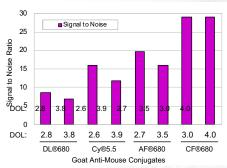


Figure 2. Intracellular staining of Jurkat cells was performed using mouse anti-human CD3 antibody or isotype control followed by goat anti-mouse secondary antibody conjugates. Fluorescence was analyzed on a BD FACSCalibur flow cytometer in the FL4 channel. Bars represent the signal-tonoise ratio of CD3-positive fluorescence to isotype control.

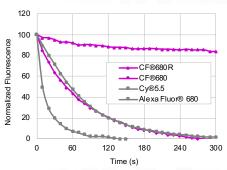


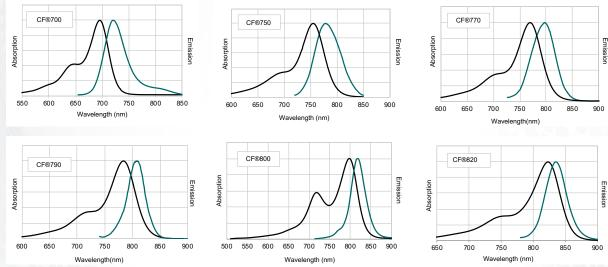
Figure 3. Relative photostability of far-red dye conjugates. Jurkat cells were stained with mouse anti-CD3 followed by the indicated goat anti-mouse IgG conjugates. Cells were continuously exposed to mercury arc lamp excitation with a Cy®5 filter set. Images were captured every 10 seconds for 5 minutes and fluorescence intensity was normalized to time 0.

CF®700 to CF®820

Unrivaled near-infrared dyes

Technical Summary

| | Dye | Ex/Em (nm) | Extinction coefficient | MW | Laser line | Spectrally similar to |
|---|--------|------------|------------------------|-------|------------|--------------------------------------|
| À | CF®700 | 695/720 | 240,000 | ~2315 | 633-685 nm | Alexa Fluor® 700, DyLight® 700 |
| | CF®750 | 755/777 | 250,000 | ~3009 | 633-685 nm | Alexa Fluor® 750, Cy®7, DyLight® 750 |
| | CF®770 | 770/797 | 220,000 | ~3138 | 785 nm | DyLight® 800, IRDye® 800CW |
| | CF®790 | 784/806 | 210,000 | ~3267 | 785 nm | Alexa Fluor® 790 |
| | CF®800 | 797/816 | 210,000 | ~3334 | 785 nm | Indocyanine Green |
| | CF®820 | 822/835 | 253,000 | ~2553 | 785 nm | Unique near-IR dye |
| | | | | | | |





Features

- Exceptionally bright and stable
- Patented PEGDye[™] technology for superior performance (see p. 3)
- Ideal for in vivo imaging
- Compatible with LI-COR® Odyssey®
- Superior signal-to-noise for conjugates
- CF®750 validated in STORM and photoacoustic imaging (see p. 21)

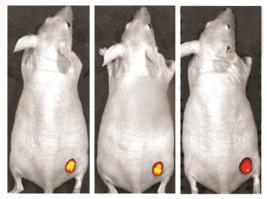


Figure 2. Tumors in mice were imaged using an IVIS® imaging system (Caliper Life Sciences) 24 hours, 48 hours, and 96 hours after IV injection of Avastin conjugated to CF®750. Images courtesy of Caliper Life Sciences.

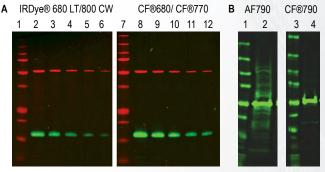


Figure 3. Near-IR western blotting with CF® Dyes compared to spectrally similar dyes. A. Two-fold dilutions of HeLa cell lysate containing from 2 ug to 0.125 ug total protein per lane were separated by SDS-PAGE, transferred to a PVDF membrane, and probed with mouse anti-tubulin and rabbit anti-COXIV antibodies. Secondary detection was performed with either IRDye® 680LT goat anti-mouse (red) and IRDye® 800CW goat anti-rabbit (green) (LI-COR®; lanes 1-6) or CF®680 goat anti-mouse (red) and CF®770 goat anti-rabbit (green) (lanes 7-12) at the same final concentrations. Membranes were scanned using an Odyssey® infrared imaging system. Quantitation showed approximately 1.5- to 2-fold higher fluorescence intensity of CF® Dye secondary antibodies compared to IRDye® secondary antibodies. B. Western blots of HeLa cell lysate (lanes 2 and 4) were probed with mouse anti-tubulin antibody followed by goat anti-mouse conjugated to Alexa Fluor® 790 (AF790, left) or CF®790 (right). CF®790 does not introduce excessive negative charge to antibody conjugates, which can increase non-specific binding.

CF® Dyes for Super-Resolution Imaging

Recent publications comparing synthetic dyes for super-resolution imaging have shown CF® Dyes give the best performance for multiple methods. The superior brightness, photostability, and photochemical switching properties of certain CF® Dyes are ideal for 3-D SIM, 3-D STORM, and other super-resolution and single molecule imaging techniques. Biotium's CF®405M has been found to be the brightest and most photostable short wavelength fluorescent dye for SIM. Six CF® Dyes spanning the visible red, far-red, and near-infrared spectra have been validated for STORM; including three color imaging with CF®568, CF®647, and CF®680.¹ See page 21 for a list of CF® Dyes validated in super-resolution and 2-photon imaging and other super-resolution techniques.

 Wide-field microscopy
 STORM

 Image: Wide-field microscopy
 CFº647

 Image: Wide-field microscopy
 CFº660C

 Image: Wide-field microscopy
 CFº680

Figure 1. Comparison of conventional wide-field microscopy (left) with STORM (right) using CF® Dye conjugates. Fixed cells were stained with mouse anti-tubulin antibody followed by CF® Dye conjugated anti-mouse secondary antibody (top row: CF®647, middle row: CF®660C, bottom row: CF®680). For STORM, samples were sealed in buffer that contained 5% (w/v) glucose, 100 mM cysteamine, 0.8 mg/mL glucose oxidase, and 40 µg/mL catalase, in Tris-HCI (pH 7.5). Samples were imaged on a Nikon Ti-Eclipse w/ PFS microscope with a CFI Plan Apo Lambda 100x oil objective. Dye molecules were photoswitched and imaged using a 647 nm laser; a 405 nm laser was used to assist dye reactivation to the emitting state. Emission was collected with an Andor iXon Ultra 897 EMCCD camera for a total of 100,000 frames per image at a frame rate of 110 Hz. Dr. Sam Kenny and Professor Ke Xu, College of Chemistry, University of California, Berkeley.

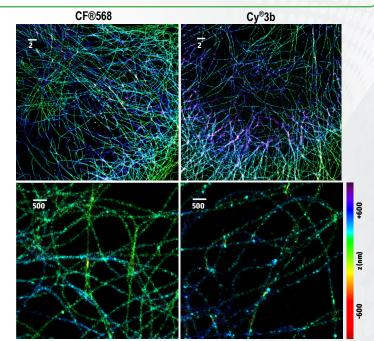


Figure 2. CF®568 (left) produces better images than Cy®3b (right) in 3-D STORM microscopy. Fixed cells were stained with mouse anti-tubulin antibody followed by dyeconjugated anti-mouse secondary antibodies. See Figure 1 for imaging conditions. Dye molecules were photoswitched and imaged using a 560 nm laser; a 405 nm laser was used to assist dye reactivation to the emitting state. Dr. Sam Kenny and Professor Ke Xu, College of Chemistry, University of California, Berkeley.

See page 4 for dye applications at a glance.

Secondary Antibodies, Single Label for STORM

1 mg/mL in PBS. 50% glycerol, 2 mg/ml BSA, 0.05% sodium azide; unit sizes: 0.5 mL, 50 uL

| • | ••• | • | | | | |
|-------------|---------------------------------------|---|---|---|-----------------------|----------------------|
| Conjugate | Donkey anti- goat | Donkey anti- guinea pig | Donkey anti- mouse (min x rat) | Donkey anti- rabbit | Goat anti- mouse | Goat anti- rabbit |
| Min x react | Ch, GP, Hs, Hu, Ms, Rb, Rt, SHm | Bv, Ch, Gt, Hs, Hu, Ms, Rb, Sh, SHm | Bv, Ch, Gt, GP, Hs, Hu, Rb, Rt, Sh, SHm | Bv, Ch, Gt, GP, Hs, Hu, Ms, Sh, SHm | Bv, Hs, Hu, Rb, Sw | Hu, Ms, Rt |
| CF®535ST | | | 20823 | 20824 | 20821 | 20822 |
| CF®568 | 20836 | 20838 | 20802 | 20803 | 20800 | 20801 |
| CF®594ST | | | 20806 | 20807 | 20804 | 20805 |
| CF®647 | 20829 | 20837 | 20810 | 20811 | 20808 | 20809 |
| CF®660C | | | 20815 | 20816 | 20812 | 20813 |
| CF®680 | | | 20819 | 20820 | 20817 | 20818 |
| CF®750 | | | 20827 | 20828 | 20825 | 20826 |
| | | | | | | |

Bv: bovine; Ch: chicken; Gt: goat; GP: guinea pig; Hs: horse; Hu: human; Ms: mouse; Rb: rabbit; Sh: sheep; SHm: Syrian hamster; Sw: swine; Rt: rat

Single-Label Secondary Antibody Conjugates for STORM

Secondary antibodies with a low degree of labeling (DOL, or number of dye molecules per antibody) have been reported to be optimal for STORM.² We offer single-label secondary antibody conjugates of our STORM-compatible dyes with an average DOL of one dye per antibody.

