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

Label your IgM antibody with one of our superior CF® Dyes in 15-30 minutes without a purification step.



Product attributes

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

Mix-n-Stain™ CF® Dye IgM Antibody Labeling Kits allow you to label 25 ug or 100 ug of your IgM antibody in 15-30 minutes, with minimal hands-on time and no purification.

Finally! Simple & Efficient Labeling of IgM Antibodies

IgM antibodies are difficult to label with fluorescent dyes because the conditions typically used for labeling IgG generally do not result in bright conjugates with good antibody reactivity for IgM. Biotium has resolved this issue by optimizing the reaction conditions for labeling IgM antibodies with Mix-n-StainTM CF® Dye IgM Antibody Labeling Kits.

Simply mix your purified IgM antibody with the reaction buffer and pre-measured dye provided, followed by a brief incubation. Any free dye or label in solution is no longer reactive at the end of the labeling, so the conjugate is ready for staining without further purification. The resulting conjugate is produced with optimal degree of labeling (DOL) for cell staining. Mix-n-StainTM labeling is covalent, therefore labeled antibodies can be used for multiplex staining without the transfer of dye between proteins.

Superior CF® Dyes

These kits feature a choice of 8 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-StainTM CF® Dye IgM Antibody Labeling Kits are very simple to use. But before you begin, you must check that your antibody meets the compatibility requirements for labeling, and choose the right labeling protocol. Download the Protocol to help you determine if your IgM antibody is compatible with the labeling protocol, or if it requires further purification.

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

Check out our main line of Mix-n-Stain™ CF® Dye Antibody Labeling Kits and Mix-n-Stain™ Maxi Antibody Labeling Kits designed for rapid and efficient labeling of IgG antibodies. We also offer Mix-n-Stain™ Nanobody Labeling Kits for labeling single-domain Nanobodies®. 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 IgM Antibody Labeling Kits

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF®405L	413/547 nm	25 ug	<u>92558</u>	CF®405L Features
		100 ug	<u>92559</u>	
<u>CF®405M</u>	416/452 nm	25 ug	<u>92560</u>	CF®405M Features
		100 ug	<u>92561</u>	
<u>CF®488A</u>	490/516 nm	25 ug	<u>92562</u>	CF®488A Features
		100 ug	<u>92563</u>	
<u>CF®555</u>	554/568 nm	25 ug	<u>92564</u>	CF®555 Features
		100 ug	<u>92565</u>	
<u>CF®568</u>	562/583 nm	25 ug	<u>92566</u>	CF®568 Features
		100 ug	<u>92567</u>	
<u>CF®594</u>	593/615 nm	25 ug	<u>92568</u>	CF®594 Features
		100 ug	<u>92569</u>	
<u>CF®640R</u>	642/663 nm	25 ug	<u>92570</u>	CF®640R Features
		100 ug	<u>92571</u>	
<u>CF®647</u>	652/668 nm	25 ug	<u>92572</u>	CF®647 Features
		100 ug	<u>92573</u>	
FITC	498/517 nm	25 ug	<u>92574</u>	
		100 ug	<u>92575</u>	

Note: Before selecting a kit, please download the Product Protocol to see Kit Compatibility and Protocol Selection guidelines.

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 lgG 20-50 ug lgG 50-100 ug lgG	~ 15 min.	Rapid, simple labeling
				 No purification
Mix-n-Stain™ FITC Antibody Labeling Kits	FITC			 Tolerates BSA & other additives
Mix-n-Stain™ Cyanine Dye Antibody Labeling Kits	Cyanine 555 Cyanine 647			Dyes for super-resolution, spectral
Mix-n-Stain™ Biotin Antibody Labeling Kits	Biotin			flow, & NIR detection
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 Kits	CF® Dyes FITC	25 ug lgM 100 ug lgM	15-30 min.	Rapid, simple labeling
				No purification
Mix-n-Stain™ STORM CF® Dye Antibody Labeling Kits	CF® Dyes	50 ug IgG	~ 30 min.	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.	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.	 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 lgG	~ 2 hours	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	 Dyes, buffers, & spin vials for labeling + purification

Nanobody is a registered trademark of Ablynx, a Sanofi company.

References

Download a list of curated CF® Dye references.

This datasheet was generated on November 12, 2025 at 07:08:27 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/mix-n-stain-cf-dye-igm-antibody-labeling-kits/