

# 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**

Call us: 800-304-5357

### **Product Description**

Mix-n-Stain™ CF® Dye Antibody Labeling Kits allow you to label ≤5 ug up to 100 ug of your lgG 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<sup>TM</sup> 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 <a href="Protocol for a">Protocol for a</a> labeling check list to help you select the right kit size and labeling protocol for your antibody. Also see our <a href="Kit Compatibility and Protocol Selection Flowchart">Kit Compatibility and Protocol Selection Flowchart</a>.

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<sup>™</sup> CF® Dye 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			
<u>CF®405S</u>	404/431 nm	5-20 ug	<u>92271</u>	CF®405S Features
20-50 ug	<u>92251</u>			
50-100 ug	<u>92231</u>			
CF®405M	408/452 nm	5-20 ug	<u>92272</u>	CF®405M Features
20-50 ug	<u>92252</u>			
50-100 ug	92232			
CF®405L	395/545 nm	5-20 ug	<u>92303</u>	CF®405L Features
20-50 ug	<u>92304</u>			
50-100 ug	<u>92305</u>			
3x(5-20 ug)	<u>92445</u>			
<u>CF®430</u>	426/498 nm	5-20 ug	<u>92316</u>	CF®430 Features
20-50 ug	<u>92317</u>			
50-100 ug	<u>92318</u>			
<u>CF®440</u>	440/515 nm	5-20 ug	<u>92319</u>	CF®440 Features
20-50 ug	<u>92320</u>			
50-100 ug	<u>92321</u>	5.00	00000	0594505
CF®450	450/538 nm	5-20 ug	<u>92322</u>	CF®450 Features
20-50 ug	<u>92323</u>			
50-100 ug	92324	F 00	00070	OF@4004 F+::
<u>CF®488A</u>	490/515 nm	5-20 ug	<u>92273</u>	CF®488A Features
20-50 ug	<u>92253</u>			
50-100 ug	<u>92233</u>			
3x(5-20 ug)	92446 F16/F49.pm	E 20 ug	<u>92331</u>	CF®514 Features
<u>CF®514</u>	516/548 nm	5-20 ug	<u>92331</u>	CF®314 Features
20-50 ug 50-100 ug	<u>92332</u> <u>92333</u>			
CF®532	527/558 nm	5-20 ug	<u>92289</u>	CF®532 Features
20-50 ug	92290	5 20 ag	<u>JLLUU</u>	OT GOOD I CALUICS
50-100 ug	<u>92291</u>			
<u>CF®543</u>	541/560 nm	5-20 ug	<u>92287</u>	CF®543 Features
20-50 ug	92267	- <del></del>		
50-100 ug	<u>92247</u>			
<u>CF®555</u>	555/565 nm	5-20 ug	92274	CF®555 Features
20-50 ug	92254	Ÿ		
50-100 ug	92234			
CF®568	562/583 nm	5-20 ug	<u>92275</u>	CF®568 Features
20-50 ug	<u>92255</u>			
50-100 ug	<u>92235</u>			
3x(5-20 ug)	92447			
CF®570	568/591 nm	5-20 ug	<u>92334</u>	CF®570 Features
20-50 ug	<u>92335</u>			
50-100 ug	<u>92336</u>			
<u>CF®583</u>	583/606 nm	5-20 ug	92337	CF®583 Features
20-50 ug	92338			
50-100 ug	92339			
<u>CF®583R</u>	586/609 nm	5-20 ug	<u>92455</u>	CF®583R Features
20-50 ug	<u>92456</u>			
50-100 ug	<u>92457</u>			
CF®594	593/614 nm	5-20 ug	<u>92276</u>	CF®594 Features
20-50 ug	<u>92256</u>			
50-100 ug	<u>92236</u>			
3x(5-20 ug)	92448 507/010 mm	F 00	00404	OF@F07D F+
<u>CF®597R</u>	597/619 nm	5-20 ug	<u>92461</u>	CF®597R Features
20-50 ug	<u>92462</u>			
50-100 ug	92463 620/650 pm	5.20 ug	02277	CE@622 Footures
<u>CF®633</u>	630/650 nm	5-20 ug	<u>92277</u>	CF®633 Features
20-50 ug	<u>92257</u> <u>92237</u>			
50-100 ug CF®640R	92237 642/662 nm	5-20 ug	<u>92278</u>	CF®640R Features
20-50 ug	92258	J-ZU uy	<u>32210</u>	OFWOHUN FEATURES
50-100 ug	<u>92236</u> <u>92245</u>			
CF®647	650/665 nm	5-20 ug	<u>92279</u>	CF®647 Features
<u>5. 8071</u>	GOO/GOO HIII	5 20 ag	<u>JLLI J</u>	OT GOTT I GALAIGS

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

<sup>\*</sup> Download the <u>Product Protocol</u> for a labeling check list to help you select the right kit size and labeling protocol for your antibody. Also see our <u>Kit Compatibility and Protocol Selection Flowchart</u>.

# **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	~ 15 min.	Rapid, simple labeling
		50-100 ug lgG		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 <u>Dye</u> Antibody Labeling Kits	Cyanine 555 Cyanine 647			
Mix-n-Stain™ Biotin Antibody Labeling Kits	Biotin			
Mix-n-Stain™ Digoxigenin Antibody Labeling Kits	Digoxygenin (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 lgG	~ 30 min.	
Mix-n-Stain™ CF® Dye IaM Antibody Labeling	CF® Dyes FITC	25 ug lgM 100 ug lgM	15-30 min.	<ul> <li>Rapid, simple labeling</li> </ul>
Kits	1110	100 ug igivi		<ul> <li>No purification</li> </ul>
Mix-n-Stain™ STORM CF® Dye Antibody Labeling Kits	CF® Dyes	50 ug IgG	~ 30 min.	Optimized to provide low 1-2.5 DOL
Mix-n-Stain™	CF® Dyes	5-20 ug Nanobody®	~ 30 min.	<ul><li>No purification</li><li>Optimized for</li></ul>
Nanobody Labeling Kits	or & byco	20-50 ug Nanobody®	00 mm.	Nanobodies®
				<ul> <li>Tolerates BSA &amp; other additives</li> </ul>
<u>Mix-n-Stain™</u> <u>Nanobody Thiol Labeling</u> <u>Kits</u>	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> <li>Label SNAP®, CLIP™, HaloTag® &amp; other small ligands</li> </ul>
				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 lgG	~ 2 hours	Dyes, buffers, & spin vials for labeling + purification
ımayıny				<ul> <li>Sterile syringes, filters and storage vials for in vivo use</li> </ul>
CF® Dye & Biotin SE Protein Labeling Kits	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.

## References

Download a list of  $\underline{\mathsf{CF}}$  Dye and  $\underline{\mathsf{Mix}}$ -n-Stain $^{\mathsf{TM}}$  references.

This datasheet was generated on November 21, 2025 at 01:16:20 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/mix-n-stain-cf-dye-antibody-labeling-kit/">https://biotium.com/product/mix-n-stain-cf-dye-antibody-labeling-kit/</a>