See page 26-29 for our full selection of antibody conjugates.

CF® Dyes for Super-Resolution Imaging

& Other specialized applications

| | Abs/Em | Extinction | | |
|----------|------------|-------------|--|--|
| CF® Dye | maxima | coefficient | Application | References |
| CF®405S | 404/431 nm | 33,000 | SIM | Demmerle, J. et al. (2017). <u>Nature Protocols 12, 988–1010.</u> Essig, K. et al. (2017). <u>Immunity https://doi.org/10.1016/j.immuni.2017.11.008</u> |
| CF®405M | 408/452 nm | 41,000 | SIM, STED | Kraus, F. et al. (2017). <u>Nat Protoc 12, 1011-1028. doi:nprot.2017.020 (SIM)</u> Markaki, Y. et al. (2013). <u>Methods Mol Biol 950, 43-64</u> . (SIM) Miron, E. et. al. (2016). In: Mark C. Leake (ed.), <u>Methods in Molecular Biology. vol. 1431, 127-140</u> . (SIM) Ohgomori, T. et al. (2017). <u>Eur J Neurosci 46, 2001-2014. doi:10.1111/ejn.13650</u> (SIM) Zhang, R. et al. (2019). bioRxiv <u>doi: https://doi.org/10.1101/586461</u> (STED) |
| CF®488A | 490/515 nm | 70,000 | STED, STORM, TIRF, 2-Photon | Angelov, B. & Angelova, A. (2017). <u>Nanoscale 9, 9797-9804. doi:10.1039/c7nr03454g</u> (STED) Mercier, L. et al. (2016). <u>Intravital 5, e1168553</u> (2-Photon) Zanetti-Domingues, L.C. et al. (2013). <u>PLoS ONE 8(9): e74200.</u> (TIRF) Collaborator communication (STORM); contact tech support through our website for more information. |
| CF®535ST | 535/568 nm | 95,000 | STORM | Collaborator communication; contact tech support through our website for more information. |
| CF®555 | 555/565 nm | 150,000 | Multicolor STORM | Lehmann, M. et al. (2015). J Biophotonics DOI 10.1002/jbio.201500119 |
| CF®568 | 562/583 nm | 100,000 | Multicolor STORM, SIM, TIRF | Gong, YN. et al. (2017). <u>Cell Cycle, 1-13. doi:10.1080/15384101.2017.1371889</u> (STORM) Gorur, A. et al. (2017). <u>J Cell Biol 216. 1745-1759. doi:10.1083/icb.201702135</u> (STORM) Heller, J. (2017). <u>OM&P 3, 48-58. doi:doi:10.20388/omp2017.002.0045</u> (STORM) Jorgans, D.M. et al. (2017). <u>J Cell Sci 2017 130: 177-189. doi: 10.1242/jcs.190967</u> (STORM) Karanasios, E. et al. (2016). <u>Nat Commun 7: 12420. DOI: 10.1038/ncomms12420</u> (STORM) Kraus, F. et al. (2017). <u>Nat Protoc 12, 1011-1028</u> doi:nprot.2017.020 (SIM) Lehmann, M. et al. (2015). <u>J Biophotonics DOI 10.1002/jbio.201500119</u> (STORM) Lim, A. et al. (2017). <u>Mol Biol Cell doi:mbc.E16-12-0820</u> (SIM) Turkowyd, B. et al. (2016). <u>Anal Bioanal Chem DOI 10.1007/s00216-016-9781-8</u> (STORM) Zanetti-Domingues, L.C. et al. (2013). <u>PLoS ONE 8(9): e74200.</u> (TIRF) Zhang, M. et al. (2015). <u>eLife 2015;10.7554/eLife.11205</u> (STORM) |
| CF®594 | 593/614 nm | 115,000 | 2-Photon | Wagner, M.C. (2016). Am J Physiol Renal Physiol310: F1089–F1102. (2-Photon) |
| CF®594ST | 593/614 nm | 115,000 | STORM | Collaborator communication; contact tech support through our website for more information. |
| CF®633 | 630/650 nm | 100,000 | FIONA, gSHRImP, Single molecule tracking, TIRF | Bosch, P. J. et al. (2014). <u>Biophys J 107, 803-814.</u> (TIRF) Huang, T. et al. (2018). <u>Biophysical Journal 114, 301–310.</u> (Single Molecule Tracking) Kim, H. J., and Selvin, P. R. (2013). <u>SpringerReference Encyclopedia of Biophysics.</u> (FIONA) Simonson, P. D. et al. (2011). <u>Nano Lett 11, 5090-5096. DOI:10.1021/nl203560r</u> (gSHRImP) Zanetti-Domingues, L.C. et al. (2013). <u>PLoS ONE 8(9): e74200.</u> (TIRF) Zhang, R. et al. (2017). <u>eLife 2017:6:e30959.</u> (TIRF) |
| CF®640R | 642/662 nm | 105,000 | FLIMP, SIM, TIRF | Bosch, P. J. et al. (2014). <u>Biophys J 107, 803-814.</u> (TIRF) Loh, L. N. (2017). MBio 8, <u>doi:mBio.02030-16</u> (SIM) Martin-Fernandez, M. L. et al. (2013). <u>J Microsc 252, 16-22.</u> (TIRF) Needham, S.R. et al. (2015). <u>Biochem Soc Trans 43, 309–314.</u> (FLImP) Needham, S.R. et al. (2016). <u>Nat Commun 7, 13307. doi:ncomms13307</u> (FLImP) Zanetti-Domingues, L.C. et al. (2013). <u>PLoS ONE 8(9): e74200.</u> (TIRF) Zanetti-Domingues, L.C. et al. (2015). <u>Prog Biophys Mol Biol. doi:S0079-6107(15)00047-4</u> (FLImP) |
| CF®647 | 650/665 nm | 240,000 | Multicolor STORM | Gong, YN. et al. (2017). <u>Cell Cycle, 1-13. doi:10.1080/15384101.2017.1371889</u> Lehmann, M. et al. (2015). <u>J Biophotonics DOI 10.1002/jbio.201500119</u> Olivier, N. et al. (2013). <u>Biomed Opt Express 4, 885-899.</u> Turkowyd, B. et al. (2016). <u>Anal Bioanal Chem DOI 10.1007/s00216-016-9781-8</u> |
| CF®660C | 667/685 nm | 200,000 | Multicolor STORM | Turkowyd, B. et al. (2016). <u>Anal Bioanal Chem DOI 10.1007/s00216-016-9781-8</u> Zhang, Z., et al. (2015). <u>Nature Methods doi:10.1038/nmeth.3528.</u> |
| CF®680 | 681/698 nm | 210,000 | Dual-Color 3D SMLM, Multicolor STORM | Früh, S.M. et al. (2015). <u>Nature Communications 6, 7275.</u> (STORM) Glebov, O.O. et al. (2017). <u>Cell Rep 18, 2715-2728. doi:S2211-1247(17)30279-6 (STORM)</u> Gorur, A. et al. (2017). <u>J Cell Biol 216, 1745-1759. doi:10.1083/jcb.201702135</u> (STORM) Lehmann, M. et al. (2015). J <u>Biophotonics DOI 10.1002/jbio.201500119</u> (STORM) Platonova, E. et al. (2015). <u>ACS Chem. Biol.10(6),1411–1416.</u> (STORM) Platonova, E. et al. (2015). <u>Methods doi: http://dx.doi.org/10.1016/j.ymeth.2015.06.018.</u> (STORM) Salvador-Gallego, R. et al. (2016). <u>EMBO J. DOI 10.15252/embj.201593384.</u> (STORM) Shrestha, R. L. et al. (2017). <u>Nat Commun 8, 150. doi:10.1038/s41467-017-00209-z (STORM)</u> Turkowyd, B. et al. (2016). <u>Anal Bioanal Chem DOI 10.1007/s00216-016-9781-8</u> (SMLM) Winterflood, C.M. et al. (2015). <u>Biophys J. 109, 3–6.</u> (SMLM) Zhang, Z., et al. (2015). <u>Nature Methods doi:10.1038/nmeth.3528.</u> (STORM) |

