



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

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

Label your IgG antibody with one of our superior CF® Dyes in as little as 15 minutes without a purification step. Labeling tolerates many common buffer components including BSA and ascites.



Product attributes

Product Description

Mix-n-Stain™ CF® Dye Antibody Labeling Kits allow you to label ≤ 5 ug up to 100 ug of your IgG antibody in as little as 15 minutes, with minimal hands-on time and no purification.

Note: This kit is for labeling IgG antibodies only. For labeling IgM antibodies, please see our [Mix-n-Stain™ CF® Dye IgM Antibody Labeling Kits](#).

Unrivaled Convenience

Mix-n-Stain™ labeling tolerates common antibody buffer formulations. An ultrafiltration centrifuge vial is provided to quickly remove interfering substances like glycerol, or to concentrate your antibody. With a slight modification in protocol, antibodies can be labeled in the presence of BSA, gelatin, or ascites fluid. Because CF® Dyes are highly water soluble, the presence of other proteins like BSA or gelatin in the labeling reaction has minimal effect on background fluorescence, because any labeled non-antibody proteins readily wash away during immunofluorescence staining.

Superior CF® Dyes

These kits feature a choice of 31 of Biotium's next-generation CF® Dyes, which were designed for advantages in brightness and photostability compared to other commercially available fluorescent dyes. Learn more about [CF® Dyes](#).

Choose the Right Labeling Kit for Your Antibody

Mix-n-Stain™ Antibody Labeling Kits are very simple to use (see the workflow overview above). But before you begin, you must check that your antibody meets the compatibility requirements for labeling, and choose the right labeling protocol. Download the updated and expanded [Product Protocol](#) for a labeling check list to help you select the right kit size and labeling protocol for your antibody. Also see our [Kit Compatibility and Protocol Selection Flowchart](#).

See the product table below for a quick reference of the maximum excitation and emission of each CF® Dye, or see our [CF® Dye Brochure](#) for more detailed dye information.

Other Labeling Kits

We now offer [Mix-n-Stain™ Maxi Antibody Labeling Kits](#) for labeling 1 mg of IgG antibody with CF® Dyes, as well as [Mix-n-Stain™ CF® Dye IgM Antibody Labeling Kits](#) and [Mix-n-Stain™ Nanobody Labeling Kits](#). Mix-n-Stain™ kits also are available for labeling antibodies with biotin, DNP, or DIG. See our full selection of [Mix-n-Stain™ Antibody Labeling Kits](#).

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

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF@350	347/448 nm	5-20 ug	92270	CF@350 Features
20-50 ug	92250			
50-100 ug	92230			
CF@405S	404/431 nm	5-20 ug	92271	CF@405S Features
20-50 ug	92251			
50-100 ug	92231			
CF@405M	408/452 nm	5-20 ug	92272	CF@405M Features
20-50 ug	92252			
50-100 ug	92232			
CF@405L	395/545 nm	5-20 ug	92303	CF@405L Features
20-50 ug	92304			
50-100 ug	92305			
3x(5-20 ug)	92445			
CF@430	426/498 nm	5-20 ug	92316	CF@430 Features
20-50 ug	92317			
50-100 ug	92318			
CF@440	440/515 nm	5-20 ug	92319	CF@440 Features
20-50 ug	92320			
50-100 ug	92321			
CF@450	450/538 nm	5-20 ug	92322	CF@450 Features
20-50 ug	92323			
50-100 ug	92324			
CF@488A	490/515 nm	5-20 ug	92273	CF@488A Features
20-50 ug	92253			
50-100 ug	92233			
3x(5-20 ug)	92446			
CF@514	516/548 nm	5-20 ug	92331	CF@514 Features
20-50 ug	92332			
50-100 ug	92333			
CF@532	527/558 nm	5-20 ug	92289	CF@532 Features
20-50 ug	92290			
50-100 ug	92291			
CF@543	541/560 nm	5-20 ug	92287	CF@543 Features
20-50 ug	92267			
50-100 ug	92247			
CF@555	555/565 nm	5-20 ug	92274	CF@555 Features
20-50 ug	92254			
50-100 ug	92234			
CF@568	562/583 nm	5-20 ug	92275	CF@568 Features
20-50 ug	92255			
50-100 ug	92235			
3x(5-20 ug)	92447			
CF@570	568/591 nm	5-20 ug	92334	CF@570 Features
20-50 ug	92335			
50-100 ug	92336			
CF@583	583/606 nm	5-20 ug	92337	CF@583 Features
20-50 ug	92338			
50-100 ug	92339			
CF@583R	586/609 nm	5-20 ug	92455	CF@583R Features
20-50 ug	92456			
50-100 ug	92457			
CF@594	593/614 nm	5-20 ug	92276	CF@594 Features
20-50 ug	92256			
50-100 ug	92236			
3x(5-20 ug)	92448			
CF@597R	597/619 nm	5-20 ug	92461	CF@597R Features
20-50 ug	92462			
50-100 ug	92463			
CF@633	630/650 nm	5-20 ug	92277	CF@633 Features
20-50 ug	92257			
50-100 ug	92237			
CF@640R	642/662 nm	5-20 ug	92278	CF@640R Features
20-50 ug	92258			
50-100 ug	92245			
CF@647	650/665 nm	5-20 ug	92279	CF@647 Features

