

*This is intended as a guide only for protein/antibody labeling; optimal conjugation protocols may vary depending on the target being labeled.*

## Conjugation Protocol for Amine-Reactive Dyes

Succinimidyl esters (NHS esters) are amine-reactive reagents that can be conjugated to (non-protonated) aliphatic amine groups. The primary reactive species for protein amine-conjugation are the  $\epsilon$ -amino groups of lysine residues. To avoid protonating these groups it is important to perform the reaction at a slightly basic pH. In addition, buffers containing primary amines should be avoided, since they will compete for conjugation with the NHS ester.

*Please note that NHS esters can be moisture sensitive, so handle accordingly. Where possible, handle and store fluorescent dyes in the dark.*

**Note:** A typical molar ratio for labeling Janelia Fluor®, SE dye:protein is 15:1, however this should be optimized for each specific protein. We recommend trying three different molar ratios of dye:protein to develop a protein-specific protocol for future use.

### Reagents

- Prepare a 10 mM stock solution of NHS ester dye in anhydrous DMSO or DMF. Briefly vortex.
- Prepare protein/antibody to be conjugated at approximately 3.0 mg/mL (minimum recommended protein concentration is 2.0 mg/mL†) in either sodium borate (50 mM, pH 8.5) or carbonate buffer (100 mM, pH 8-8.5). N.B. for carbonate buffer, we recommend adding 75 mg/mL Sodium Bicarbonate solution as indicated below to adjust the pH of the starting protein/antibody PBS solution to approximately 8.25.

*Protein/Antibody Volume ( $\mu$ L)  $\times$  0.112 = Volume Sodium Bicarbonate required ( $\mu$ L).*

† N.B. This is critical for achieving the optimal degrees of labeling (DOLs) for Janelia Fluor® dyes.

## Conjugation

- Add dye solution to protein solution at the appropriate molar ratio (see note above) while stirring/gently vortexing.
- Incubate at room temperature for 60 minutes in the dark†.
- Quench the reaction by adding Tris-HCl or Glycine (pH 7.4, 50-100 mM final concentration), incubate with stirring for 10-15 minutes at room temperature (this step is optional).

† N.B. Increasing the incubation time to 18 hours in the dark can increase the DOL with Janelia Fluor® dyes.

## Purification

- Remove excess dye with a Zeba™ Spin desalting column (Thermo), with appropriate MWCO, or a PD MiniTrap™ G-25 (GE Healthcare), following the manufacturer's instructions.

## Degree of labeling calculation

- Dilute the protein-dye conjugate to approximately 0.1 mg/mL and measure the absorbance at 280 nm (protein  $A_{280}$ ) and at the maximum absorbance wavelength ( $A_{max}$ ) for the fluorescent dye used (please refer to the individual product descriptions for the max.  $\lambda_{abs}$  values).
- Calculate the corrected  $A_{280}$  ( $A_{280C}$ ) using the following equation:

$$A_{280C} = A_{280} - (A_{max} \times CF)$$

*Please refer to the individual product descriptions for the correction factor (CF) values for the fluorescent dye used.*

- Calculate the final protein concentration (**[protein]** in mg/mL) using the corrected  $A_{280C}$  value calculated above, the extinction coefficient ( $\epsilon$ ) for your protein and the Beer-Lambert Law equation:

$$A_{280} = \epsilon \times [protein] \times l$$

- Calculate the final F:P ratio (Degree of Labeling), where  $\epsilon_{dye}$  is the extinction coefficient for the fluorescent dye used, *please refer to the individual product descriptions for this value.*

$$F:P = \frac{A_{max} \times MW_{protein}}{[protein] \times \epsilon_{dye}}$$

*Zeba is a trademark of Thermo Scientific and MiniTrap is a trademark of GE Healthcare*