FIONA: Fluorescence Imaging with One Nanometer Accuracy; FLImP: Fluorophore localization imaging with photobleaching; SHRImP: Single-molecule High-Resolution Imaging with Photobleaching; SIM: Structured Illumination Microscopy; STED: Stimulated Emission Depletion; STORM: Stochastical Optical Reconstruction Microscopy; TIRF: Total Internal Reflection Fluorescence

Mix-n-Stain™ Antibody Labeling Kits

Mix-n-Stain™ CF® Dye & Hapten Antibody Labeling Kits

- Labeling in 15 minutes with minimal hands-on time & no purification
- · Covalent conjugation, suitable for multiplex staining
- Choice of small-scale labeling sizes to conserve precious antibodies
- Reaction tolerates common antibody buffers & stabilizers
- Modified protocol for antibodies with excess BSA/gelatin or ascites

Mix-n-Stain™ Labeled Antibodies Perform Better Than Lightning-Link® Labeled Antibodies

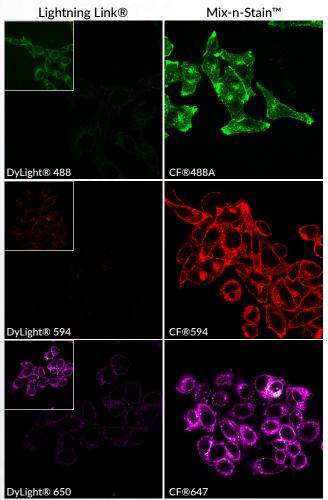


Figure 2. Mouse anti-transferrin receptor antibody from BD Biosciences (endosome and plasma membrane marker) was labeled using Lightning-Link® Rapid Conjugation Kits with the indicated DyLight® dyes (left) or Mix-n-Stain™ CF® Dye Antibody Labeling Kits (right). CF® Dye conjugates shows brighter signal and more specific staining compared to the spectrally similar DyLight® conjugates using the same laser and gain settings. The insets show the DyLight® conjugates imaged with higher gain settings to demonstrate the presence of cells in the field of view.



Figure 1. Mix-n-Stain[™] CF® Dye labeling protocol. Simply mix your antibody with the reaction buffer and pre-measured dye, and incubate 15 minutes for a ready-to-use conjugate covalently labeled with one of our bright & photostable CF® Dyes, biotin, or other label.

Mix-n-Stain™ CF® Dye or Hapten Antibody Labeling Kits

| | in a blain ci bycor napten intibouy Labernig (115 | | | | | |
|--------------|---|-------------------------|--------------------------|---------------------------|---------------------------------------|--|
| Dye or Label | Ex/Em (nm) | 1 x 5-20 ug labeling | 1 x 20-50 ug labeling | 1 x 50-100 ug labeling | Mix-n-Stain™ Maxi 1 mg labeling | |
| CF®350 | 347/448 | 92270 | 92250 | 92230 | 92420 | |
| CF®405S | 404/431 | 92271 | 92251 | 92231 | 92421 | |
| CF®405M | 408/452 | 92272 | 92252 | 92232 | 92404 | |
| CF®405L | 395/495 | 92303 | 92304 | 92305 | | |
| CF®430 | 426/498 | 92316 | 92317 | 92318 | | |
| CF®440 | 440/515 | 92319 | 92320 | 92321 | | |
| CF®450 | 405/460 | 92322 | 92323 | 92324 | | |
| CF®488A | 490/515 | 92273 | 92253 | 92233 | 92405 | |
| CF®514 | 516/548 | 92331 | 92332 | 92333 | | |
| CF®532 | 527/558 | 92289 | 92290 | 92291 | | |
| CF®543 | 541/560 | 92287 | 92267 | 92247 | | |
| CF®555 | 555/565 | 92274 | 92254 | 92234 | 92406 | |
| CF®568 | 562/583 | 92275 | 92255 | 92235 | 92407 | |
| CF®570 | 568/591 | 92334 | 92335 | 92336 | | |
| CF®583 | 586/609 | 92337 | 92338 | 92339 | | |
| CF®594 | 593/614 | 92276 | 92256 | 92236 | 92408 | |
| CF®633 | 630/650 | 92277 | 92257 | 92237 | 92409 | |
| CF®640R | 642/662 | 92278 | 92258 | 92245 | | |
| CF®647 | 650/665 | 92279 | 92259 | 92238 | 92410 | |
| CF®660C | 667/685 | 92280 | 92260 | 92239 | | |
| CF®660R | 663/682 | 92281 | 92261 | 92243 | | |
| CF®680 | 681/698 | 92282 | 92262 | 92240 | 92422 | |
| CF®680R | 680/701 | 92283 | 92263 | 92246 | | |
| CF®700 | 695/720 | 92425 | 92426 | 92427 | | |
| CF®750 | 755/777 | 92284 | 92264 | 92241 | 92423 | |
| CF®770 | 770/797 | 92285 | 92265 | 92242 | 92424 | |
| CF®790 | 784/806 | 92288 | 92268 | 92248 | | |
| CF®800 | 797/816 | 92428 | 92429 | 92430 | | |
| CF®820 | 822/835 | 92431 | 92432 | 92433 | | |
| FITC | 494/519 | 92294 | 92295 | 92296 | 92411 | |
| Cyanine 555 | 555/565 | 92412 | 92413 | 92414 | 92415 | |
| Cyanine 647 | 650/665 | 92416 | 92417 | 92418 | 92419 | |
| Biotin | N/A | 92286 | 92266 | 92244 | | |
| Digoxigenin | N/A | 92328 | 92329 | 92330 | | |
| DNP | N/A | 92325 | 92326 | 92327 | | |

Antibody, Protein, and Ligand Labeling Kits

Mix-n-Stain™ Enzyme or Fluorescent Protein Antibody Labeling Kits

- · Easy conjugation in just a few hours with no special equipment required
- Choose AP, HRP, or GOx enzyme conjugation
- Easy labeling with R-PE, APC, PerCP, or tandem dyes for flow cytometry

Mix-n-Stain™ Enzyme Antibody Labeling Kits

| Conjugation | 1 x 10-20 ug labeling | 1 x 25-50 ug labeling | 1 x 50-100 ug labeling | 1 x 1 mg labeling |
|------------------------------|--------------------------|--------------------------|---------------------------|----------------------|
| Horseradish peroxidase (HRP) | 92300 | 92301 | 92302 | 92437 |
| Alkaline phosphatase (AP) | | 92314 | 92315 | |
| Glucose oxidase (GOx) | | 92312 | 92313 | |

Mix-n-Stain™ Fluorescent Protein Antibody Labeling Kits

| Conjugation | Ex/Em (nm) | 1 x 25-50 ug labeling | 1 x 50-100 ug labeling | 1 x 1 mg labeling |
|--------------|-------------|--------------------------|---------------------------|----------------------|
| R-PE | 496,564/578 | 92298 | 92299 | |
| R-PE-CF®647T | 496/665 | 92340 | 92341 | 92346 |
| R-PE-CF®583R | 496/609 | 92442 | 92443 | |
| APC | 650/660 | 92306 | 92307 | |
| Per-CP | 482/678 | 92308 | 92309 | |
| APC-CF®750T | 650/780 | 92310 | 92311 | |

CF® Dye SE and VivoBrite™ Protein Labeling Kits

- Everything you need to label and purify 3 x 1 mg antibody
- VivoBrite™ kits feature superior near-IR CF® Dyes for *in vivo* imaging, and 0.2 um sterile mini-syringe filters

CF® Dye or Biotin SE Protein Labeling Kits

| | | - | |
|--------------|------------|----------|-------|
| Dye or Label | Ex/Em (nm) | Cat. No. | NIR S |
| CF®350 | 347/448 | 92210 | |
| CF®405S | 404/431 | 92211 | CF |
| CF®405M | 408/452 | 92212 | CF |
| CF®405L | 395/495 | 92228 | CF |
| CF®488A | 490/515 | 92213 | CF |
| CF®532 | 527/558 | 92208 | |
| CF®543 | 541/560 | 92209 | |
| CF®555 | 555/565 | 92214 | |
| CF®568 | 562/583 | 92215 | |
| CF®594 | 593/614 | 92216 | |
| CF®633 | 630/650 | 92217 | |
| CF®640R | 642/662 | 92225 | |
| CF®647 | 650/665 | 92218 | |
| CF®660C | 667/685 | 92219 | |
| CF®660R | 663/682 | 92223 | |
| CF®680 | 681/698 | 92220 | |
| CF®680R | 680/701 | 92226 | |
| CF®750 | 755/777 | 92221 | |
| CF®770 | 770/797 | 92222 | |
| Biotin | N/A | 92224 | |

VivoBrite[™] Antibody Labeling Kits for NIR Small Animal *In Vivo* Imaging

| Dye | Ex/Em (nm) | Cat. No. |
|--------|------------|----------|
| CF®680 | 681/698 | 92160 |
| CF®750 | 755/777 | 92161 |
| CF®770 | 770/797 | 92162 |
| CF®790 | 784/806 | 92163 |

