Labeling Antibodies with mFluorTM Dye Succinimidyl Esters

Introduction

mFluorTM dyes are a series of excellent fluorescent labeling dyes that span the full visible spectrum. All the mFluorTM dyes have excellent water solubility and large Stokes Shifts. Their hydrophilic property minimizes the use of organic solvents. The mFluorTM dyes have been predominantly used to label antibodies for multicolor flow cytometry applications.

Succinimidyl (NHS) esters are proven to be the best reagents for amine modifications because the amide bonds that are formed are essentially identical to, and as stable as the natural peptide bonds. These reagents are generally stable and show good reactivity and selectivity with aliphatic amines. There are few factors that need be considered when succinimidyl esters compounds are used for conjugation reaction: 1). *Solvents:* For the most part, reactive dyes should be dissolved in anhydrous dimethylformamide (DMF) or dimethylsulfoxide (DMSO). 2). *Reaction pH:* The labeling reactions of amines with succinimidyl esters are strongly pH dependent. Amine-reactive reagents react with non-protonated aliphatic amine groups, including the terminal amines of proteins and the \varepsilon-amino groups of lysines. Thus amine acylation reactions are usually carried out above pH 7.5. Protein modifications by succinimidyl esters can typically be done at pH 8.5-9.5. 3). *Reaction Buffers:* Buffers that contain free amines such as Tris and glycine and thiol compounds must be avoided when using an amine-reactive reagent. Ammonium salts (such as ammonium sulfate and ammonium acetate) that are widely used for protein precipitation must also be removed (such as via dialysis) before performing dye conjugations. 4). *Reaction Temperature:* Most conjugations are done at room temperature. However, either elevated or reduced temperature may be required for a particular labeling reaction.



Storage and Handling

Upon receipt, mFluorTM dyes should be stored at <-15 °C, and kept from light and moisture. The reconstituted DMSO stock solution of mFluorTM dyes can be stored at <-15 °C for less than two weeks. The protein conjugate should be stored at > 0.5 mg/mL in the presence of a carrier protein (e.g., 0.1% bovine serum albumin). The conjugate solution could be stored at 4 °C for two months without significant change when stored in the presence of 2 mM sodium azide and kept from light. For longer storage, the protein conjugates could be lyophilized or divided into single-used aliquots and stored at \leq -60 °C, and protected from light.

Sample Labeling Protocol

Note: This labeling protocol was developed for the conjugate of Goat anti-mouse IgG with $mFluor^{TM}$ 450 SE. You might need further optimization for your particular proteins.

1. Prepare protein stock solution (Solution A):

Mix 100 μ L of a reaction buffer (e.g., 1 M sodium carbonate solution or 1 M phosphate buffer with pH ~9.0) with 900 μ L of the target protein solution (e.g. antibody, protein concentration >2 mg/ml if possible) to give 1 mL protein labeling stock solution.

Note 1: The pH of the protein solution (Solution A) should be 8.5 \pm 0.5. If the pH of the protein solution is lower than 8.0, adjust the pH to the range of 8.0-9.0 using 1 M sodium bicarbonate solution or 1 M pH 9.0 phosphate buffer.

Note 2: The protein should be dissolved in 1X phosphate buffered saline (PBS), pH 7.2-7.4. If the protein is dissolved in Tris or glycine buffer, it must be dialyzed against 1X PBS, pH 7.2-7.4, to remove free amines or ammonium salts (such as ammonium sulfate and ammonium acetate) that are widely used for protein precipitation.

Note 3: Impure antibodies or antibodies stabilized with bovine serum albumin (BSA) or gelatin will not be labeled well. The presence of sodium azide or thimerosal might also interfere with the conjugation reaction. Sodium azide or thimerosal can be removed by dialysis or spin column for optimal labeling results.

Note 4: The conjugation efficiency is significantly reduced if the protein concentration is less than 2 mg/mL. For optimal labeling efficiency the final protein concentration range of 2-10 mg/mL is recommended.

2. Prepare dye stock solution (Solution B):

Add anhydrous DMSO into the vial of mFluor™ dye SE to make a 10-20 mM stock solution. Mix well by pipetting or vortex.

Note: Prepare the dye stock solution (Solution B) before starting the conjugation. Use promptly. Extended storage of the dye stock solution may reduce the dye activity. Solution B can be stored in freezer for two weeks when kept from light and moisture. Avoid freeze-thaw cycles.

3. Determine the optimal dye/protein ratio (optional):

Note: Each protein requires distinct dye/protein ratio, which also depends on the properties of dyes. Over labeling of a protein could detrimentally affects its binding affinity while the protein conjugates of low dye/protein ratio gives reduced sensitivity. We recommend you experimentally determine the best dye/protein ratio by repeating Steps 4 and 5 using a serial different amount of labeling dye solutions. In general 4-6 dyes/protein are recommended for most of dye-protein conjugates.

