

Technical Note

Biotinylate antibody from stock which contains carrier protein (Recommended procedure for use with FortéBio streptavidin biosensors)

Materials required:

- Antibody to be biotinylated (>200µg antibody in the presence of carrier protein; lyophilized or in solution)
- NAb Protein A or G Purification Kit (Pierce cat # 45200 or 45201)
- EZ-Link NHS-LC-LC-Biotin (Pierce cat # 21343)
- Dimethylformamide (DMF)
- 1x PBS
- Slide-A-Lyzer, 10000 MWCO (Pierce cat# 66383)
- De-salt Dextran Columns (Pierce cat #43230)- optional

Time required:

Antibody isolation: 1.5 hour hands-on; 1 day + overnight total time

Biotinylation: 1 hour hands-on; 1 day + overnight total time

Protocol:

1. Antibody isolation

- 1.1. **NOTE:** For mIgG2a, mIgG2b and IgG3 use Protein A Spin Purification Kit. For mIgG1 use the Protein G Spin Purification Kit.
- 1.2. Mix resin gently to make slurry; use cut pipet tip to transfer 200 µL of the slurry to spin cup column placed in a collection tube. Tip should be cut to provide a wider opening for pipeting the resin.
- 1.3. Add 300 µL of the appropriate Binding buffer from Purification Kit to resin. Mix gently. Centrifuge 1 min at ~5000 g. Discard solution from collection tube.
- 1.4. Add 400 µL of Binding buffer to resin. Mix gently. Centrifuge 1 min at ~5000 g. Discard solution from collection tube.
- 1.5. Repeat wash 1 more time.
- 1.6. If the antibody is lyophilized, add 300 µL of the binding buffer.
- 1.7. If the antibody is in solution, add an equal volume of binding buffer (eg 250 µL of antibody plus 250 µL binding buffer)
- 1.8. Add antibody solution to resin. Cap cup and incubate 30 minute with gentle shaking.
- 1.9. Uncap cup and centrifuge in collection tube 1 minute.

- 1.10. Move cup to a new collection tube. Add 400 μ L binding buffer and mix briefly. Centrifuge 1 minute.
- 1.11. Repeat previous wash step 2 more times.
- 1.12. Transfer to a new collection tube. Add 400 μ L of Elution Buffer. Cap cup and mix gently for 5 minutes.
- 1.13. Uncap and centrifuge 1 minute. Transfer cup to a new collection tube.
- 1.14. Repeat steps 2 more times for a total of 3 elutions. Elutions should be neutralized with 40 μ L of 1M sodium phosphate (pH 8).
- 1.15. A denaturing, non-reducing gel should be run of the washes and elutions to determine efficiency of isolation and location of the antibody (**Figure 1**).

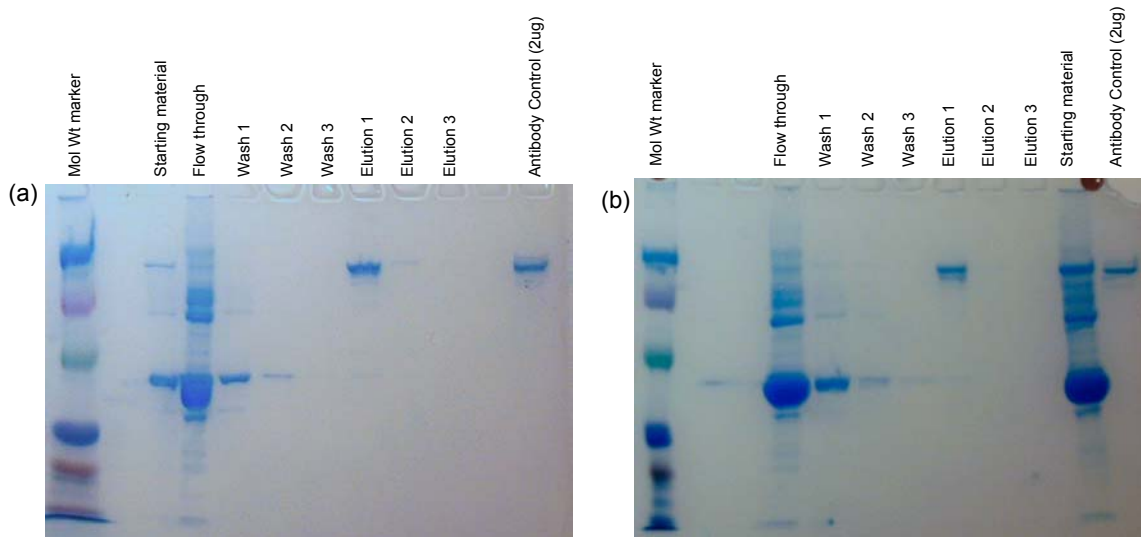


Figure 1: Denaturing, non-reducing gels of fractions from antibody isolation. **(a)** using Protein G spin purification with mIgG1. **(b)** using Protein A spin purification with mIgG2a