Mix-n-Stain™ Small Ligand Labeling Kits

- For labeling small molecules with primary amines
- Label SNAP-Tag®, CLIP-Tag™, HALO-Tag®, or TMP-tag ligands
- 30 minute labeling with minimal hands-on time and no purification
- Choose from 10 CF® Dye colors for surface targets, or 5 CF® Dye colors for intracellular targets

Mix-n-Stain™ Small Ligand Labeling Kits

| | | - | |
|---------|------------|----------------------|-----------------------|
| Dye | Ex/Em (nm) | Cell surface targets | Intracellular targets |
| CF®405M | 408/452 | 92362 | |
| CF®408 | 408/450 | | 92356 |
| CF®488A | 490/515 | 92350 | |
| CF®500 | 500/510 | | 92357 |
| CF®540 | 540/570 | | 92358 |
| CF®555 | 555/565 | | 92364 |
| CF®568 | 562/583 | 92351 | |
| CF®594 | 593/614 | 92352 | |
| CF®633 | 630/650 | 92353 | |
| CF®640R | 642/662 | 92354 | |
| CF®647 | 650/665 | 92359 | |
| CF®650 | 650/670 | | 92363 |
| CF®660C | 667/685 | 92360 | |
| CF®680 | 681/698 | 92361 | |
| CF®680R | 680/701 | 92355 | |
| | | | |

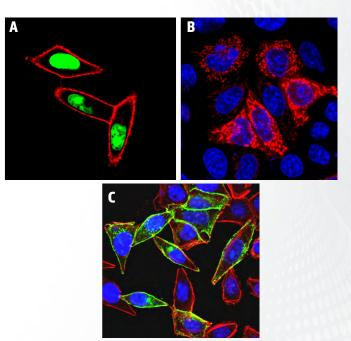


Figure 1. Versatility of Mix-n-Stain [™]-labeled ligands for multicolor live cell imaging. (A) CF®500-labeled CLIP-Tag[™] ligand was used to detect nuclear protein H2B (green), and CF®568-labeled SNAP-Tag® ligand was used to detect cell surface protein ADRβ2. (B) CF®540-labeled CLIP-Tag[™] ligand was used to detect mitochondrial protein Cox8A in living cells (red); nuclei were stained with Hoechst 33342 (blue). (C) Three color imaging in fixed cells. CF®488A-labeled CLIP-Tag[™] ligand was used to stain cell surface protein NK1R (green). Cells were then fixed and stained with CF®633 phalloidin (red) and mounted with EverBrite[™] mounting medium with DAPI (blue).

CF® Dye: Reactive Dyes

A wide selection of colors and functional groups for dye conjugation

| Reacts withAzides, picolyl azidesActivated carboxylic acidsAldehydes & ketonesAlkynes, BCN (Cu-catalyzed)Azides (Cu-free)Polar tracer1ThiolsAlkynes (low [Cu])Primary amines lysine residues | HRP substrate |
|--|---------------|
| Size 0.5 mg 1 mg 1 mg 0.5 mg 0.5 mg 1 mg 1 mg 0.5 mg 1 umol | 0.5 mg |
| CF®350 92035 92050 92151 92020 92109 | 92170 |
| CF®405S 92036 92055 92113 92183 92030 92110 | 92197 |
| CF®405M 92093 92056 92092 92114 92021 92111 | 96057 |
| CF®405L 92046 92112 | 92198 |
| CF®430 96063 92118 92117 | 96053 |
| CF®440 96070 ³ 96064 92124 92123 | |
| CF®450 96012 96011 | |
| CF®488A 92086 92037 92051 92080 92075 92152 92022 92097 92187 92120 | 92171 |
| CF®503R 96026 ³ 96079 96078 | |
| CF®514 92103 | 92199 |
| CF®532 92180 92045 92104 | 96066 |
| CF®543 92181 92044 92098 92105 | 92172 |
| CF®550R 92087 92038 92081 92153 96704 96073 | 96077 |
| CF®568 92088 92039 92057 92082 92076 92154 92024 92188 92131 | 92173 |
| CF®570 96015 96014 | |
| CF®583 96017 96016 | |
| CF®594 92089 92040 92052 92083 92077 92158 92025 92099 92189 92132 | 92174 |
| CF®620R 92033 92106 | 92194 |
| CF®633 92041 92053 92156 92026 92133 | |
| CF®640R 92091 92043 92058 92085 92078 92157 92034 92096 92190 92108 | 92175 |
| CF®647 92090 92042 92084 92136 92027 92191 92135 | 96022 |
| CF®650 96027 ³ | |
| CF®660C 92095 92094 92028 96001 92137 | |
| CF®660R 96004 96010 92059 92182 96024 92031 96002 92134 | 92195 |
| CF®680 96005 92119 92029 96003 92139 | |
| CF®680R 96006 92054 96000 92079 96025 92032 96007 92107 | 92196 |
| CF®700 96067 | |
| CF®750 92102 96062 92142 | 96052 |
| CF®770 92065 92192 92150 | |
| CF®790 921554 | |
| CF®800 92128 921274 | |
| CF®820 960684 | |

¹ For conjugation to aldehyde or ketone groups, we recommend using CF® Dye aminooxy forms.

²See page 30 for Tyramide Amplification Kits and Ready-to-Use Tyramide Amplification Buffer.

³ Membrane-permeant, compatible with intracellular copper-free reaction with azide.

⁴ Size: 0.25 umol

Don't see what you're looking for?

We regularly add new CF® Dye products to our catalog according to customer demand. Be sure to check our website for updates. If you are looking for a CF® Dye product not listed in our catalog, please contact tech support through our website. We may be able to add it as a new product, or perform a custom synthesis for you.

Visit www.biotium.com to see our full selection of reactive biotin reagents, traditional reactive dyes, cyanine dyes, and sets of size- and chargematched dyes for DIGE.

CF® Dye Bioconjugates

Bioconjugate Applications

| Conjugate | Application |
|-----------------------------|---|
| Annexin V | Phosphatidylserine probe; apoptotic cell surface marker Available in solution with azide, or lyophilized, azide-free for real-time imaging |
| a-Bungarotoxin (BTX) | Acetylcholine receptor probe; neuromuscular junction stain |
| Bovine serum albumin (BSA) | Fluid-phase endocytosis tracer; in vivo blood flow tracer |
| Cholera Toxin Subunit B | GM1 receptor probe; lipid raft, endocytic vesicle, neuronal tracing |
| Concanavalin A (Con A) | Lectin; binds α -D-mannosyl and α -D-glucosyl groups, stains yeast cell wall |
| Dextran amine, anionic | Fixable fluid-phase endocytosis tracer |
| Nucleotide conjugates | Fluorescent DNA or RNA probe synthesis; TUNEL apoptosis assay |
| Phalloidin | Filamentous actin probe |
| Peanut agglutinin (PNA) | Lectin; specific for terminal b-galactose |
| Streptavidin | Detection of biotinylated probes |
| Transferrin (human) | Recycling endosome tracer |
| Wheat germ agglutinin (WGA) | Lectin, binds N-acetyl-D-glucosamine and sialic acid; Fluorescent bacterial Gram stain, stains yeast bud scars |

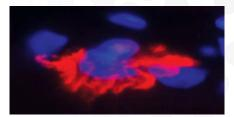


Figure 1. Frozen section of rat skeletal muscle stained with CF®633 a-bungarotoxin (magenta) to detect nicotinic acetylcholine receptors at the neuromuscular junction. Nuclei are stained with DAPI (blue).

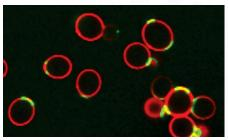


Figure 2. S. cerevisiae yeast stained with CF®488A WGA and CF®594 ConA. ConA (red) stains the cell wall, while WGA (green) preferentially stains bud scars.