3.1 Use 10:1 molar ratio of Solution B (dye)/Solution A (protein) as the starting point: Add 5 µl of the dye stock solution (Solution B, assuming the dye stock solution is 10 mM) into the vial of the protein solution (95 µl of Solution A) with effective shaking. The concentration of the protein is ~0.05 mM assuming the protein concentration is 10 mg/mL and the molecular weight of the protein is ~200KD.

Note: The concentration of the DMSO in the protein solution should be <10%.

- 3.2 Run conjugation reaction (see Step 4 below).
- 3.3 Repeat #3.2 with the molar ratios of Solution B/Solution A at 5:1; 15:1 and 20:1 respectively.
- 3.4 Purify the desired conjugates using premade spin columns.
- 3.5 Calculate the dye/protein ratio (DOS) for the above 4 conjugates (see next page).
- 3.6 Run your functional tests of the above 4 conjugates to determine the best dye/protein ratio to scale up your labeling reaction.

4. Run conjugation reaction:

4.1 Add the appropriate amount of dye stock solution (Solution B) into the vial of the protein solution (Solution A) with effective shaking.

Note: The best molar ratio of Solution B/Solution is determined from Step 3.6. If Step 3 is skipped, we recommend using 10:1 molar ratio of Solution B (dye)/Solution A (protein).

4.2 Continue to rotate or shake the reaction mixture at room temperature for 30-60 minutes.

5. Purify the conjugation

The following protocol is an example of dye-protein conjugate purification by using a Sephadex G-25 column.

- 5.1 Prepare Sephadex G-25 column according to the manufacture instruction.
- 5.2 Load the reaction mixture (directly from Step 4) to the top of the Sephadex G-25 column.
- 5.3 Add PBS (pH 7.2-7.4) as soon as the sample runs just below the top resin surface.
- 5.4 Add more PBS (pH 7.2-7.4) to the desired sample to complete the column purification. Combine the fractions that contain the desired dye-protein conjugate.

Note 1: For immediate use, the dye-protein conjugate need be diluted with staining buffer, and aliquoted for multiple uses.

Note 2: For longer term storage, dye-protein conjugate solution need be concentrated or freeze dried (see below).

Characterize the Desired Dye-Protein Conjugate

The Degree of Substitution (DOS) is the most important factor for characterizing dye-labeled protein. Proteins of lower DOS usually have weaker fluorescence intensity, but proteins of higher DOS (e.g. DOS > 6) tend to have reduced fluorescence too. The optimal DOS for most antibodies is recommended between 2 and 10 depending on the properties of dye and protein. For effective labeling, the degree of substitution should be controlled to have 4-10 moles of mFluorTM Violet 450 SE to one mole of antibody. The following steps are used to determine the DOS of mFluorTM Violet 450 SE labeled proteins.

Last Updated: February 2019

1. Measure absorption:

To measure the absorption spectrum of a dye-protein conjugate, it is recommended to keep the sample concentration in the range of $1-10 \mu M$ depending on the extinction coefficient of the dye.

2. Read OD (absorbance) at 280 nm and dye maximum absorption ($\lambda_{max} = 403$ nm for mFluorTM Violet 450 dyes):

For most spectrophotometers, the sample (from the column fractions) need be diluted with de-ionized water so that the OD values are in the range of 0.1 to 0.9. The O.D. (absorbance) at 280 nm is the maximum absorption of protein while 403 nm is the maximum absorption of mFluorTM Violet 450 SE. To obtain accurate DOS, make sure that the conjugate is free of the non-conjugated dye.

3. Calculate DOS using the following equations:

- 3.1 Calculate protein concentration [Protein] = $\frac{\text{A280- (OD @ Dye Maximum Absorrtion X CF@280nm)}}{\text{Protein Extinction Coefficient}} \times \text{dilution factor}$
- 3.2 Calculate dye concentration [Dye] = $\frac{\text{OD @ Maximum Absortion}}{\text{Dye Extinction Coefficient}} \times \text{dilution factor}$
- 3.3 Calculate the degree of labeling DOS = [Dye]/[Protein] = $[^{D}OD_{403} \times ^{P} \epsilon_{280}]$ /[35000×(A₂₈₀-0.238A₄₀₃)]

[Dye] is the dye concentration, and can be readily calculated from the Bee-Lambert Law: $A=\epsilon_{dye}CL$. [Protein] is the protein concentration. This value can be either estimated by the weight (added to the reaction) if the conjugation efficiency is high enough (preferably > 70%) or more accurately calculated by the Beer-Lambert Law: $A=\epsilon_{protein}CL$. For example, IgG has the ϵ value to be 203,000 cm⁻¹M⁻¹. $^{P}\epsilon_{280}$ = protein molar extinction coefficient at 280 nm (e. g. the molar extinction coefficient of IgG is 203,000 cm⁻¹M⁻¹). CF (dye absorption correction factor at 280 nm) = $OD_{280}/OD_{403} = 0.238$ for mFluorTM Violet dye SE. 35,000 cm⁻¹M⁻¹ is the molar extinction coefficient of mFluorTM Violet 450 SE.

