



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357) Email: [btinfo@biotium.com](mailto:btinfo@biotium.com)

## 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  $\leq 5$  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
<a href="#">CF@350</a>	347/448 nm	5-20 ug	<a href="#">92270</a>	<a href="#">CF@350 Features</a>
		20-50 ug	<a href="#">92250</a>	
		50-100 ug	<a href="#">92230</a>	
<a href="#">CF@405S</a>	404/431 nm	5-20 ug	<a href="#">92271</a>	<a href="#">CF@405S Features</a>
		20-50 ug	<a href="#">92251</a>	
		50-100 ug	<a href="#">92231</a>	
<a href="#">CF@405M</a>	408/452 nm	5-20 ug	<a href="#">92272</a>	<a href="#">CF@405M Features</a>
		20-50 ug	<a href="#">92252</a>	
		50-100 ug	<a href="#">92232</a>	
<a href="#">CF@405L</a>	395/545 nm	5-20 ug	<a href="#">92303</a>	<a href="#">CF@405L Features</a>
		20-50 ug	<a href="#">92304</a>	
		50-100 ug	<a href="#">92305</a>	
<a href="#">CF@430</a>	426/498 nm	3x(5-20 ug)	<a href="#">92445</a>	<a href="#">CF@430 Features</a>
		5-20 ug	<a href="#">92316</a>	
		20-50 ug	<a href="#">92317</a>	
		50-100 ug	<a href="#">92318</a>	
<a href="#">CF@440</a>	440/515 nm	5-20 ug	<a href="#">92319</a>	<a href="#">CF@440 Features</a>
		20-50 ug	<a href="#">92320</a>	
		50-100 ug	<a href="#">92321</a>	
<a href="#">CF@450</a>	450/538 nm	5-20 ug	<a href="#">92322</a>	<a href="#">CF@450 Features</a>
		20-50 ug	<a href="#">92323</a>	
		50-100 ug	<a href="#">92324</a>	
<a href="#">CF@488A</a>	490/515 nm	5-20 ug	<a href="#">92273</a>	<a href="#">CF@488A Features</a>
		20-50 ug	<a href="#">92253</a>	
		50-100 ug	<a href="#">92233</a>	
		3x(5-20 ug)	<a href="#">92446</a>	
<a href="#">CF@514</a>	516/548 nm	5-20 ug	<a href="#">92331</a>	<a href="#">CF@514 Features</a>
		20-50 ug	<a href="#">92332</a>	
		50-100 ug	<a href="#">92333</a>	
<a href="#">CF@532</a>	527/558 nm	5-20 ug	<a href="#">92289</a>	<a href="#">CF@532 Features</a>
		20-50 ug	<a href="#">92290</a>	
		50-100 ug	<a href="#">92291</a>	
<a href="#">CF@543</a>	541/560 nm	5-20 ug	<a href="#">92287</a>	<a href="#">CF@543 Features</a>
		20-50 ug	<a href="#">92267</a>	
		50-100 ug	<a href="#">92247</a>	
<a href="#">CF@555</a>	555/565 nm	5-20 ug	<a href="#">92274</a>	<a href="#">CF@555 Features</a>
		20-50 ug	<a href="#">92254</a>	
		50-100 ug	<a href="#">92234</a>	
<a href="#">CF@568</a>	562/583 nm	5-20 ug	<a href="#">92275</a>	<a href="#">CF@568 Features</a>
		20-50 ug	<a href="#">92255</a>	
		50-100 ug	<a href="#">92235</a>	
		3x(5-20 ug)	<a href="#">92447</a>	
<a href="#">CF@570</a>	568/591 nm	5-20 ug	<a href="#">92334</a>	<a href="#">CF@570 Features</a>
		20-50 ug	<a href="#">92335</a>	
		50-100 ug	<a href="#">92336</a>	
<a href="#">CF@583</a>	583/606 nm	5-20 ug	<a href="#">92337</a>	<a href="#">CF@583 Features</a>
		20-50 ug	<a href="#">92338</a>	
		50-100 ug	<a href="#">92339</a>	
<a href="#">CF@583R</a>	586/609 nm	5-20 ug	<a href="#">92455</a>	<a href="#">CF@583R Features</a>
		20-50 ug	<a href="#">92456</a>	
		50-100 ug	<a href="#">92457</a>	
<a href="#">CF@594</a>	593/614 nm	5-20 ug	<a href="#">92276</a>	<a href="#">CF@594 Features</a>
		20-50 ug	<a href="#">92256</a>	
		50-100 ug	<a href="#">92236</a>	
		3x(5-20 ug)	<a href="#">92448</a>	
<a href="#">CF@597R</a>	597/619 nm	5-20 ug	<a href="#">92461</a>	<a href="#">CF@597R Features</a>
		20-50 ug	<a href="#">92462</a>	
		50-100 ug	<a href="#">92463</a>	
<a href="#">CF@633</a>	630/650 nm	5-20 ug	<a href="#">92277</a>	<a href="#">CF@633 Features</a>
		20-50 ug	<a href="#">92257</a>	
		50-100 ug	<a href="#">92237</a>	
<a href="#">CF@640R</a>	642/662 nm	5-20 ug	<a href="#">92278</a>	<a href="#">CF@640R Features</a>
		20-50 ug	<a href="#">92258</a>	
		50-100 ug	<a href="#">92245</a>	