CF® Dye Bioconjugates

| Dye | Annexin V | Annexin V, azide-free | a- BTX | BSA | Cholera Toxin B | Con A | Dextran 3.5K | Dextran 10K | Dextran 40K | Dextran 70K | Dextran 150K | Dextran 250K | Phalloidin | PNA | Streptavidin | Transferrin | WGA |
|---------|-----------|--------------------------|--------|-------|--------------------|-------|-----------------|----------------|----------------|----------------|-----------------|-----------------|------------|-------|--------------|-------------|-------|
| CF®350 | 29012 | 29012R-5ug | | | | 29015 | 80137 | | | | | | 00049 | | 29031 | | 29021 |
| CF®405S | | | 00002 | | | 29075 | | | | | | | | | 29032 | | 29027 |
| CF®405M | 29009 | 29009R-5ug | | | | 29074 | | | | | | | 00034 | | 29033 | | 29028 |
| CF®405L | | | | | | | | | | | | | | | 29056 | | |
| CF®430 | | | | | | | | | | | | | 00054 | | 29065 | | |
| CF®440 | | | | | | | | | | | | | 00055 | | 29066 | | |
| CF®450 | 29083 | 29083R-5ug | | | | | | | | | | | | | | | |
| CF®488A | 29005 | 29005R-5ug | 00005 | 20289 | 00070 | 29016 | | 80110 | 80126 | 80117 | 80131 | 80134 | 00042 | 29060 | 29034 | 00081 | 29022 |
| CF®514 | | | | | | | | | | | | | | | 29081 | | |
| CF®532 | | | | | 00074 | | | | | | | | 00051 | | 29030 | | 29064 |
| CF®543 | | | 00026 | | 00075 | | | 80111 | | | | | 00043 | | 29043 | 00082 | |
| CF®555 | 29004 | 29004R-5ug | 00018 | | | | | 80112 | | | | | 00040 | | 29038 | | 29076 |
| CF®568 | 29010 | 29010R-5ug | 00006 | | 00071 | | | 80113 | | | | | 00044 | 29061 | 29035 | 00083 | 29077 |
| CF®583R | | 29085R-5ug | | | | | | | | | | | 00064 | | | | |
| CF®594 | 29011 | 29011R-5ug | 00007 | 20290 | 00072 | 29017 | | 80114 | | | | | 00045 | 29062 | 29036 | 00084 | 29023 |
| CF®620R | | | | | 00076 | | | | | | | | | | | | |
| CF®633 | 29008 | 29008R-5ug | 00009 | | 00077 | 29018 | | | | | | | 00046 | | 29037 | | 29024 |
| CF®640R | 29014 | 29014R-5ug | 00004 | 20291 | 00073 | 29019 | | 80115 | | | | | 00050 | 29063 | 29041 | 00085 | 29026 |
| CF®647 | 29003 | 29003R-5ug | | | | | | | | | | | 00041 | | 29039 | | |
| CF®660C | | | | | | | | | | | | | 00052 | | | | |
| CF®660R | 29069 | 29069R-5ug | | | 00078 | | | | | | | | 00047 | | 29040 | | |
| CF®680 | | 29007 | | 20292 | | 29020 | | 80118 | 80127 | 80129 | 80132 | 80135 | 00053 | | | | 29029 |
| CF®680R | | 29070 | 00003 | | 00079 | | | 80116 | | | | | 00048 | | 29072 | 00086 | 29025 |
| CF®700 | | 29082 | | | | | | | | | | | | | | | |
| CF®750 | | 29006 | | | | 29080 | | 80119 | 80128 | 80130 | 80133 | 80136 | | | | 00087 | |
| CF®770 | | 29046 | | | | 29058 | | 80120 | 80122 | 80123 | 80124 | 80125 | | | | | 29059 |
| CF®790 | | 29047 | | | | | | 80121 | | | | | | | | | |
| CF®800 | | 29078 | | | | | | -1-1-1-1 | | | | line and the | | | | | |

Visit www.biotium.com to see our selection of apoptosis staining kits, bacterial Gram stain kits, and phalloidin conjugates of biotin and traditional dyes.

Nucleotide Conjugates

| Nucleotide | CF®405S | CF®405M | CF®488A | CF®532 | CF®543 | CF®555 | CF®568 | CF®594 | CF®640R | CF®647 | CF®660R | CF®680R |
|------------|---------|---------|---------|--------|--------|--------|--------|--------|---------|--------|---------|---------|
| dCTP | | | 40067 | 40057 | 40058 | 40027 | 40055 | 40056 | 40066 | 40028 | 40068 | |
| ddCTP | | | | | | 40031 | | | | | | |
| UTP | | | | | | | | | 40032 | | | |
| dUTP | 40004 | 40100 | 40008 | | 40002 | | 40005 | 40006 | 40007 | | | 40003 |

Visit www.biotium.com to see our CF® Dye TUNEL staining kits, plus a selection of nucleotide conjugates of biotin, traditional dyes, and amino-allyl nucleotides.

Primary Antibody Conjugates

Features

- More than 1000 monoclonal antibodies
- Growing selection of recombinant monoclonal mAbs & monoclonal rabbit antibodies
- Validated in IHC and other applications
- Select mAbs verified as monospecific in HuProt[™] human protein array
- Choice of 13 bright and photostable CF® Dyes
- Also available with R-PE, APC, PerCP, HRP, AP, or biotin
- Matched isotype controls for mouse and rabbit monoclonal antibodies
- Purified antibodies available BSA-free, 1 mg/mL, and ready-to-use for Mix-n-Stain[™] labeling or other conjugation

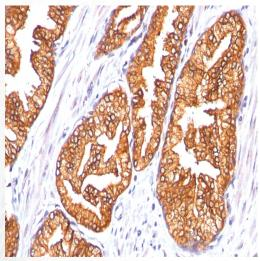
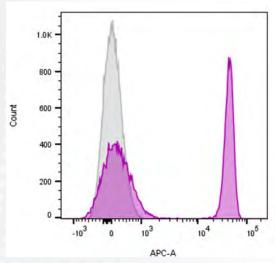
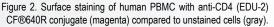


Figure 1. IHC staining of human prostate carcinoma with anti-ODC1 clone ODC1/485.





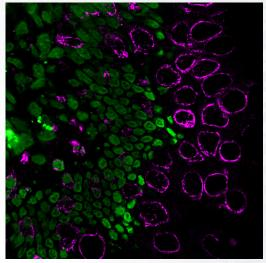


Figure 3. Immunofluorescence staining of rat jejunum with CF \otimes 488A mouse anti-Histone H1 (nuclei, green) and CF \otimes 647 mouse anti-Pan Cytokeratin (microfilaments, magenta).

Your Choice of Size and Format

| Format | Concentration | Size |
|---------------------------------|---------------|---------------|
| CF® Dye conjugates (13 colors) | 0.1 mg/mL | 100 or 500 uL |
| Biotin, HRP, or AP conjugates | 0.1 mg/mL | 100 or 500 uL |
| R-PE, APC, or Per-CP conjugates | 0.1 mg/mL | 250 uL |
| Purified, with BSA | 0.1 mg/mL | 100 or 500 uL |
| Purified, BSA-free | 1 mg/mL | 50 uL |

Your Choice of 13 Bright and Photostable CF® Dyes

| CF® Dye | Ex/Em (nm) | Features |
|---------|------------|---|
| CF®405S | 404/431 | Better fit for the 450/50 flow cytometer channel than Alexa Fluor® 405 |
| CF®405M | 408/452 | More photostable than Pacific Blue [™], with less green spill-over Compatible with super-resolution imaging by SIM |
| CF®488A | 490/515 | Less non-specific binding and spill-over than Alexa Fluor® 488 Very photostable and pH-insensitive Compatible with super-resolution imaging by TIRF |
| CF®543 | 541/560 | Brighter than Alexa Fluor® 546 |
| CF®555 | 555/565 | Brighter than Cy®3 Validated in multicolor super-resolution imaging by STORM |
| CF®568 | 562/583 | Optimized for the 568 nm line of the Ar-Kr mixed-gas Brighter and more photostable than Alexa Fluor® 568 Compatible with TIRF and multicolor STORM |
| CF®594 | 593/614 | Brighter than Texas Red® or Alexa Fluor® 594 Extremely photostable |
| CF®640R | 642/662 | Most photostable Cy®5-like dye with excellent brightness Compatible with TIRF and FLImP super-resolution techniques |
| CF®647 | 650/665 | Brighter than Cy®5 Compatible with super-resolution imaging by STORM |
| CF®660R | 663/682 | Brighter than Alexa Fluor® 660, remarkably photostable |
| CF®680 | 681/698 | Brighter than Cy®5.5, Alexa Fluor® 680, or IRDye® 680LT Validated in STORM and 3D super-resolution imaging Compatible with LI-COR® Odyssey® System |
| CF®680R | 680/701 | The most photostable 680 excitable dye Compatible with LI-COR® Odyssey® System |
| CF®770 | 770/797 | Exceptionally bright and stable Compatible with LI-COR® Odyssey® System Replacement for DyLight® 800 or IRDye® 800CW |