References

- 1. Hermanson GT (1996). *Biocojugate Techniques*, Academic Press, New York.
- 2. Haugland RP (1995). Coupling of monoclonal antibodies with fluorophores. *Methods Mol Biol* **45**, 205-21.
- 3. Brinkley M (1992). A brief survey of methods for preparing protein conjugates with dyes, haptens, and cross-linking reagents. *Bioconjug Chem* **3**, 2-13.

Last Updated: February 2019

Appendix 1. mFluor™ Fluorescence-Labeling Dye Selection Guide Chart

mFluor TM Dye	Ex(nm)	Em (nm)	Features and Benefits	Ordering Information	
mFluor™ Blue 570	553 nm	564 nm	 Alternative to RPE Higher labeling yield Photostable 	#1160 (SE, NH ₂ -reactive)	
mFluor™ Green 620	522 nm	617 nm	Optimized for Green Laser Excitation Strong fluorescence Photostable	#1165 (SE, NH ₂ -reactive)	
mFluor™ Red 700	650 nm	694 nm	Alternative to Alexa Fluor®700-APC Tandem Higher labeling yield Photostable	#1190 (SE, NH ₂ -reactive)	
mFluor™ Red 780	651nm	768 nm	Alternative to Alexa Fluor® Alexa Fluor®700-APC Tandem Higher labeling yield Photostable	#1191 (SE, NH ₂ -reactive)	
mFluor™ UV 375	351nm	388 nm	Alternative to BD Biosciences' BUV395 Readily conjugated to antibodies	#1135 (SE, NH ₂ -reactive)	
mFluor™ UV 460	357nm	456 nm	Readily conjugated to antibodies	#1136 (SE, NH ₂ -reactive)	
mFluor™ Violet 450	405 nm	445 nm	Alternative to Pacific Blue TM Brighter on some antibodies More water-soluble	#1150 (SE, NH ₂ -reactive) #1100 (labeling kit)	
mFluor™ Violet 500	433 nm	497 nm	Alternative to Pacific Green™ and BD Horizon™ V500 ■ Brighter on some antibodies ■ More water-soluble	#1149 (SE, NH ₂ -reactive)	
mFluor™ Violet 510	411 nm	504 nm	Alternative to AmCyan and Pacific Green™ ■ pH-insensitive fluorescence ■ Photostable	#1151 (SE, NH ₂ -reactive) #1110 (labeling kit)	
mFluor TM Violet 540	401 nm	532 nm	Alternative to Pacific Orange™ ■ pH-insensitive fluorescence ■ Photostable	#1152 (SE, NH ₂ -reactive) #1114 (labeling kit)	
mFluor™ Yellow 630	554 nm	625 nm	Optimized for Yellow Laser Excitation Strong fluorescence Photostable	#1170 (SE, NH ₂ -reactive)	

Appendix 2. Spectral Properties of mFluorTM Fluorescent Labeling Dyes

Dye	Abs _{max} (nm)	Em _{max} (nm)	ε 1	Φ2	CF at 260 nm ³	CF at 280 nm ⁴
mFluor Red 780	631	768	70000	0.034^{6}	0.101	0.116
mFluor Red 700	657	694	250000	0.029^6	0.135	0.127
mFluor Violet 450	406	445	25000	0.92	0.338	0.078
mFluor Violet 510	409	504	30000	0.86	0.464	0.366
mFluor Blue 570	552	564	120000	0.08^{6}	0.228	0.179
mFluor Green 620	521	617	50000	0.06*	0.895	0.569
mFluor Violet 500	426	497	35000	0.81	0.769	0.365
mFluor Violet 540	400	532	15000	0.64	1.392	0.529
mFluor Yellow 630	546	625	110000	0.016	0.283	0.413
mFluor UV 375	354	388	35000	0.94	0.099	0.138
mFluor UV 460	362	456	15000	0.86	0.35	0.134

Note 1.1. Extinction coefficient at their maximum absorption wavelength. The units of extinction coefficient are $cm^{-1}M^{-1}$. 2. FQY = fluorescence quantum yield in aqueous buffer (pH 7.2). 3. CF at 260 nm is the correction factor used for eliminating the dye contribution to the absorbance at 260 nm. 4. CF at 280 nm is the correction factor used for eliminating the dye contribution to the absorbance at 280 nm (for peptides and protein labeling). 5. Fluorescence intensity is significantly increased upon coupling to proteins. 6. Fluorescence quantum yield may increase upon conjugation to antibodies.

Note 2. Pacific BlueTM, Pacific GreenTM and Pacific OrangeTM are the trademarks of Molecular Probes. mFluorTM is the trademark of AAT Bioquest.