2. Biotinylation

- 2.1. Using a MWCO 10000 dialyzer, dialyze the elution fractions containing the antibody against PBS at 4C (elution 1 in both examples shown in **Figure 1**).
- 2.2. **NOTE:** Since the elution buffer contains primary amines, dialysis must be extensive. Six changes at 1:1000 (> 3 hours between each change) are recommended. Use of a desalting column is not recommended for this step as the procedure is not efficient enough to remove all the reactive amines from the buffer.
- 2.3. Recover the dialyzed sample from the dialyzer and transfer to a new eppendorf tube.
- 2.4. Prepare a 10mM Biotin reagent solution: add 2.0 mg NHS-LC-LC-Biotin reagent in 350 μ L of DMF. Mix to dissolve.
- 2.5. Calculate the volume of 10 mM biotin reagent needed based on the mass of antibody started with prior to antibody isolation. For most antibodies, a

5:1 molar coupling ratio (moles NHS-LCLC-biotin: moles antibody at start of procedure).

$$\frac{\text{mg protein}}{\text{MW (mg/mmol)}} \times \frac{5 \text{ mmol biotin}}{1 \text{ mmol protein}} \times \frac{1000\text{mL}}{10 \text{ mmol biotin}} \times \frac{1000 \mu\text{L}}{1 \text{ mL}} = \mu\text{L of 10mM solution of biotin reagent}$$

- 2.6. To each sample, add the appropriate **volume (μL)** of NHS-LC-LC-Biotin reagent as calculated. Mix immediately
- 2.7. Incubate 30 minutes at Room Temperature.
- 2.8. Stop the reaction by removing the excess biotin reagent by either dialysis (recommended) or desalting column. Dialysis should involve 4 changes of PBS at 1:1000.
- 2.9. **NOTE:** Take care to remove all free biotin in order to most efficiently bind the biotinylated protein to the streptavidin sensor surface.
- 2.10. Biotinylated antibodies in PBS can be stored at 4C. See **Figure 2** for example immobilization of biotinylated antibody onto streptavidin SBC biosensor.

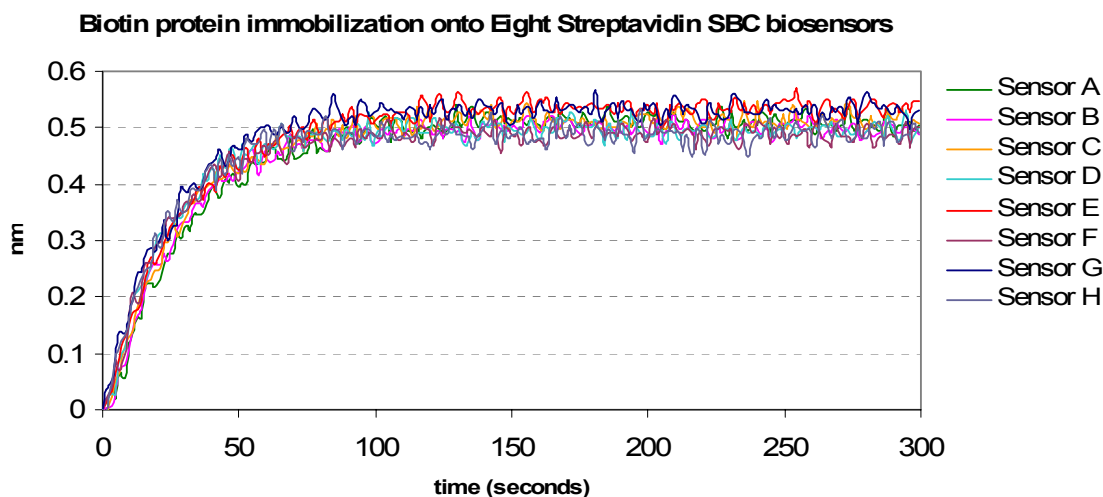


Figure 2: Example data for the parallel immobilization of a biotinylated protein onto 8 streptavidin SBC biosensors.