



Glowing products for science

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

## Mix-n-Stain™ Maxi Antibody Labeling Kits, 1 mg Labeling



The simplicity and convenience of Mix-n-Stain™ labeling, for conjugation of 1 mg IgG with our CF® Dyes or other dyes.

### Product attributes

Labeling scale	1 mg
Purification method	Purification not required after labeling

### Product Description

Mix-n-Stain™ Maxi Antibody Labeling Kits allow you to rapidly label 1 mg of IgG with one of our next-generation CF® Dyes or other label.

#### Unrivaled Convenience

Mix-n-Stain™ Maxi Antibody Labeling Kits allow you to rapidly label 1 mg of IgG with one of our next-generation CF® Dyes or other label. Simply mix your antibody with the reaction buffer and optimally formulated dye provided, followed by a 30 minute incubation and addition of the provided stop buffer. Any remaining free dye or other label is no longer reactive at the end of the labeling, so the conjugate is ready for staining without further purification. Note the labeling protocol and buffer requirements for the Maxi kits differ from our original micro-scale Mix-n-Stain™ kits, see the [Reaction Compatibility Chart](#) below.

After the reaction, the antibody will be labeled with an average of 4-6 dye molecules per antibody molecule. Mix-n-Stain™ labeling is covalent, so labeled antibodies can be used for multiplex staining without transfer of dyes/labels between antibodies.

#### Superior CF® Dyes

These kits feature a choice of Biotium's next-generation CF® Dyes, which were designed for advantages in brightness and photostability compared to Alexa Fluor®, DyLight®, and other fluorescent dyes. Learn more about [CF® Dyes](#). We also offer kits for labeling with Cyanine 555, Cyanine 647, or fluorescein/FITC. See [Ordering Information](#) below.

#### Choose the Right Kit for Your Application

These kits are optimized for a single labeling reaction with 1 mg of IgG. Mix-n-Stain™ kits are not designed to be used for multiple labeling reactions. For smaller-scale conjugation, see our original [Mix-n-Stain™ CF® Dye Antibody Labeling Kits](#); original Mix-n-Stain™ kits also have greater tolerance for antibody buffer additives, and provide special protocols for labeling antibodies in buffer with BSA or in ascites fluid. The kits are not recommended for labeling IgM antibodies, which may denature under the reaction conditions.

Mix-n-Stain™ kits can be used to label other proteins, however, the degree of labeling (DOL, or number of dye molecules per antibody) may not be optimal, and the kits are not designed to allow for DOL measurement after labeling. For performing reactions to optimize DOL for your antibody or application, we recommend using our [CF® Dye SE Protein Labeling Kits](#). The kits include everything you need to perform controlled conjugation reactions and post-labeling purification, as well as detailed instructions and correction factors for determining DOL for CF® Dyes.

Learn about all of our [Mix-n-Stain™ and protein labeling options](#), or see our [Frequently Asked Questions \(FAQs\)](#) for more information.

#### Reaction Compatibility for Mix-n-Stain™ Maxi Antibody Labeling Kits, 1 mg Labeling

Component	Compatibility
Sodium azide	Compatible
Glycerol ≤ 10%:	Compatible
Glycerol > 10%:	Compatible after ultrafiltration*
Tris ≤20 mM	Compatible
Tris >20 mM	Compatible after ultrafiltration*
Glycine	Compatible after ultrafiltration*
BSA or gelatin	Not compatible; purify antibody**
Serum, cell culture supernatant, or ascites fluid	Not compatible; purify antibody**

\* Ultrafiltration vials are sold separately (catalog no. 22004).

\*\* Protocol/reagents for antibody purification not provided.

### Mix-n-Stain™ Maxi Ordering Information

Conjugation	Ex/Em	Labeling size	Catalog No.	Dye Features
<a href="#">CF@350</a>	347/448 nm	1 mg	<a href="#">92420</a>	<a href="#">CF@350 Features</a>
<a href="#">CF@405S</a>	404/431 nm	1 mg	<a href="#">92421</a>	<a href="#">CF@405S Features</a>
<a href="#">CF@405M</a>	408/452 nm	1 mg	<a href="#">92404</a>	<a href="#">CF@405M Features</a>
<a href="#">CF@405L</a>	395/545 nm	1 mg	<a href="#">92454</a>	<a href="#">CF@405L Features</a>
<a href="#">CF@488A</a>	490/515 nm	1 mg	<a href="#">92405</a>	<a href="#">CF@488A Features</a>
<a href="#">CF@555</a>	555/565 nm	1 mg	<a href="#">92406</a>	<a href="#">CF@555 Features</a>
<a href="#">CF@568</a>	562/583 nm	1 mg	<a href="#">92407</a>	<a href="#">CF@568 Features</a>
<a href="#">CF@594</a>	593/614 nm	1 mg	<a href="#">92408</a>	<a href="#">CF@594 Features</a>
<a href="#">CF@633</a>	630/650 nm	1 mg	<a href="#">92409</a>	<a href="#">CF@633 Features</a>
<a href="#">CF@647</a>	650/665 nm	1 mg	<a href="#">92410</a>	<a href="#">CF@647 Features</a>
<a href="#">CF@680</a>	681/698 nm	1 mg	<a href="#">92422</a>	<a href="#">CF@680 Features</a>
<a href="#">CF@750</a>	755/777 nm	1 mg	<a href="#">92423</a>	<a href="#">CF@750 Features</a>
<a href="#">CF@770</a>	770/797 nm	1 mg	<a href="#">92424</a>	<a href="#">CF@770 Features</a>
<a href="#">Fluorescein (FITC)</a>	494/518 nm	1 mg	<a href="#">92411</a>	
<a href="#">Cyanine 555</a>	555/565 nm	1 mg	<a href="#">92415</a>	
<a href="#">Cyanine 647</a>	650/665 nm	1 mg	<a href="#">92419</a>	