CF® Dye Anti-Tag and Secondary Antibody Conjugates

Anti-GFP, Anti-Hapten, and Anti-Epitope Tag Antibody Conjugates

In PBS, 50% glycerol, 2 mg/ml BSA, 0.05% sodium azide

| Conjugate | Goat anti- GST 1 mg/mL 0.1 mL | Mouse monoclonal anti-biotin 2 mg/mL 50 uL or 0.25 mL | Mouse monoclonal anti-fluorescein 2 mg/mL 50 uL or 0.25 mL | Mouse monoclonal anti-GFP 1 mg/mL 0.1 mL | Mouse monoclonal anti-6X His tag 1 mg/mL 50 uL | Rabbit anti- HA tag 1 mg/mL 50 uL | Rabbit anti- RFP 1 mg/mL 0.1 mL | Rabbit anti- V5 tag 1 mg/mL 0.1 mL |
|-----------|--|--|---|---|---|--|--|---|
| CF®405S | | 20203 | | | | | | |
| CF®405M | | | 20214 | | | | | |
| CF®488A | 20424 | 20204 | 20210 | 20215 | 20228 | 20238 | 20421 | 20440 |
| CF®543 | | | | | | | 20476 | 20441 |
| CF®568 | | | | | | | 20477 | |
| CF®588 | | | | 20480 | | | | 20441 |
| CF®594 | 20425 | 20205 | 20211 | 20216 | 20229 | 20239 | 20422 | 20442 |
| CF®633 | | 20206 | 20212 | 20217 | | | | |
| CF®640R | 20426 | 20207 | 20213 | 20218 | 20237 | 20237 | 20423 | 20443 |
| CF®647 | | | | | | 20486 | | |
| CF®660R | | | 20399 | 20481 | | | | |
| CF®680R | | | | 20482 | 20359 | | 20478 | |
| CF®750 | | 20501 | | 20220 | 20360 | | | |

Secondary Antibodies, Whole IgG (H+L), Not Cross-Adsorbed

2 mg/mL in PBS, 50% glycerol, 2 mg/ml BSA, 0.05% sodium azide, or preservative-free lyophilized form

Unit size: 0.5 mL, 50 uL, or 1 mg (lyophilized)

| Conjugate | Chicken anti-goat | Chicken anti-mouse | Chicken anti-rabbit | Goat anti- guinea pig | Goat anti- Ilama | Goat anti- mouse | Goat anti- rabbit | Goat anti- swine | Llama anti- mouse | Llama anti- rabbit | Rabbit anti- chicken | Rabbit anti- goat | Rabbit anti- guinea pig |
|-----------|----------------------|-----------------------|------------------------|--------------------------|---------------------|---------------------|----------------------|---------------------|----------------------|-----------------------|-------------------------|----------------------|----------------------------|
| CF®350 | 20364 | 20331 | 20332 | 20198 | | 20140 | 20141 | | | | | | |
| CF®405S | | | | | 20844 | 20080 | 20082 | | | | | | |
| CF®405M | | | | | | 20180 | 20181 | | | | | | |
| CF®405L | | | | | | 20408 | 20409 | | | | | | |
| CF®488A | 20225 | 20208 | 20209 | 20017 | 20845 | 20010 | 20012 | 20028 | 20454 | 20449 | 20079 | 20021 | |
| CF®514 | | | | | | 20386 | 20387 | | | | | | |
| CF®532 | | | | | | 20365 | 20366 | | | | | | |
| CF®543 | 20333 | 20334 | 20335 | 20317 | 20846 | 20306 | 20309 | 20324 | | | 20312 | 20315 | 20336 |
| CF®555 | | | | 20036 | 20847 | 20030 | 20033 | 20236 | | | | 20031 | |
| CF®568 | 20337 | 20338 | 20339 | 20108 | 20848 | 20100 | 20102 | 20091 | 20455 | 20450 | | 20107 | |
| CF®594 | 20226 | 20221 | 20223 | 20118 | 20849 | 20110 | 20112 | 20160 | 20456 | 20451 | 20164 | 20117 | |
| CF®633 | 20227 | 20222 | 20224 | 20129 | | 20120 | 20122 | 20138 | | | 20165 | 20128 | |
| CF®640R | | | | 20085 | 20850 | 20197 | 20202 | 20089 | 20457 | 20452 | | 20090 | |
| CF®647 | | | | 20041 | 20851 | 20040 | 20043 | 20286 | 20458 | 20453 | | 20049 | |
| CF®660C | | | | | 20852 | 20050 | 20053 | | | | | | |
| CF®660R | | | | | 20853 | 20054 | 20055 | | | | | | |
| CF®680 | | | | | 20855 | | | | | | | 20068 | 20243 |
| CF®750 | | | | | 20856 | 20070 | 20073 | | | | | | |
| CF®770 | | | | | | | | | | | | | 20244 |
| CF®790 | | | | | | 20378 | 20379 | | | | | | |

Don't see what you're looking for?

We regularly add new CF® Dye conjugates to our catalog according to customer demand. Be sure to check our website for updates. If you are looking for a CF® Dye product not listed in our catalog, please contact tech support through our website and let us know. We may be able to add it as a new product, or perform a custom conjugation for you.

Visit www.biotium.com to see our full selection of secondary antibodies, including conjugates of biotin, HRP, R-PE, and APC.

CF® Dye Secondary Antibody Conjugates

Highly cross-adsorbed for multiple labeling

Drop-n-Stain™ Secondary Antibodies, Whole IgG (H+L), Highly Cross-Adsorbed

5 mL solution in convenient dropper bottle format for quick and easy immunofluorescence staining

| Conjugate | Donkey anti-mouse (min x rat) | Donkey anti-rabbit | Goat anti-mouse | Goat anti-rabbit |
|-------------|--|--|-------------------|------------------|
| Min x react | Bv, Ch, Gt, GP, Hs, Hu, Rt, Rb, Sh, SHm | Bv, Ch, Gt, GP, Hs, Hu, Ms, Rt, Sh, SHm | Bv, Hs, Hu, Rb,Sw | Hu, Ms, Rt |
| CF®488A | 20952 | 20950 | 20956 | 20954 |
| CF®543 | 20967 | 20966 | 20969 | 20968 |
| CF®594 | 20953 | 20951 | 20957 | 20955 |
| CF®640R | 20963 | 20962 | 20965 | 20964 |

Bv: bovine; Ch: chicken; Gt: goat; GP: guinea pig; Hs: horse; Hu: human; Ms: mouse; Rb: rabbit; Sh: sheep; SHm: Syrian hamster; Sw: swine; Rt: rat

Secondary Antibodies, Whole IgG (H+L), Highly Cross-Adsorbed

2 mg/mL in PBS. 50% glycerol, 2 mg/ml BSA, 0.05% sodium azide, or preservative-free lyophilized form

CF®350 through CF®660R unit sizes: 0.5 mL, 50 uL, or 1 mg (lyophilized); near-IR conjugates available in 0.25 mL or 50 uL sizes

| Conjugate | Bovine anti-goat | Donkey anti- chicken | Donkey anti- goat | Donkey anti- guinea pig | Donkey anti- human | Donkey anti- mouse (min x rat) | Donkey anti- rabbit | Donkey anti- rat | Donkey anti- sheep |
|-------------|---|---|------------------------------------|---|---|---|---|--|------------------------------------|
| Min x react | Bv, Ch, GP, Hs, Hu, Ms, Rb, Rt, SHm | Bv, Gt, GP, Hs, Hu, Ms, Rb, Rt, Sh, SHm | Ch, GP, Hs, Hu, Ms, Rb, Rt, SHm | Bv, Ch, Gt, Hs, Hu, Ms, Rb, Sh, SHm | Bv, Ch, GP, Gt, Hs, Ms, Rb, Rt, Sh, SHm | Bv, Ch, Gt, GP, Hs, Hu, Rb, Rt, Sh, SHm | Bv, Ch, Gt, GP, Hs, Hu, Ms, Sh, SHm | Bv, Ch, GP, Gt, Hs, Hu, Ms, Rb, Sh, SHm | Ch, GP, Hs, Hu, Ms, Rb, Rt, SHm |
| CF®350 | | 20275 | 20142 | | | 20350 | 20351 | 20361 | 20148 |
| CF®405S | | | 20416 | 20356 | | | 20420 | 20419 | |
| CF®405M | | | 20398 | 20376 | | | | | |
| CF®430 | | | | | | 20461 | 20462 | | |
| CF®488A | 20293 | 20166 | 20016 | 20169 | 20074 | 20014 | 20015 | 20027 | 20024 |
| CF@514 | | | | | | 20483 | | | |
| CF®543 | 20313 | 20310 | 20314 | 20316 | 20318 | 20305 | 20308 | 20320 | 20322 |
| CF®555 | | | 20039 | 20276 | | 20037 | 20038 | | 20234 |
| CF®568 | 20294 | | 20106 | 20377 | | 20105 | 20098 | 20092 | 20095 |
| CF®594 | 20295 | 20167 | 20116 | 20170 | 20075 | 20115 | 20152 | 20159 | 20156 |
| CF®633 | 20296 | 20168 | 20127 | 20171 | 20076 | 20124 | 20125 | 20137 | 20134 |
| CF®640R | 20297 | | 20179 | | | 20177 | 20178 | 20199 | 20083 |
| CF®647 | | | 20048 | | | 20046 | 20047 | 20843 | 20284 |
| CF®660C | | | 20051 | 20372 | | | | | |
| CF®660R | | | 20391 | | | 20388 | 20389 | 20390 | |
| CF®680 | | | 20060 | 20241 | 20278 | | 20418 | 20417 | 20062 |
| CF®680R | | | 20196 | | | 20194 | 20195 | | |
| CF®750 | | | 20362 | | | | 20298 | 20857 | |
| CF®770 | | | 20277 | 20242 | | | 20484 | | |
| CF®790 | | | 20345 | | 20279 | 20363 | 20344 | | |
| CF@800 | | | 20834 | | | 20835 | | | |

Bv: bovine; Ch: chicken; Gt: goat; GP: guinea pig; Hs: horse; Hu: human; Ms: mouse; Rb: rabbit; Sh: sheep; SHm: Syrian hamster; Sw: swine; Rt: rat

Don't see what you're looking for?

