CF® Dyes Quick Reference

CF® dye	λ _{Ex} (nm)	λ _{Em} (nm)	Excitation*	Replacement for	Features and applications
CF®350	347	448	UV	Alexa Fluor® 350, AMCA, DyLight® 350	Brightest blue fluorescent conjugates for 350 nm excitation Highly water-soluble and pH insensitive
CF®405S	404	431	405 nm	Alexa Fluor® 405, Cascade Blue®, DyLight® 405	Better compatibility with common instruments
CF®405M	408	452	405 nm	BD Horizon™ V450, eFluor® 450, Pacific Blue®	More photostable than Pacific Blue® dye with less green spill-over Excellent choice for super-resolution imaging by SIM
CF®405L	395	545	405 nm	Pacific Orange®	405 nm excitable orange fluorescent dye for multicolor detection
CF®430	426	498	405 nm	Pacific Green®, BD Horizon™ V500, Krome Orange™	Photostable 405 nm excitable green dye Perfect match for the CFP filter set
CF®440	440	515	405 nm	Alexa Fluor® 430	Photostable 405 nm excitable green dye
CF®450	450	538	405 nm	Unique dye	Green dye with unique spectral properties
CF®488A	490	515	488 nm	ATTO 488, Alexa Fluor® 488, Cy®2, DyLight® 488, FAM, FITC, Fluorescein	Less non-specific binding and less red spill-over than Alexa Fluor® 488 Very photostable Compatible with super-resolution imaging by TIRF
CF®514	516	548	488 nm	Alexa Fluor® 514	Green dye that can be separated from CF®488A by spectral unmixing
CF®532	527	558	532 nm	Alexa Fluor® 532, ATTO 532	Significantly brighter than Alexa Fluor® 532
CF®535ST	535	568	532 nm	Unique dye for STORM	Orange dye designed for STORM super-resolution microscopy
CF®543	541	560	532, 543, or 546 nm	Alexa Fluor® 546, Tetramethylrhodamine (TAMRA)	Brighter than Alexa Fluor® 546
CF®555	555	565	532, 543, 546,, 555, or 568 nm	Alexa Fluor® 555, ATTO 550, Cy®3, DyLight® 549, TRITC	\bullet Brighter than Cy®3 Validated in multicolor super-resolution imaging by STORM
CF®568	562	583	532, 543, 546, 555, or 568 nm	Alexa Fluor® 568, ATTO 565, Rhodamine Red	 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®570	568	591	532, 543, 546, 555, or 568 nm	Alexa Fluor® 568, ATTO 565, DY-560, Rhodamine Red	Yields brighter conjugates compared to spectrally similar dyes
CF®583	583	606	532, 543, 546, 555, or 568 nm	Cy®3.5	Yields brighter conjugates compared to spectrally similar dyes
CF®594	593	614	532, 543, 546, 555, or 568 nm	Alexa Fluor® 594, ATTO 594, DyLight® 594, Texas Red®	Yields the brightest conjugates among spectrally similar dyesExtremely photostable
CF®594ST	593	614	532, 543, 546, 555, or 568 nm	Unique dye for STORM	Specifically designed for super-resolution imaging by STORM
CF®620R	617	639	633 or 635 nm	LightCycler® Red 640	Highly fluorescent dye with unique spectral properties
CF®633	630	650	633 or 635 nm	Alexa Fluor® 633, Alexa Fluor® 647, Cy®5, DyLight® 633	Yields the brightest antibody conjugates among spectrally similar dyes Far more photostable than Alexa Fluor® 647 Compatible with super-resolution TIRF, FIONA, and gSHRImP
CF®640R	642	662	633, 635, or 640 nm	Alexa Fluor® 647, ATTO 647N, Cy®5, DyLight® 649	Has the best photostability among dyes with Cy®5-like spectra Yields highly fluorescent protein conjugates Compatible with TIRF and FLIMP super-resolution techniques
CF®647	650	665	633, 635, or 640 nm	Alexa Fluor® 647, ATTO 647N, Cy®5, DyLight® 649	Brighter than Cy®5 Compatible with multicolor super-resolution imaging by STORM
CF®660C	667	685	633, 635, or 640 nm	Alexa Fluor® 660	Much brighter and more photostable than Alexa Fluor® 660 Compatible with multicolor super-resolution imaging by STORM
CF®660R	663	682	633, 635, or 640 nm	Alexa Fluor® 660	Brighter than Alexa Fluor® 660 The most photostable 660 nm dye
CF®680	681	698	680 or 685 nm	Alexa Fluor® 680, Cy®5.5, DyLight® 680, IRDye® 680LT	The brightest among spectrally similar 680 nm dyes Validated in multicolor STORM and 2-color 3D super-resolution imaging Compatible with LI-COR® Odyssey® System
CF®680R	680	701	680 or 685 nm	Alexa Fluor® 680, Cy®5.5, DyLight® 680, IRDye® 680LT	 The most photostable 680 nm dye Suitable for labeling nucleic acids and small biomolecules Compatible with LI-COR® Odyssey® System Compatible with STED and single molecule spectroscopy
CF®750	755	777	680 or 685 nm	Alexa Fluor® 750, Cy®7, DyLight® 750, IRDye® 750	Exceptionally bright and stableHighly water soluble without bearing excessive chargeValidated in super-resolution imaging by STORM
CF®770	770	797	785 nm	DyLight® 800, IRDye® 800CW, ZW800-1	Exceptionally bright and stable Compatible with LI-COR® Odyssey® System
CF®790	784	806	785 nm	Alexa Fluor® 790	Exceptionally bright and stable Highly water soluble without bearing excessive charge