## Antibody & Protein Labeling Kits

Product Name	Label Options	Labeling Scale	Labeling Time	Features
<a href="#">Mix-n-Stain™ CF@ Dye Antibody Labeling Kits</a>	CF@ Dyes	≤5-20 ug IgG 20-50 ug IgG 50-100 ug IgG	~ 15 min.	<ul style="list-style-type: none"> <li>• Rapid, simple labeling</li> <li>• No purification</li> <li>• Tolerates BSA &amp; other additives</li> <li>• Dyes for super-resolution, spectral flow, &amp; NIR detection</li> </ul>
<a href="#">Mix-n-Stain™ FITC Antibody Labeling Kits</a>	FITC			
<a href="#">Mix-n-Stain™ Cyanine Dye Antibody Labeling Kits</a>	Cyanine 555 Cyanine 647			
<a href="#">Mix-n-Stain™ Biotin Antibody Labeling Kits</a>	Biotin			
<a href="#">Mix-n-Stain™ Digoxigenin Antibody Labeling Kits</a>	Digoxigenin (DIG)			
<a href="#">Mix-n-Stain™ DNP Antibody Labeling Kits</a>	DNP			
<a href="#">Mix-n-Stain™ Maxi Antibody Labeling Kits</a>	CF@ Dyes Cyanine Dyes	1 mg IgG	~ 30 min.	
<a href="#">Mix-n-Stain™ CF@ Dye IgM Antibody Labeling Kits</a>	CF@ Dyes FITC	25 ug IgM 100 ug IgM	15-30 min.	<ul style="list-style-type: none"> <li>• Rapid, simple labeling</li> <li>• No purification</li> </ul>
<a href="#">Mix-n-Stain™ STORM CF@ Dye Antibody Labeling Kits</a>	CF@ Dyes	50 ug IgG	~ 30 min.	<ul style="list-style-type: none"> <li>• Optimized to provide low 1-2.5 DOL</li> <li>• No purification</li> </ul>

Product Name	Label Options	Labeling Scale	Labeling Time	Features
<a href="#">Mix-n-Stain™ Nanobody Labeling Kits</a>	CF® Dyes	5-20 ug Nanobody® 20-50 ug Nanobody®	~ 30 min.	<ul style="list-style-type: none"> <li>• Optimized for Nanobodies®</li> <li>• Tolerates BSA &amp; other additives</li> </ul>
<a href="#">Mix-n-Stain™ Nanobody Thiol Labeling Kits</a>	CF® Dyes	~ 2 hours	• Optimized for Nanobodies® with single exposed cysteine residue	
<a href="#">Mix-n-Stain™ Small Ligand Labeling Kits</a>	CF® Dyes	0.1 umol small ligand	~ 30 min.	<ul style="list-style-type: none"> <li>• Label SNAP®, CLIP™, HaloTag® &amp; other small ligands</li> <li>• Dye options for surface or intracellular targets</li> </ul>
<a href="#">VivoBrite™ Rapid Antibody Labeling Kits for Small Animal In Vivo Imaging</a>	Near-IR CF® Dyes	3 x 1 mg IgG	~ 2 hours	<ul style="list-style-type: none"> <li>• Dyes, buffers, &amp; spin vials for labeling + purification</li> <li>• Sterile syringes, filters and storage vials for in vivo use</li> </ul>
<a href="#">CF® Dye &amp; Biotin SE Protein Labeling Kits</a>	CF® Dyes Biotin	3 x 1 mg protein	~ 2 hours	• Dyes, buffers, & spin vials for labeling + purification

Alexa Fluor and DyLight are registered trademarks of Thermo Fisher Scientific; SNAP-tag and CLIP-tag are trademarks or registered trademarks of New England BioLabs; HaloTag is a registered trademark of Promega Corporation.

## References

Download a list of curated [CF® Dye references](#).

This datasheet was generated on June 11, 2026 at 06:44:22 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/mix-n-stain-antibody-labeling-kits-1-mg-labeling/>