We regularly add new CF® Dye conjugates to our catalog according to customer demand. Be sure to check our website for updates. If you are looking for a CF® Dye product not listed in our catalog, please contact tech support through our website and let us know. We may be able to add it as a new product, or perform a custom conjugation for you.

Visit www.biotium.com to see our full selection of secondary antibodies, including conjugates of biotin, HRP, R-PE, and APC.

CF® Dye Secondary Antibody Conjugates Highly cross-adsorbed, F(ab'), fragments, and isotype-specific secondary antibodies

Secondary Antibodies, Whole IgG (H+L) , Highly Cross-Adsorbed (continued from p. 28)

2 mg/mL in PBS. 50% glycerol, 2 mg/ml BSA, 0.05% sodium azide, or preservative-free lyophilized form CF®350 through CF®660R unit sizes: 0.5 mL, 50 uL, or 1 mg (lyophilized); CF®680 through CF®790 available in 0.25 mL or 50 uL sizes

| Conjugate | Goat anti- chicken | Goat anti- guinea pig | Goat anti- human | Goat anti- mouse | Goat anti-mouse (min x rat) | Goat anti- rabbit | Goat anti- rat | Rabbit anti-human | Rabbit anti- mouse | Rabbit anti- rat | Rabbit anti-sheep |
|-------------|--|---|---------------------|-----------------------|---|----------------------|-------------------|----------------------|-----------------------|---------------------|----------------------|
| Min x react | Bv, Gt, GP, Hs, Hu, Ms, Rb, Rt, Sh, SHm | Bv, Ck, Gt, Hs, Hu, Ms, Rb, Rt, SHm, Shp | Bv, Hs, Ms | Bv, Hs, Hu, Rb, Sw | Bv, Ch, Gt, GP Hs Hu Rb Rt, Sh, SHm | Hu, Ms, Rt | Bv, Hs, Hu, Rb | Ms | Hu | Hu | Hu |
| CF®350 | | | | 20143 | | 20144 | 20147 | | 20149 | | |
| CF®405S | | 20488 | | | 20830 | | | | | | |
| CF®405M | 20375 | 20487 | | 20182 | 20340 | 20373 | 20374 | | | | |
| CF®430 | | | | 20459 | | 20460 | | | | | |
| CF®488A | 20020 | 20489 | 20022 | 20018 | 20302 | 20019 | 20023 | 20071 | 20026 | 20025 | 20172 |
| CF®532 | | | | 20468 | | 20469 | | | | | |
| CF®543 | 20311 | 20492 | 20319 | 20299 | 20328 | 20300 | 20321 | | 20307 | | 20323 |
| CF®555 | 20034 | 20491 | 20320 | 20231 | | 20232 | 20233 | | 20235 | | |
| CF®568 | 20104 | 20492 | 20097 | 20101 | 20301 | 20103 | 20096 | | 20093 | 20094 | |
| CF®594 | 20114 | 20493 | 20154 | 20111 | 20303 | 20113 | 20155 | 20072 | 20158 | 20157 | 20173 |
| CF®633 | 20126 | | 20132 | 20121 | 20341 | 20123 | 20133 | 20066 | 20136 | 20135 | 20174 |
| CF®640R | 20084 | 20494 | 20081 | 20175 | 20304 | 20176 | 20088 | | 20200 | 20201 | |
| CF®647 | 20044 | 20495 | 20280 | 20281 | | 20282 | 20283 | | 20285 | | |
| CF®660C | 20371 | 20497 | | 20052 | 20368 | 20369 | 20370 | | | | |
| CF®660R | | 20496 | | | | | | | | | |
| CF®680 | | 20499 | 20287 | 20065 | | 20067 | 20069 | | 20061 | | |
| CF®680R | | 20498 | | 20192 | | 20193 | | | | | |
| CF®750 | | | | 20463 | | | | | | | |
| CF®770 | | 20500 | 20288 | 20077 | | 20078 | 20383 | | | | |
| CF®790 | | | | 20342 | | 20343 | | | | | |

Bv: bovine; Ch: chicken; Gt: goat; GP: guinea pig; Hs: horse; Hu: human; Ms: mouse; Rb: rabbit; Sh: sheep; SHm: Syrian hamster; Sw: swine; Rt: rat

Secondary Antibodies, F(ab'), Fragments

Conjugate

CF®350

CF®488A

CF®543

CF®555

CF®568

CF®594

CF®633

CF®640R

CF®647

CF®680

2 mg/mĹ, unit size: 0.25 mL or 50 uL

Goat anti-

mouse, F(ab'),

20145

20011

20329

20032

20109

20119

20130

20086

20042

20063

Goat anti-rabbit,

F(ab')₂

20146

20013

20330

20035

20099

20153

20131

20087

20045

20064

| Goat Anti-Mouse |
|-----------------------------|
| Isotype-Specific Antibodies |

2 mg/mL, unit size: 0.25 mL or 50 uL

| Conjugate | Goat anti- mouse lgG1 | Goat anti- mouse IgG2a | Goat anti-mouse IgG2b | Goat anti- mouse IgM |
|-------------|--------------------------|---------------------------|-----------------------------|-------------------------|
| Min x react | Bv, Hu, Rb | Bv, Hu, Rb | Bv, Hu, Rb | Bv, Hu, Rb |
| CF®350 | 20245 | 20255 | 20265 | |
| CF®405S | 20380 | 20381 | 20382 | |
| CF®488A | 20246 | 20256 | 20266 | 20840 |
| CF®543 | 20325 | 20356 | 20326 | |
| CF®555 | 20247 | 20257 | 20267 | 20485 |
| CF®568 | 20248 | 20258 | 20268 | |
| CF®594 | 20249 | 20259 | 20269 | |
| CF®633 | 20250 | 20260 | 20270 | |
| CF®640R | 20251 | 20261 | 20271 | |
| CF®647 | 20252 | 20262 | 20272 | |
| CF®680 | 20253 | 20263 | 20273 | 20384 |
| CF®680R | | 20842 | | 20841 |
| CF®750 | | | 20430 | |
| CF®770 | 20254 | 20264 | 20274 | 20385 |

Goat Anti-Human Isotype-Specific Antibodies 2 mg/mL, unit size: 0.25 mL or 50 uL

| Conjugate | Goat anti- human IgA (alpha chain) | Goat anti- human IgM (mu chain) | |
|-----------|--|---------------------------------------|--|
| CF®488A | 20428 | 20347 | |
| CF®594 | 20429 | 20348 | |
| CF®640R | | 20349 | |
| CF®633 | 20427 | | |
| CF®647 | | 20346 | |
| CF®680 | | 20384 | |

See more highly cross-adsorbed secondaries on the previous page; see p. 20 for single-label antibody conjugates for STORM.

Tyramides & Signal Amplification Kits

Tyramide signal amplification (TSA), sometimes called catalyzed reporter deposition (CARD), is a highly sensitive method enabling the detection of low-abundance targets immunofluorescence applications. For multiplexing, TSA not only facilitates detection of low-abundance targets, but also simplifies antibody panel design since primary antibodies of choice may be used, irrespective of host species or isotype.

In TSA, horseradish peroxidase (HRP) converts a labeled tyramide substrate into a highly reactive form that can covalently bind to tyrosine residues on proteins at or near the HRPconjugate. This generates high density tyramide labeling and is the reason for the exceptional sensitivity of this system. Because the label is covalently linked to the sample, the antibodies can be stripped off without affecting signal, allowing multiple rounds of staining for multiplex detection using antibodies from the same host species.

We offer CF® Dye and other tyramide conjugates with a wide color selection, plus easy-to-use kits and reaction buffer.

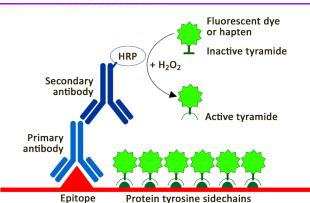


Figure 1. Illustration of tyramide signal amplification. A cell or tissue sample is labeled

with primary and secondary antibody using conventional methods. The horseradish peroxidase (HRP), conjugated to the secondary antibody, catalyzes the conversion of labeled tyramide into a reactive radical. The tyramide radical then covalently binds to nearby tyrosine residues, providing high-density labeling.