^{*}Visible and far-red dyes can be excited by a UV light source for epifluorescence microscopy.



CF® Dye Product Lines

Bioconjugates

Primary antibodies

- More than 1000 mouse & rabbit monoclonal antibodies
- Protein array-validated monospecific recombinant monoclonals
- Validated in IHC and other applications
- Choose from 13 bright and photostable CF® dyes
- Also available with R-PE, APC, PerCP, HRP, AP, or biotin
- Available purified without BSA, ready to use for Mix-n-Stain™ labeling or other conjugation
- · Offered in affordable 100 uL sizes

Anti-tag antibodies

- Antibodies against GFP, RFP, biotin, HA-tag, 6X His tag, and more
- Choose from several bright and photostable CF® dye options

Secondary antibodies

- · Wide variety of host and target species
- F(ab)'2, highly cross-adsorbed, and isotype-specific antibodies available
- · Choose from 20 bright and photostable CF® dyes
- Also available with R-PE, APC, PerCP, HRP, AP, or biotin
- Available in multiple sizes, liquid format or lyophilized

Other bioconjugates

- Streptavidin
- Phalloidin
- Annexin V
- · BSA, dextrans, transferrin, and cholera toxin
- Lectins
- Nucleotide conjugates

Reactive Dyes and Labeling Kits

Mix-n-Stain™ Antibody Labeling Kits

- · The simplest antibody labeling protocol available
- Label your antibody with your choice of more than 20 CF® dye colors, biotin, DNP, digoxygenin, or FITC in just 30 minutes, with minimal hands-on time
- No post-labeling purification required
- Labeling is covalent, suitable for multiplex staining
- Choice of small-scale labeling sizes preserves precious primary antibodies
- Kits tolerate common antibody buffer components and stabilizer proteins
- Mix-n-Stain™ kits also available for labeling antibodies with R-PE, APC, Per-CP, R-PE-CF®750T, HRP, AP, and glucose oxidase

Mix-n-Stain™ Small Ligand Labeling Kits

- · For labeling small molecules on primary amines
- Label 0.1 umol SNAP-Tag®, CLIP-Tag™, or HaloTag® ligand
- Choose from 10 CF® dye colors for surface targets, or 3 CF® dye colors for intracellular targets

CF® Dye SE and VivoBrite™ Protein Labeling Kits

- Everything you need to label and purify 3 x 1 mg antibody or other protein using standard conjugation techniques
- VivoBrite™ kits feature our superior near-IR CF® dyes for in vivo imaging

Reactive Dyes

 Our full selection of CF® dyes with a wide variety of functional groups

Related Products and Accessories

- TrueBlack® autofluorescence quencher, background suppressor, and blocking buffer for IF or western
- EverBrite™ Antifade Mounting Media
- CoverGrip™ Coverslip Sealant
- RedDot[™] and NucSpot® nuclear counterstains
- Ready-to-use, preservative-free PFA, buffers, blocking agents, Pap pens, and more

