

## Buccutite™ Rapid APC-Cy5.5 Tandem Antibody Labeling Kit \*Microscale Optimized for Labeling 100 ug Antibody Per Reaction\*

Catalog number: 1320

Unit size: 2 Labeledings

| Component                                       | Storage                                       | Amount                |
|---|---|-----------------------|
| Component A: Buccutite™ FOL-Activated APC-Cy5.5 | Refrigerate (2-4 °C), Minimize light exposure | 2 vials (lyophilized) |
| Component B: Buccutite™ MTA                     | Refrigerate (2-4 °C), Minimize light exposure | 2 vials (lyophilized) |
| Component C: Reaction Buffer                    | Refrigerate (2-4 °C), Minimize light exposure | 1 vial (20 µL)        |
| Component D: Spin Column                        | Room temperature                              | 2 columns             |

### OVERVIEW

APC-Cy5.5 is a popular color used in flow cytometry. Its primary absorption peak is at 651 nm with emission peak at ~700 nm. AAT Bioquest offers this Buccutite™ rapid labeling kit to facilitate the APC-Cy5.5 tandem conjugations to antibodies and other proteins such as streptavidin and other secondary reagents. Buccutite™ APC-Cy5.5 Tandem Conjugation Kit provides a robust and convenient method to conjugate antibodies with APC. The kit includes a preactivated APC-Cy5.5 tandem and reaction buffer. The conjugated antibody can be used in WB, ELISA and IHC applications. This kit is sufficient for 2 labeling reactions, each up to 100 ug of antibody. The best ratio for any new antibody reagent must be determined by experimentation. Our kit provides preactivated APC-Cy5.5 tandem to facilitate the APC-Cy5.5 tandem conjugations to antibodies and other proteins such as streptavidin and other secondary reagents. Our preactivated APC-Cy5.5 tandem is ready to conjugate, giving much higher yield than the conventionally tedious SMCC-based conjugation chemistry. In addition, our preactivated APC-Cy5.5 tandem is conjugated to a protein via its amino group that is abundant in proteins while SMCC chemistry targets the thiol group that has to be regenerated by the reduction of antibodies.

### AT A GLANCE

#### Protocol summary

1. Add 5 µL Reaction Buffer (Component C) into antibody (100 µL)
2. Add the antibody solution into Buccutite™ MTA vial (Component B)
3. Incubate at room temperature
4. Remove free Buccutite™ MTA by spin column
5. Mix with 50 µL Buccutite™ FOL-Activated APC-Cy5.5 (Component A)
6. Incubate at room temperature

**Important** Upon receipt, store the kit at 4 °C. When stored properly, the kit should be stable for six months. Alternatively, Component B can be stored at -20°C. Do not freeze Buccutite™ FOL-Activated APC-Cy5.5 (Component A), Reaction Buffer (Component C) and Spin Column (Component D). Warm all the components and centrifuge the vials briefly before opening, and immediately prepare the required solutions before starting your conjugation. The following SOP is an example for labeling goat anti-mouse IgG antibody.

### PREPARATION OF WORKING SOLUTION

#### Antibody working solution:

For labeling 100 µg antibody (assuming the target antibody concentration is 1 mg/mL), mix 5 µL (5% of the total reaction volume) of Reaction Buffer (Component C) with 100 µL of the target antibody solution.

**Note** If you have a different concentration, adjust the antibody volume accordingly to make ~100 µg antibody available for your labeling reaction.

**Note** The antibody should be dissolved in 1X phosphate buffered saline (PBS), pH 7.2-7.4; If the antibody is dissolved in glycine buffer, it must be dialyzed against 1X PBS, pH 7.2-7.4, or use Amicon Ultra-0.5, Ultracel-10 Membrane, 10 kDa ( Cat. # UFC501008 from Millipore) to remove free amines or ammonium salts (such as ammonium sulfate and ammonium acetate) that are widely used for antibody precipitation.

**Note** Impure antibodies or antibodies stabilized with bovine serum albumin (BSA) or gelatin will not be labeled well.

**Note** The antibody –Buccutite™ MTA reaction efficiency is significantly reduced if the antibody concentration is less than 1 mg/mL. For optimal labeling efficiency the final antibody concentration range of 1-10 mg/mL is recommended.

### SAMPLE EXPERIMENTAL PROTOCOL

#### Run Antibody-Buccutite™ MTA reaction

1. Add the antibody solution directly into the vial of Buccutite™ MTA (Component B), and mix them well by repeatedly pipetting for a few times or vortex the vial for a few seconds.
2. Keep the antibody- Buccutite™ MTA reaction mixture at room temperature for 30 - 60 minutes.

**Note** The antibody-Buccutite™ MTA reaction mixture can be rotated or shaken for longer time if desired.

#### Prepare spin column for antibody-Buccutite™ MTA purification

1. Invert the provided spin column (Component D) several times to re-suspend the settled gel and remove any bubbles.
2. Snap off the tip and place the column in a washing tube (2 mL, not provided). Remove the cap to allow the excess packing buffer to drain by gravity to the top of the gel bed. If column does not begin to flow, push cap back into column and remove it again to start the flow. Discard the drained buffer, and then place the column back into the Washing Tube. However, centrifuge immediately if the column is placed into a 12 x 75 mm test tube (not provided).
3. Centrifuge for 2 minutes in a swinging bucket centrifuge at 1,000 x g (see Centrifugation Notes section) to remove the packing buffer. Discard the buffer.
4. Apply 1-2 mL 1X PBS (pH 7.2-7.4) to the column. After each application of PBS, let the buffer drain out by gravity, or centrifuge the column for 2 minutes to remove the buffer. Discard the buffer from the collection tube. Repeat this process for 3-4 times.
5. Centrifuge for 2 minutes in a swinging bucket centrifuge at 1,000 x g (see Centrifugation Notes section) to remove the packing buffer. Discard the buffer.