Advantages of Tyramide Signal Amplification

- Detect low-abundance targets
- ICC, IHC, and FISH-compatible
- Sensitivity up to 100-fold that of conventional methods
- · Similar workflow to conventional staining methods
- Use less antibody
- Simplify primary antibody panel design for multiplexing

Tyramide Signal Amplification Kits

Everything you need for the tyramide labeling reaction

- · Biotin tyramide or one of six CF® Dye tyramides
- HRP conjugate: goat anti-mouse, goat anti-rabbit, or streptavidin
- Tyramide Amplification Buffer Plus
- BSA (for blocking buffer preparation)

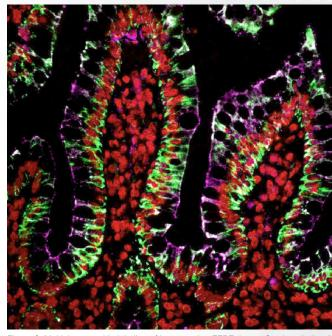


Figure 2. Multiplex tyramide labeling of human colon FFPE tissue. Cytokeratin (pan) was labeled with CF®488A tyramide (cytoskeleton, green); Histone H1 was labeled with Cyanine 555 tyramide (nuclei, red); ZO1 was labeled with CF®640R tyramide (tight junctions, magenta). All primary antibodies were from mouse; secondary antibody was HRP-conjugated goat anti-mouse. Each labeling was performed sequentially, with antibody removal by microwaving between each labeling step.

Streptavidin

HRP

33002

33005

33008

33011

33014

33017

33020

Tyramide Signal Amplification Kits

Goat anti-

rabbit HRP

33001

33004

33007

33010

33013

33016

33019

Goat anti-

mouse HRP

33000

33003

33006

33009

33012

33015

33018

Tyramide Amplification Buffer

Tyramide Amplification Buffer Plus

Tyramide

CF®488A

CF®543

CF®568

CF®594

CF®640R

CF®680R

Biotin-XX

22029

Cat. No. Product

Tyramides

| iyrannues | | |
|-------------|----------|--|
| Dye/Label | Cat. No. | |
| CF®350 | 92170 | |
| CF®405S | 92197 | |
| CF®405M | 96057 | |
| CF®405L | 92198 | |
| CF®430 | 96053 | |
| CF®488A | 92171 | |
| CF®514 | 92199 | |
| CF®532 | 96066 | |
| CF®543 | 92172 | |
| CF®550R | 96077 | |
| CF®555 | 96021 | |
| CF®568 | 92173 | |
| CF®583R | 96085 | |
| CF®594 | 92174 | |
| CF®620R | 92194 | |
| CF®640R | 92175 | |
| CF®647 | 96022 | |
| CF®660R | 92195 | |
| CF®680R | 92196 | |
| CF®750 | 96052 | |
| Biotin-XX | 92176 | |
| Fluorescein | 96018 | |
| DNP | 96019 | |
| Cyanine 555 | 96020 | |
| | | |

30 • www.biotium.com

Tyramides & Signal Amplification Kits

Background Suppressors and Accessory Products for IF/IHC/ICC

Lipofuscin autofluorescence in human cerebral cortex sections

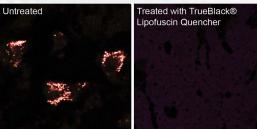


Figure 1. Left: Human brain tissue showing lipofuscin granules with bright, broad-spectrum autofluorescence that appear white in the merged image of the green, red, and far-red channels. Right: Tissue after TrueBlack® treatment, which quenches lipofuscin fluorescence.

Non-specific background from Alexa Fluor® 647 conjugate

| Non-specific background from Alexa Fluor® 047 conjugate | | |
|---|--|--|
| Gelatin blocking buffer | TrueBlack® IF Background Suppressor | |
| | | |
| S G X DRI | e e e e e e e e e e e e e e e e e e e | |
| 1843-1883 | 6 | |
| | É. | |

Figure 2. Left: Non-specific signal in HeLa cells caused by binding of negatively charged Alexa Fluor® 647 dye conjugated to secondary antibody. Right: TrueBlack® IF Background Suppressor blocks background from non-specific interactions of charged dyes with biological samples.

Our TrueBlack® line of background quenchers and blocking buffers are designed to reduce background from multiple sources, including tissue autofluorescence, non-specific antibody binding, and non-specific interactions of charged dye conjugates with cells or blotting membranes. We also offer a variety of essential accessory products for immunofluorescence staining.



Figure 3. Western detection of phospho-Erk1/2 in PDGF-stimulated NIH-3T3 cell lysate. Membranes were blocked with fish gelatin, LI-COR® Odyssey® TBS Blocking Buffer, or TrueBlack® WB Blocking Buffer. Rabbit anti-pErk1/2 and CF®680R donkey anti-rabbit antibodies were used for detection. TrueBlack® WB Blocking Buffer gave lower background fluorescence and highest specificity.

| Cat. No. | Product | Features | | |
|----------|--|--|--|--|
| 23012 | TrueBlack® IF Background Suppressor System (Permeabilizing) | Suppresses background from non-specific antibody binding and charged fluorescent dyes More efficient than Image-iT® FX, block and permeabilize in just 10 minutes Non-mammalian blocking agents for broad secondary antibody compatibility For immunofluorescence on cells or tissue sections | | |
| 23013 | TrueBlack® WB Blocking Buffer Kit | Blocks as well or better than Odyssey® Blocking Buffer, at a lower price Reduces non-specific protein bands and background from charged dyes Compatible with PVDF and nitrocellulose membranes For visible and near-IR fluorescent westerns | | |
| 23007 | TrueBlack® Lipofuscin Autofluorescence Quencher | Eliminates lipofuscin autofluorescence with less background than Sudan Black B Reduces background from other sources Can be used before or after IF staining | | |
| 23001 | EverBrite™ Mounting Medium | Excellent protection from photobleaching | | |
| 23002 | EverBrite™ Mounting Medium with DAPI | Compatible with a wide variety of dyes, including Cy®3, Cy®5, and Alexa Fluor® 647 Available in wet-set or hard-set formulations Available with or without DAPI | | |
| 23003 | EverBrite™ Hardset Mounting Medium | | | |
| 23004 | EverBrite™ Hardset Mounting Medium with DAPI | | | |
| 23008 | Drop-n-Stain EverBrite™ Mounting Medium | • EverBrite™ antifade medium in a convenient dropper bottle for pipette-free mounting | | |
| 23009 | Drop-n-Stain EverBrite™ Mounting Medium with DAPI | | | |
| 23005 | CoverGrip™ Coverslip Sealant | Replaces nail polish for coverslip sealing Won't mix with aqueous mounting media | | |
| 40061 | RedDot™2 Far Red Nuclear Counterstain | Far-red nuclear dye for the Cy®5 channel More specific than Draq7™ | | |
| 40083 | NucSpot® 470 Nuclear Stain, | Green fluorescent nuclear-specific counterstain for fixed cells or tissues | | |
| 22005 | Mini Super HT Pap Pen 2.5 mm tip, ~400 uses | Create hydrophobic barriers around tissue sections | | |
| 22006 | Super ^{HT} Pap Pen 4 mm tip, ~800 uses | Heat-stable to 120°C Insoluble in aqueous buffers, detergents, alcohol and acetone; can be removed with xylene | | |
| 22023 | Paraformaldehyde, 4% in PBS, Ready-to-Use Fixative | Ready-to-use fixation buffer Methanol-free, prepared from EM grade paraformaldehyde No glass ampoules to break, store in the original bottle | | |
| 22020 | 10X Phosphate Buffered Saline (PBS) 4L Cubitainer® | Convenient buffers and blocking agents for immunofluorescence or western | | |
| 22010 | 10X Fish Gelatin Blocking Agent | | | |
| 22014 | 30% Bovine Serum Albumin Solution | | | |
| | | | | |

Alexa Fluor and Image-iT are registered trademarks of Thermo Fisher Scientific; Cubitainer is a registered trademark of Cubitainer is a registered trademark of Hedwin Corporation; Cy dye is a registered trademarks of GE Healthcare; Draq7 is a trademark of BioStatus Ltd; Odyssey is a registered trademark of LI-COR, Inc.

About Us

At Biotium, we are dedicated to developing cutting-edge fluorescent dyes and life science assays. Innovation is at the very heart of what we do every day. Our efforts have resulted in a growing number of unique and industry-leading fluorescence-based technologies for a wide array of molecular and cellular biology applications. Our products are available in the U.S. through our website, and worldwide through our extensive network of domestic and international distributors.

CF® Dyes and Mix-n-Stain[™] antibody labeling technology are covered by granted and pending U.S. and international patents. We license our technologies to number of international biotechnology companies, and collaborate with academic laboratories to develop new tools for the ever-changing needs of the research community. We welcome inquiries about licensing the use of our dyes, technologies, or trademarks; email us at btinfo@biotium.com.

Biotium implements a Quality System, certified by QAS according to Standard QAS ISO 9001:2015.

Biotium, Inc. 46117 Landing Parkway Fremont, CA 94538 USA 800-304-5357 www.biotium.com