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
20-50 ug	92259			
50-100 ug	92238			
3x(5-20 ug)	92449			
CF@660R	663/682 nm	5-20 ug	92281	CF@660R Features
20-50 ug	92261			
50-100 ug	92243			
CF@660C	667/685 nm	5-20 ug	92280	CF@660C Features
20-50 ug	92260			
50-100 ug	92239			
CF@680	681/698 nm	5-20 ug	92282	CF@680 Features
20-50 ug	92262			
50-100 ug	92240			
CF@680R	680/701 nm	5-20 ug	92283	CF@680R Features
20-50 ug	92263			
50-100 ug	92246			
CF@700	695/720 nm	5-20 ug	92425	CF@700 Features
20-50 ug	92426			
50-100 ug	92427			
CF@710	712/736 nm	5-20 ug	92579	CF@710 Features
20-50 ug	92580			
50-100 ug	92581			
CF@725	729/750 nm	5-20 ug	92582	CF@725 Features
20-50 ug	92583			
50-100 ug	92584			
CF@740	742/767 nm	5-20 ug	92576	CF@740 Features
20-50 ug	92577			
50-100 ug	92578			
CF@750	755/779 nm	5-20 ug	92284	CF@750 Features
20-50 ug	92264			
50-100 ug	92241			
CF@770	770/797 nm	5-20 ug	92285	CF@770 Features
20-50 ug	92265			
50-100 ug	92242			
CF@790	784/806 nm	5-20 ug	92288	CF@790 Features
20-50 ug	92268			
50-100 ug	92248			
CF@800	797/816 nm	5-20 ug	92428	CF@800 Features
20-50 ug	92429			
50-100 ug	92430			
CF@820	822/835 nm	5-20 ug	92431	CF@820 Features
20-50 ug	92432			
50-100 ug	92433			

* Download the [Product Protocol](#) for a labeling check list to help you select the right kit size and labeling protocol for your antibody. Also see our [Kit Compatibility and Protocol Selection Flowchart](#).

Antibody & Protein Labeling Kits

Product Name	Label Options	Labeling Scale	Labeling Time	Features
Mix-n-Stain™ CF@ Dye Antibody Labeling Kits	CF@ Dyes	≤5-20 ug IgG 20-50 ug IgG 50-100 ug IgG	~ 15 min.	<ul style="list-style-type: none"> • Rapid, simple labeling • No purification • Tolerates BSA & other additives • Dyes for super-resolution, spectral flow, & NIR detection
Mix-n-Stain™ FITC Antibody Labeling Kits	FITC			
Mix-n-Stain™ Cyanine Dye Antibody Labeling Kits	Cyanine 555 Cyanine 647			
Mix-n-Stain™ Biotin Antibody Labeling Kits	Biotin			
Mix-n-Stain™ Digoxigenin Antibody Labeling Kits	Digoxigenin (DIG)			
Mix-n-Stain™ DNP Antibody Labeling Kits	DNP			

Product Name	Label Options	Labeling Scale	Labeling Time	Features
Mix-n-Stain™ Maxi Antibody Labeling Kits	CF® Dyes Cyanine Dyes	1 mg IgG	~ 30 min.	
Mix-n-Stain™ CF® Dye IgM Antibody Labeling Kits	CF® Dyes FITC	25 ug IgM 100 ug IgM	15-30 min.	<ul style="list-style-type: none"> • Rapid, simple labeling • No purification
Mix-n-Stain™ STORM CF® Dye Antibody Labeling Kits	CF® Dyes	50 ug IgG	~ 30 min.	<ul style="list-style-type: none"> • Optimized to provide low 1-2.5 DOL • No purification
Mix-n-Stain™ Nanobody Labeling Kits	CF® Dyes	5-20 ug Nanobody® 20-50 ug Nanobody®	~ 30 min.	<ul style="list-style-type: none"> • Optimized for Nanobodies® • Tolerates BSA & other additives
Mix-n-Stain™ Nanobody Thiol Labeling Kits	CF® Dyes	~ 2 hours	• Optimized for Nanobodies® with single exposed cysteine residue	
Mix-n-Stain™ Small Ligand Labeling Kits	CF® Dyes	0.1 umol small ligand	~ 30 min.	<ul style="list-style-type: none"> • Label SNAP®, CLIP™, HaloTag® & other small ligands • Dye options for surface or intracellular targets
VivoBrite™ Rapid Antibody Labeling Kits for Small Animal In Vivo Imaging	Near-IR CF® Dyes	3 x 1 mg IgG	~ 2 hours	<ul style="list-style-type: none"> • Dyes, buffers, & spin vials for labeling + purification • Sterile syringes, filters and storage vials for in vivo use
CF® Dye & Biotin SE Protein Labeling Kits	CF® Dyes Biotin	3 x 1 mg protein	~ 2 hours	<ul style="list-style-type: none"> • Dyes, buffers, & spin vials for labeling + purification

Alexa Fluor and DyLight are registered trademarks of Thermo Fisher Scientific.

References

Download a list of [CF® Dye and Mix-n-Stain™ references](#).

This datasheet was generated on June 5, 2026 at 01:55:40 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/mix-n-stain-cf-dye-antibody-labeling-kit/>