

Product Information

Mix-n-Stain™ CF™405M Antibody Labeling Kits

Catalog Numbers **MX405MS20**, **MX405MS50**, and **MX405MS100**

Storage Temperature -20°C

TECHNICAL BULLETIN

Product Description

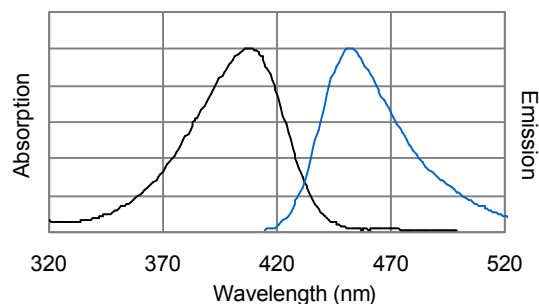
The kits contain everything needed to rapidly label an antibody. Select the labeling kit suitable for the amount of antibody to be labeled. The labeling procedure is simply mixing of the antibody with the optimally formulated dye in the reaction buffer provided, followed by a brief incubation. The resulting solution is ready for staining without further purification. The kit is suitable for labeling commercially available primary antibodies, either directly or after a simple antibody clean-up step. After labeling, the dye is covalently linked to the antibody with a degree of labeling of 4–6 dye molecules per antibody molecule. Multiple antibodies labeled with different dyes using Mix-n-Stain™ kits can be used for multicolor fluorescence staining without transfer of dyes between antibodies.

CF™405M is a blue fluorescent dye spectrally similar to Pacific Blue™ dye. CF405M is as bright as Pacific Blue dye but is significantly more photostable. An additional advantage of CF405M is it has less spillover in the 525/50 green channel than Pacific Blue dye due to narrower emission peak, making CF405M a superior choice in multicolor detection applications. CF405S is also offered, another 405 nm-excitable dye with an even shorter emission wavelength.

Note: Labeled secondary antibodies will bind to primary antibodies labeled using Mix-n-Stain kits; therefore, if multiple primary antibodies from the same species are to be used for multicolor immunofluorescence staining, a secondary antibody cannot be used to distinguish an unlabeled primary antibody from a Mix-n-Stain labeled primary antibody. Biotin Mix-n-Stain labeling kits are offered for secondary detection using CF dye-labeled streptavidin or CF dye-labeled monoclonal mouse anti-biotin.

Figure 1.

Absorption and emission spectra of CF405M conjugated to goat anti-mouse IgG in PBS.



CF405M dye properties:

Abs/Em Maxima: 408/450 nm

Flow cytometry laser line: 360 nm, 405 nm, or 407 nm

Microscopy laser line: 405 nm

Direct replacement for: BD Horizon™ V450, eFluor® 450, and Pacific Blue

Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

Storage/Stability

Store the kit at -20°C .

Procedure

Procedures

A. Antibody Preparation

Check the compatibility of the antibody formulation with the compatibility table (see Table 1). An antibody solution free of stabilizers produces better labeling results. However, low levels of stabilizers can be tolerated in the standard protocol.

Non-protein stabilizers can be removed using the ultrafiltration vial provided.

Table 1.

Mix-n-Stain Compatibility with Common Antibody Storage Components

Component	Compatibility
Sodium Azide	compatible
Glycerol	≤10%: proceed to Procedure C >10%: proceed to Procedure B
Tris	≤20 mM: proceed to Procedure C >20 mM: proceed to Procedure B
Glycine	Proceed to Procedure B
BSA or gelatin	≤4× IgG by weight: proceed to Procedure C >4× IgG by weight: proceed to Procedure D
Ascites fluid	Proceed to Procedure D
Serum	Not compatible; purify IgG
Hybridoma supernatant	Not compatible; purify IgG

Antibodies can be labeled in the presence of excess protein stabilizers or in ascites fluid by selecting a kit size that matches the amount of total protein in the labeling reaction (IgG plus stabilizer or amount of protein in ascites fluid) and using the modified Mix-n-Stain labeling procedure (Procedure D). The modified Mix-n-Stain protocol can be used to label antibody amounts that fall below the range of the kit by adding stabilizer protein to the antibody to bring the total protein amount within the kit range (see Procedure C).

One may experience higher background after staining if the antibody is labeled in the presence of BSA or gelatin. Background staining can be reduced by using blocking and wash solutions containing at least 1% BSA or gelatin, respectively. Alternatively, protein stabilizers (BSA or gelatin) can be removed using an antibody clean-up kit.

For optimal labeling, use the antibody at a concentration of 0.5–1 mg/mL. If necessary, adjust the antibody concentration with 1× PBS. The ultrafiltration protocol (Procedure B) can be performed to concentrate more dilute solutions of antibody.

B. Ultrafiltration

Notes: Before beginning, it is important to use Table 1 to determine whether the antibody requires ultrafiltration before labeling. If ultrafiltration is not required, proceed to the appropriate labeling procedure as determined by Table 1.

