

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

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-Stain™ 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-Stain™ 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-Stain™ 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 [Product 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 100 ug	413/547 nm 92559	25 ug	92558	CF@405L Features
CF@405M 100 ug	416/452 nm 92561	25 ug	92560	CF@405M Features
CF@488A 100 ug	490/516 nm 92563	25 ug	92562	CF@488A Features
CF@555 100 ug	554/568 nm 92565	25 ug	92564	CF@555 Features
CF@568 100 ug	562/583 nm 92567	25 ug	92566	CF@568 Features
CF@594 100 ug	593/615 nm 92569	25 ug	92568	CF@594 Features
CF@640R 100 ug	642/663 nm 92571	25 ug	92570	CF@640R Features
CF@647 100 ug	652/668 nm 92573	25 ug	92572	CF@647 Features
FITC 100 ug	498/517 nm 92575	25 ug	92574	

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 IgG 20-50 ug IgG 50-100 ug IgG	~ 15 min.	<ul style="list-style-type: none"> • Rapid, simple labeling • 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 Dye Antibody Labeling Kits	Cyanine 555 Cyanine 647			
Mix-n-Stain™ Biotin Antibody Labeling Kits	Biotin			
Mix-n-Stain™ Digoxigenin Antibody Labeling Kits	Digoxigenin (DIG)			
Mix-n-Stain™ DNP Antibody Labeling Kits	DNP			
Mix-n-Stain™ Maxi Antibody Labeling Kits	CF@ Dyes Cyanine Dyes	1 mg IgG	~ 30 min.	
Mix-n-Stain™ CF@ Dye IgM Antibody Labeling Kits	CF@ Dyes FITC	25 ug IgM 100 ug IgM	15-30 min.	<ul style="list-style-type: none"> • Rapid, simple labeling • No purification
Mix-n-Stain™ STORM CF@ Dye Antibody Labeling Kits	CF@ Dyes	50 ug IgG	~ 30 min.	<ul style="list-style-type: none"> • Optimized to provide low 1-2.5 DOL • No purification • Optimized for Nanobodies® • Tolerates BSA & other additives
Mix-n-Stain™ Nanobody Labeling Kits	CF@ Dyes	5-20 ug Nanobody® 20-50 ug Nanobody®	~ 30 min.	<ul style="list-style-type: none"> • Optimized for Nanobodies® with single exposed cysteine residue
Mix-n-Stain™ Nanobody Thiol Labeling Kits	CF@ Dyes	~ 2 hours		
Mix-n-Stain™ Small Ligand Labeling Kits	CF@ Dyes	0.1 umol small ligand	~ 30 min.	<ul style="list-style-type: none"> • 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 IgG	~ 2 hours	<ul style="list-style-type: none"> • 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	<ul style="list-style-type: none"> • 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 June 5, 2026 at 08:16:02 PM. 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/>