



Glowing products for science

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

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

Label your antibody with one of our CF® Dyes specialized for STORM super-resolution microscopy. Labeling takes just 30 minutes without a purification step, and yields an antibody conjugate with a low degree of labeling optimal for STORM.



Product attributes

Product Description

Mix-n-Stain™ STORM CF® Dye Antibody Labeling Kits allow you to label 50 ug of your antibody with one of Biotium's STORM CF® Dyes to produce an antibody conjugate with a low 1-2.5 DOL (degree of labeling) optimal for super-resolution STORM (Stochastic Optical Reconstruction Microscopy). Labeling takes just 30 minutes, with minimal hands-on time and no purification.

Superior CF® Dyes for STORM Imaging

These kits feature a choice of 8 CF® Dyes developed with superior brightness and unique photoswitching properties that are ideal for super-resolution imaging by STORM. In addition, these specialized Mix-n-Stain™ STORM CF® Dye Antibody Labeling Kits are designed to offer a low degree of labeling (DOL, or dye molecules per antibody molecule) of 1-2.5, which is reported to be optimal for STORM [\(1\)](#). Learn more about [CF® Dyes for STORM](#) or our [other fluorescent reagents validated for super-resolution microscopy](#).

Discover More with Green-Excitable CF® Dyes Developed for Multi-Color STORM

High-quality STORM imaging relies heavily on bright fluorescent dyes with excellent photoswitching properties. While there are several commonly used red-excited dyes suitable for STORM imaging such as CF®647, CF®660C and CF®680, generating comparable dyes for other excitation sources to produce quality multi-color STORM images remains a challenge. Biotium scientists have met this challenge by developing novel chemical modifications for green-excited fluorescent dyes to offer optimal performance for STORM super-resolution imaging. This includes CF®583R and CF®597R which were developed specifically for STORM imaging and display exceptional performance for multi-color STORM [\(2\)](#).

See the product table below for a quick reference of the maximum excitation and emission of each CF® Dye. Select the links in the Dye Features column to learn more about each CF® Dye and potential applications.

Other Labeling Kits

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

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

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF®505	505/519 nm	50 ug	92549	
CF®568	562/584 nm	50 ug	92550	CF®568 Features
CF®583R	585/609 nm	50 ug	92551	CF®583R Features
CF®597R	597/619 nm	50 ug	92553	CF®597R Features
CF®647	652/668 nm	50 ug	92554	CF®647 Features
CF®660C	667/685 nm	50 ug	92555	CF®660C Features
CF®680	681/698 nm	50 ug	92556	CF®680 Features
CF®750	755/779 nm	50 ug	92557	CF®750 Features

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

Product Name	Label Options	Labeling Scale	Labeling Time	Features
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			
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

CF is a registered trademark of Biotium, Inc.

References

1. Proc. SPIE 9331, (2015), DOI: 10.1117/12.2083209
2. Angew. Chem. Int. Ed. 9331, e202113612, (2021), DOI: 10.1002/ANIE.202113612

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

This datasheet was generated on December 29, 2025 at 10:57:20 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/mix-n-stain-storm-cf-dye-antibody-labeling-kits/>