The ultrafiltration column membrane has a molecular mass cut-off of 10,000 Da. Therefore, molecules smaller than 10 kDa will flow through the membrane and molecules larger than 10 kDa, including IgG antibodies, will be retained on the upper surface of the membrane. Take care not to touch the membrane with pipette tips, which could tear or puncture the membrane, resulting in loss of antibody.

Ultrafiltration vial capacities:

Maximum sample volume: 500 µL

Final concentrate volume: 15 µL

Filtrate receiver volume: 500 µL

Hold-up volume (membrane/support): <5 µL

1. Add an appropriate amount of antibody to the ultrafiltration vial being careful not to touch the membrane. Spin the solution at $14,000 \times g$ in a microcentrifuge for a few minutes until the liquid is removed. Discard the liquid in the collection vessel.
2. Rinse the original antibody vial with an equal volume of 1× PBS and add to the ultrafiltration vial. Spin the vial at $14,000 \times g$ until the liquid is removed. Resuspend the antibody to a concentration of 0.5–1 mg/mL in an appropriate volume of 1× PBS and transfer to a clean vial.
3. If using the modified Mix-n-Stain labeling procedure (Procedure D), save the ultrafiltration vial to concentrate the antibody after labeling.

C. Standard Mix-n-Stain Labeling

Note: Before beginning, use Table 1 to select the appropriate labeling protocol for the antibody.

1. Warm up the Mix-n-Stain Reaction Buffer vial and the Mix-n-Stain Storage Buffer vial to room temperature before use. Centrifuge the vials briefly to collect the solutions at the bottom of the vials.
2. Dilute the 10× Mix-n-Stain Reaction Buffer to 1× with the antibody solution. Mix by pipetting up and down a few times.
3. Transfer the entire solution from step 2 to the vial containing the CF dye (no need to weigh the dye). Vortex the vial for a few seconds.
4. Incubate the vial in the dark for 30 minutes. The solution is now ready to use. The concentration of the CF dye-labeled antibody conjugate is approximately the amount of the starting antibody divided by the total volume (i.e., ~100% labeling yield).
5. Dilute the labeled antibody solution with the provided Storage Buffer. Simply transfer the entire labeled antibody solution into the Storage Buffer and store at 2–8 °C. Recalculate the concentration of the antibody solution.
Notes: Storage buffer contains 2 mM sodium azide.

Alternatively, the labeled antibody solution can be aliquoted and stored at –20 °C. Without repeated freeze-thaws, the labeled antibody solution remains active for at least 6 months.

D. Modified Mix-n-Stain Labeling

Note: Before beginning, use Table 1 to select the appropriate labeling protocol for the antibody.

1. Use the antibody solution at a concentration of 0.5–1 mg/mL total protein (IgG plus stabilizer protein) for optimal labeling, using 1× PBS to dilute the solution if necessary. Make sure the amount of total protein (μg) in the labeling reaction falls within the range of the kit. To label an amount of IgG that falls below the lower limit of the kit, add BSA to bring to the total protein concentration (IgG + BSA) within the range of the kit and proceed with labeling based on total protein concentration.

2. Warm up the Mix-n-Stain Reaction Buffer vial and the Mix-n-Stain Storage Buffer vial to room temperature before use. Centrifuge the vials briefly to collect the solutions at the bottom of the vials.
3. Dilute the 10× Mix-n-Stain Reaction Buffer to 1× with the antibody solution. Mix by pipetting up and down a few times.
4. Transfer the entire solution from step 3 to the vial containing the CF dye (no need to weigh the dye). Vortex the vial for a few seconds.
5. Incubate the vial in the dark for 30 minutes.
Optional: One can transfer the entire labeling reaction to the tube of antibody storage buffer provided. However, this may result in a highly dilute IgG solution, which may not be practical for subsequent use. To transfer the antibody to storage buffer without additional dilution, follow steps 6–8.
Note: Storage buffer contains 2 mM sodium azide.
6. Transfer the labeling reaction to the membrane of the ultrafiltration vial provided (or saved from Procedure B). Centrifuge the vial at 14,000 × g until all of the liquid has filtered into the receiving vial as described in Procedure B.
7. Resuspend the labeled antibody in antibody storage buffer at the desired final concentration of IgG. Carefully pipette the storage buffer up and down over the upper surface of the membrane to recover and resuspend the antibody.
Note: Storage buffer contains 2 mM sodium azide.
8. Transfer the recovered antibody solution to a fresh microcentrifuge tube. The antibody is now ready to use for staining. The labeled antibody is stable for at least 6 months when stored at 2–8 °C, protected from light. Alternatively, the antibody can be stored in single use aliquots at –20 °C for longer term storage. Without repeated freeze-thaws, the labeled antibody solution remains active for at least 6 months.

Mix-n-Stain and CF are trademarks of Biotium.
BD Horizon is a trademark of BD Biosciences.
eFluor is a registered trademark of eBioscience.
Pacific Blue is a trademark of Molecular Probes, Inc.

AKN,MAM 03/13-1