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
<a href="#">CF@647</a>	650/665 nm	5-20 ug	<a href="#">92279</a>	<a href="#">CF@647 Features</a>
		20-50 ug	<a href="#">92259</a>	
		50-100 ug	<a href="#">92238</a>	
		3x(5-20 ug)	<a href="#">92449</a>	
<a href="#">CF@660R</a>	663/682 nm	5-20 ug	<a href="#">92281</a>	<a href="#">CF@660R Features</a>
		20-50 ug	<a href="#">92261</a>	
		50-100 ug	<a href="#">92243</a>	
<a href="#">CF@660C</a>	667/685 nm	5-20 ug	<a href="#">92280</a>	<a href="#">CF@660C Features</a>
		20-50 ug	<a href="#">92260</a>	
		50-100 ug	<a href="#">92239</a>	
<a href="#">CF@680</a>	681/698 nm	5-20 ug	<a href="#">92282</a>	<a href="#">CF@680 Features</a>
		20-50 ug	<a href="#">92262</a>	
		50-100 ug	<a href="#">92240</a>	
<a href="#">CF@680R</a>	680/701 nm	5-20 ug	<a href="#">92283</a>	<a href="#">CF@680R Features</a>
		20-50 ug	<a href="#">92263</a>	
		50-100 ug	<a href="#">92246</a>	
<a href="#">CF@700</a>	695/720 nm	5-20 ug	<a href="#">92425</a>	<a href="#">CF@700 Features</a>
		20-50 ug	<a href="#">92426</a>	
		50-100 ug	<a href="#">92427</a>	
<a href="#">CF@710</a>	712/736 nm	5-20 ug	<a href="#">92579</a>	<a href="#">CF@710 Features</a>
		20-50 ug	<a href="#">92580</a>	
		50-100 ug	<a href="#">92581</a>	
<a href="#">CF@725</a>	729/750 nm	5-20 ug	<a href="#">92582</a>	<a href="#">CF@725 Features</a>
		20-50 ug	<a href="#">92583</a>	
		50-100 ug	<a href="#">92584</a>	
<a href="#">CF@740</a>	742/767 nm	5-20 ug	<a href="#">92576</a>	<a href="#">CF@740 Features</a>
		20-50 ug	<a href="#">92577</a>	
		50-100 ug	<a href="#">92578</a>	
<a href="#">CF@750</a>	755/779 nm	5-20 ug	<a href="#">92284</a>	<a href="#">CF@750 Features</a>
		20-50 ug	<a href="#">92264</a>	
		50-100 ug	<a href="#">92241</a>	
<a href="#">CF@770</a>	770/797 nm	5-20 ug	<a href="#">92285</a>	<a href="#">CF@770 Features</a>
		20-50 ug	<a href="#">92265</a>	
		50-100 ug	<a href="#">92242</a>	
<a href="#">CF@790</a>	784/806 nm	5-20 ug	<a href="#">92288</a>	<a href="#">CF@790 Features</a>
		20-50 ug	<a href="#">92268</a>	
		50-100 ug	<a href="#">92248</a>	
<a href="#">CF@800</a>	797/816 nm	5-20 ug	<a href="#">92428</a>	<a href="#">CF@800 Features</a>
		20-50 ug	<a href="#">92429</a>	
		50-100 ug	<a href="#">92430</a>	
<a href="#">CF@820</a>	822/835 nm	5-20 ug	<a href="#">92431</a>	<a href="#">CF@820 Features</a>
		20-50 ug	<a href="#">92432</a>	
		50-100 ug	<a href="#">92433</a>	

\* 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
<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> </ul>
<a href="#">Mix-n-Stain™ FITC Antibody Labeling Kits</a>	FITC			<ul style="list-style-type: none"> <li>• Tolerates BSA &amp; other additives</li> </ul>
<a href="#">Mix-n-Stain™ Cyanine Dye Antibody Labeling Kits</a>	Cyanine 555 Cyanine 647			<ul style="list-style-type: none"> <li>• Dyes for super-resolution, spectral flow, &amp; NIR detection</li> </ul>
<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>

Product Name	Label Options	Labeling Scale	Labeling Time	Features
<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>
<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™ 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	<ul style="list-style-type: none"> <li>• Dyes, buffers, &amp; spin vials for labeling + purification</li> </ul>

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 April 1, 2025 at 11:46:00 PM. Visit product page to check for updated information before use.

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