# **CF® Dye & Biotin SE Protein Labeling Kits**

#### **Product attributes**

Call us: 800-304-5357

Kits for succinimidyl ester labeling antibodies or other proteins with Biotium's next-generation CF® Dyes or biotin.  $3 \times 1$  mg labeling reactions per kit.

#### **Product Description**

This kit contains reagents for labeling antibodies or other proteins with a fluorescent CF® dye succinimidyl ester (CF® dye SE) or biotin succinimidyl ester (biotin SE). Learn more about next-generation CF® dyes with superior brightness, photostability, and water solubility compared to other fluorescent dyes.

**Note:** Conjugates of blue-fluorescent dyes like CF®350, CF®405S and CF®405M are not recommended for detecting low abundance targets and may be challenging to use in tissue specimens. Blue dyes have lower fluorescence and photostability, and cells and tissue have high autofluorescence in blue wavelengths, resulting in lower signal to noise compared to other colors.

See our full line of Antibody & Protein Labeling Kits.

### **CF® Dye & Biotin SE Protein Labeling Kits**

SE Protein Labeling Kit	Ex/Em	Size	Catalog No.	Dye Features
CF®350	347/448 nm	3 x 1 mg labelings	92210	CF®350 Features
CF®405S	404/431 nm	3 x 1 mg labelings	<u>92211</u>	CF®405S Features
CF®405M	408/452 nm	3 x 1 mg labelings	<u>92212</u>	CF®405M Features
CF®405L	395/545 nm	3 x 1 mg labelings	<u>92228</u>	CF®405L Features
<u>CF®488A</u>	490/515 nm	3 x 1 mg labelings	<u>92213</u>	CF®488A Features
<u>CF®532</u>	527/558 nm	3 x 1 mg labelings	<u>92208</u>	CF®532 Features
CF®543	541/560 nm	3 x 1 mg labelings	<u>92209</u>	CF®543 Features
<u>CF®555</u>	555/565 nm	3 x 1 mg labelings	<u>92214</u>	CF®555 Features
<u>CF®568</u>	562/583 nm	3 x 1 mg labelings	<u>92215</u>	CF®568 Features
CF®594	593/614 nm	3 x 1 mg labelings	<u>92216</u>	CF®594 Features
CF®633	630/650 nm	3 x 1 mg labelings	<u>92217</u>	CF®633 Features
<u>CF®640R</u>	642/662 nm	3 x 1 mg labelings	<u>92225</u>	CF®640R Features
CF®647	650/665 nm	3 x 1 mg labelings	<u>92218</u>	CF®647 Features
<u>CF®660R</u>	663/682 nm	3 x 1 mg labelings	<u>92219</u>	CF®660R Features
<u>CF®660C</u>	667/685 nm	3 x 1 mg labelings	<u>92223</u>	CF®660C Features
CF®680	681/698 nm	3 x 1 mg labelings	<u>92220</u>	CF®680 Features
<u>CF®680R</u>	680/701 nm	3 x 1 mg labelings	<u>92226</u>	CF®680R Features
CF®750	755/777 nm	3 x 1 mg labelings	<u>92221</u>	CF®750 Features
CF®770	770/797 nm	3 x 1 mg labelings	<u>92222</u>	CF®770 Features
<u>Biotin</u>	N/A	3 x 1 mg labelings	92224	

## **Antibody & Protein Labeling Kits**

**Biotin** 

Antibody Labeling Kits
Mix-n-Stain<sup>TM</sup> Biotin

Antibody Labeling Kits

-	_			
Product Name	Label Options	Labeling Scale	Labeling Time	Features
Mix-n-Stain™ CF® Dye Antibody Labeling Kits	CF® Dyes	≤5-20 ug lgG 20-50 ug lgG 50-100 ug lgG	~ 15 min.	Rapid, simple labeling
				<ul> <li>No purification</li> </ul>
				<ul> <li>Tolerates BSA &amp; other additives</li> </ul>
				Dyes for super-resolution, spectral flow, & NIR detection
Mix-n-Stain™ FITC Antibody Labeling Kits	FITC			
Mix-n-Stain™ Cyanine Dye	Cyanine 555 Cyanine 647			

Product Name	Label Options	Labeling Scale	Labeling Time	Features
Mix-n-Stain™ Digoxigenin Antibody Labeling Kits	Digoxygenin (DIG)			
Mix-n-Stain™ DNP Antibody Labeling Kits	DNP			
Mix-n-Stain™ Maxi Antibody Labeling Kits	CF® Dyes Cyanine Dyes	1 mg lgG	~ 30 min.	
Mix-n-Stain™ CF® Dye IgM Antibody Labeling	CF® Dyes FITC	25 ug IgM 100 ug IgM	15-30 min.	Rapid, simple labeling
Kits				No purification
Mix-n-Stain™ STORM CF® Dye Antibody Labeling Kits	CF® Dyes	50 ug lgG	~ 30 min.	Optimized to provide low     1-2.5 DOL
M' OL TH	05e D	500 N L L	00 '	No purification
Mix-n-Stain™ Nanobody Labeling Kits	CF® Dyes	5-20 ug Nanobody® 20-50 ug Nanobody®	~ 30 min.	Optimized for Nanobodies®
				<ul> <li>Tolerates BSA &amp; other additives</li> </ul>
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> <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
magity				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	Dyes, buffers, & spin vials for labeling + purification

#### References

Download a list of  $\underline{\mathsf{CF}}\underline{\mathsf{B}}$  dye references.

This datasheet was generated on November 20, 2025 at 05:32:15 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/cf-dye-biotin-se-protein-labeling-kits/">https://biotium.com/product/cf-dye-biotin-se-protein-labeling-kits/</a>