## Purify the antibody-Buccutite™ MTA solution

1. Place the column (from Step Prepare spin column for antibody-Buccutite™ MTA purification) in a clean Collecting Tube (1.5 mL, not provided). Carefully load the sample (~105 µL, from Step Antibody-Buccutite™ MTA reaction) directly to the center of the column.
2. After loading the sample, add 5 µL of 1X PBS (pH 7.2-7.4) to make the total volume of 110 µL. Centrifuge the column for 5-6 minutes at 1,000 x g, and collect the solution that contains the desired antibody-Buccutite™ MTA solution.

## Make antibody-APC-Cy5.5 conjugation

1. Make Buccutite™ FOL-Activated APC-Cy5.5 solution by adding 50 µL ddH<sub>2</sub>O into the vial of Buccutite™ FOL-Activated APC-Cy5.5 (Component A), mix well by repeatedly pipetting for a few times or vortex the vial for a few seconds.
2. Mix whole vial of Buccutite™ FOL-Activated APC-Cy5.5 solution into the purified antibody-Buccutite™ MTA solution (from Step Purify the antibody-Buccutite™ MTA solution), mix well and rotating the mixture for 1 hour at room temperature.
3. The antibody-APC-Cy5.5 conjugate is now ready to use.

**Note** For immediate use, the antibody-APC-Cy5.5 conjugate need be diluted with the buffer of your choice.

**Note** For longer term storage, antibody-APC-Cy5.5 conjugate solution need be concentrated or freeze dried.

## Storage of Antibody-APC-Cy5.5 Conjugate

The antibody conjugate should be stored at > 0.5 mg/mL in the presence of a carrier antibody (e.g., 0.1% bovine serum albumin). The Antibody-APC-Cy5.5 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 antibody-APC-Cy5.5 conjugates could be lyophilized and stored at ≤ -20 °C.

## Centrifugation Notes

Spin column (Component D) can fit into 2 mL microcentrifuge tubes or 12 x 75 mm test tubes for sample collection during centrifugation. Use the 2 mL microtube with the columns for the initial column equilibration step.

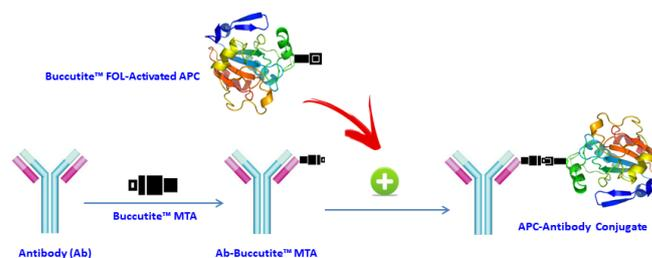
Swinging bucket centrifuges capable of generating a minimum force of 1,000 x g are suitable for Bio-Spin column use. The gravitational force created at a particular revolution speed is a function of the radius of the microcentrifuge rotor. Consult the swinging bucket centrifuge instruction manual for the information about conversion from revolutions per minute (RPM) to centrifugal or g-force. Alternatively, use the following equation to calculate the speed in RPM required to reach the gravitational force of 1,000 x g.

$RCF (x g) = (1.12 \times 10^{-5}) \times (RPM)^2 \times r$  (RCF is the relative centrifugal force, r is the radius in centimeters measured from the center of the rotor to the middle of the Bio-Spin column, and RPM is the speed of the rotor).

**Table 1.** Available fluorophores at AAT Bioquest Buccutite™ Rapid Antibody Labelling Kits

| Cat# | Labels          | Ex (nm) | Em (nm) |
|------|-----------------|---------|---------|
| 1310 | PE              | 565     | 575     |
| 1322 | PE-Cy5          | 565     | 674     |
| 1316 | PE-Cy5.5        | 565     | 700     |
| 1317 | PE-Cy7          | 565     | 780     |
| 1318 | PE-Texas Red    | 565     | 600     |
| 1311 | APC             | 651     | 662     |
| 1319 | APC-iFluor™ 700 | 651     | 713     |
| 1320 | APC-Cy5.5       | 651     | 700     |
| 1321 | APC-Cy7         | 651     | 780     |
| 1325 | PerCP           | 482     | 677     |

## EXAMPLE DATA ANALYSIS AND FIGURES



**Figure 1.**

AAT Bioquest offers this Buccutite™ rapid labeling kit to facilitate the APC-Cy5.5 tandem conjugations to antibodies and other proteins such as streptavidin and other secondary reagents. Our preactivated APC was premodified with our Buccutite™ FOL. Your antibody (or other proteins) is modified with our Buccutite™ MTA to give MTA-modified protein (such as antibody). The MTA-modified protein readily reacts with FOL-modified APC to give the desired APC-antibody conjugate in much higher yield than the SMCC chemistry. In addition, our preactivated APC reacts with MTA-modified biopolymers at much lower concentrations than the SMCC chemistry.

## DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.