

Andy Fluor™ NHS Ester

Protein Labeling Quick Reference

Quick Protocol

1. Dissolve ~10 mg of the protein in 1 mL of 0.1 M sodium bicarbonate buffer pH ~8.3.
2. Dissolve 1 µmol of the Andy Fluor™ NHS Ester in 100 µL of DMSO or DMF.
3. While stirring the protein solution, slowly add 50-100 µL of the Andy Fluor™ NHS Ester solution from step 2.
4. Incubate the reaction for 1 hour at room temperature with continuous stirring.
5. Equilibrate a 10 × 300 mm gel filtration column (Sephadex® G-25, BioGel® P-30, or equivalent) with PBS.
6. Separate the conjugate on the gel filtration column.
7. Store the conjugates under the same conditions used for the parent protein.

For detail protocol visit www.abpbio.com/product/andy-fluor-dyes/