

Product Information

CF™440 maleimide

Catalog Number: 92124

Unit Size: 1.0 μmole

Color and Form: Yellow solid.

Storage and Handling

Store CF™440 maleimide desiccated at ≤ -20°C. When stored as directed, CF™440 maleimide should be stable for at least 6 months from the time of receipt.

Technical Summary

Abs/Em Maxima: 440/515 nm (See Figure 1)

Extinction coefficient: 40,000

Molecular weight: ~839

A₂₈₀/A_{max} or CF: 0.044 (correction factor for estimating degree of protein labeling)

Direct replacement for: Alexa Fluor® 430

Solubility

Soluble in H₂O, DMF, or DMSO. For making stock solution, we recommend dissolving the dye in anhydrous DMSO (Biotium cat# 90082) at 10 mM.

Product Application

CF™440 maleimide is a green fluorescent dye with a maleimide reactive group, which can be used to label a target molecule with a thiol group. CF™440 (abs/em=440 nm/515 nm) is a superior alternative to Alexa Fluor 430 owing to its exceptional photostability.

CF™440 maleimide reacts with thiol groups to form thioether-coupled products. The reaction can take place at pH 7 in the presence of amines. Under the neutral pH condition, the maleimide group does not react with histidine or arginine.

Protocol for Labeling IgG antibodies

The protocol below is for labeling proteins. Protocols for labeling other thiol-containing molecules are similar except for the purification procedures which may need to be modified.

1. Materials Required but not Provided

- 10-100 mM phosphate (e.g., PBS), Tris or HEPES buffer with pH 7.0-7.5
- Sephadex G-25
- Anhydrous dimethylsulfoxide (DMSO, #90082) for preparing stock solution
- (optional) Tris-(2-carboxyethyl)phosphine (TCEP, #91049) for reducing disulfide binds in proteins to produce free thiol groups.
- BSA

2. Labeling Procedure

2.1 Prepare protein solution for labeling.

a) Dissolve the protein at 50-100 μM (7.5 mg/mL-15 mg/mL for IgG) in any of the mentioned buffers (See Materials section) at room temperature.

b) As an optional step, if you wish to free up more thiol groups from the disulfide bonds in the protein, you may add ~10-fold molar excess of TCEP at this stage. Incubate the reaction solution for ~30 min. The reduction reaction and the subsequent labeling reaction are best to be carried out in the presence of an inert gas (N₂ or Ar) to prevent re-formation of disulfide bonds.

2.2 Prepare dye stock solution

Let a vial of the CF™ maleimide (1 μmole) warm up to room temperature. Add 0.1 mL anhydrous DMSO (Biotium Cat# 90082) to the vial to form a 10 mM dye stock solution. Vortex the vial briefly to fully dissolve the dye, followed by brief centrifugation to concentrate the dye at the bottom of the vial. If the labeling reaction is to be carried out with a much smaller amount of protein, the dye stock solution may need to be more dilute for accurate pipetting.

Note: 1) Any left-over stock solution may be stored at -20°C for later use. If anhydrous DMSO is used for making the solution, the dye should be stable for at least one month.

2) Dye stock solution may also be prepared in de-ionized water. However, because the dye will hydrolyze slowly, the stock solution in water should only be prepared immediately before the conjugation reaction and cannot be stored for later use.

2.3 Carry out the labeling reaction

a) While stirring or vortexing the protein solution, add the dye stock to result in a dye/protein molar ratio of 10-20.

b) Continue to stir or rock the reaction solution at room temperature for 2 hour or at 4 °C overnight.

Tip: while the labeling reaction is underway, proceed to the next step (Step 2.4a) to prepare a Sephadex G-25 column.

2.4 Separate the labeled protein from the free dye

a) Prepare a Sephadex G-25 column (10 mm x 300 mm) equilibrated in PBS buffer (pH~7.4).

b) Load the reaction solution from Step 2.3b onto the column and elute the column with PBS buffer. The first band excluded from the column corresponds to the antibody conjugate.

Note: For small scale labeling reaction, you may use a ultrafiltration device, such as an ultrafiltration device (MWCO~10k) (Biotium Cat#22004), to remove the free dye from the conjugate in order to avoid overly dilute product.

3. Determination of Degree of Labeling

3.1 Determine the protein concentration

The concentration of the antibody conjugate can be calculated from the formula:

$$[\text{conjugate}](\text{mg/mL}) = \left(\frac{A_{280} - A_{\text{max}} \times C_f}{1.4} \right) \times (\text{dilution factor})$$

where [conjugate] is the concentration of the antibody conjugate collected from the column; "dilution factor" is the fold of dilution used for spectral measurement; A_{280} and A_{max} are the absorbance readings of the conjugate at 280 nm and the absorption maximum (~440 nm for CF™440), respectively; C_f is the absorbance correction factor (0.044 for CF™440); and the value 1.4 is the extinction coefficient of IgG in mL/mg.

Note: the protein solution eluted from the column may be too concentrated for accurate absorbance measurement and thus must be diluted to approximately ~0.1 mg/mL. The fold of dilution ("dilution factor") necessary can be estimated from the amount of starting antibody and the total volume of the protein solution collected from the column.

3.2 Calculate the degree of labeling (DOL)

The DOL is calculated according to the formula:

$$\text{DOL} = \frac{A_{\text{max}} \times \text{Mwt} \times (\text{dilution factor})}{\epsilon \times [\text{conjugate}]}$$

where A_{max} , "dilution factor" and [conjugate] are as defined in Step 3.1, Mwt is the molecular weight of IgG (~150,000), and ϵ is the molar extinction coefficient of CF™440 (i.e., ~40,000).

4. Storage and Handling

For long-term storage, we recommend that BSA and sodium azide be added to the conjugate solution to final concentrations of 5-10 mg/mL and 0.01-0.03%, respectively, to prevent denaturation and microbial growth. The conjugate solution should be stored at 4 °C and protected from light.

Spectral Property

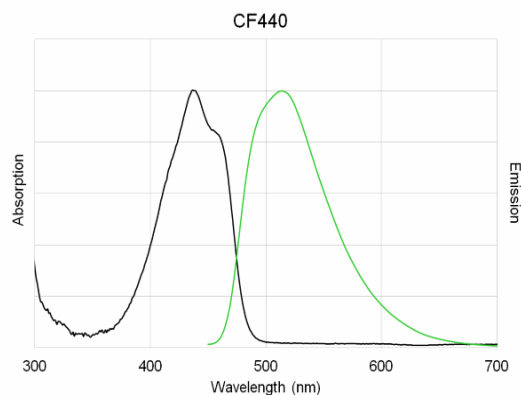


Figure 1. Absorption and emission spectra of CF™440 conjugated to goat anti-mouse IgG in PBS.

Other Related Products

A full selection of secondary antibodies, antibody labeling kits, and other bioconjugates including phalloidin, annexin V and α -bungarotoxin are available for many of our CF™ dyes. Please visit the Biotium website at www.biotium.com for details.

*CF™ dye technology is covered by pending US and international patents.
**Alexa Fluor is a registered trademark of Invitrogen.