

## Buccutite™ Rapid PE-iFluor® 710 Tandem Antibody Labeling Kit \*Microscale Optimized for Labeling 25 ug Antibody Per Reaction\*

Catalog number: 1357  
Unit size: 2 Labelings

| Component  | Storage  | Amount (Cat No. 1357) |
|--|--|-----------------------|
| Component A: Buccutite™ FOL-Activated PE-iFluor® 710 | Refrigerated (2-8 °C), Minimize light exposure | 2 Vials (Lyophilized) |
| Component B: Buccutite™ MTA                          | Freeze (< -15 °C), Minimize light exposure     | 1 Vial (Lyophilized)  |
| Component C: Reaction Buffer                         | Freeze (< -15 °C), Minimize light exposure     | 1 Vial (20 uL)        |

### OVERVIEW

The Buccutite™ Rapid PE-iFluor® 710 Tandem Antibody Labeling Kit offers a highly efficient and reproducible method for small-scale conjugation of antibodies with PE-iFluor® 710. Utilizing the advanced Buccutite™ crosslinking platform, this kit significantly simplifies the labeling workflow compared to conventional strategies such as SMCC-mediated crosslinking. The streamlined two-step protocol allows for the rapid conjugation of antibodies or proteins to PE-iFluor® 710 in under two hours, with minimal hands-on time.

Each kit includes all reagents necessary for two labeling reactions, with each reaction optimized to conjugate 25 µg of purified antibody or protein using Buccutite™ FOL-Activated PE-iFluor® 710. For optimal performance, the removal of stabilizing proteins (e.g., BSA) and avoidance of amine-containing buffers such as Tris are critical, as these components can interfere with the conjugation chemistry.

PE-iFluor® 710 is a tandem fluorophore with excitation and emission maxima at ~565 nm and ~747 nm, respectively, offering a substantial Stokes shift and high fluorescence intensity. These properties make it particularly well-suited for applications such as multicolor flow cytometry, spectral flow cytometry, and other fluorescence-based immunoassays requiring high sensitivity. However, due to its limited photostability, it is not recommended for applications involving prolonged light exposure. This kit enables direct labeling of primary antibodies, eliminating the need for secondary antibody-based detection systems. The resulting conjugates reduce experimental complexity and enhance assay sensitivity and reproducibility.

### AT A GLANCE

#### Protocol Summary

1. Add 1.25 µL Reaction Buffer (Component C) into antibody (25 µL)
2. Add 2.5 µL Buccutite™ MTA working solution
3. Incubate at room temperature for 30 - 60 minutes
4. Mix with 50 µL Buccutite™ FOL-Activated PE-iFluor® 710 working solution
5. Incubate at room temperature for 60 minutes

**Important:** Upon receipt, store the kit at 4 °C. When stored properly, the kit should be stable for six months. Alternatively Components A and B can be stored at -20 °C. Do not freeze Reaction Buffer (Component C). 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

1. To label 25 µg of antibody (assuming the target antibody concentration is 1 mg/mL), mix 1.25 µL (5% of the total reaction

volume) of Reaction Buffer (Component C) with 25 µL of the target antibody solution.

**Note:** If your antibody has a different concentration, adjust the volume to ensure approximately 25 µg of antibody is available for the 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:** Ensure the antibody is pure and not stabilized with bovine serum albumin (BSA) or gelatin, as these stabilizers significantly impair labeling efficiency.

**Note:** For optimal labeling efficiency, the antibody concentration should be in the range of 1–10 mg/mL. If the concentration is below 1 mg/mL, the Buccutite™ MTA reaction efficiency is significantly reduced.

#### Buccutite™ MTA Working Solution

1. Add 10 µL of DMSO (Not provided) to the vial of Buccutite™ MTA (Component B).

#### Buccutite™ FOL-Activated PE-iFluor® 710 Working Solution

1. Add 50 µL of ddH<sub>2</sub>O to the vial of Buccutite™ FOL-Activated PE-iFluor® 710 (Component A).

### SAMPLE EXPERIMENTAL PROTOCOL

#### Run Antibody-Buccutite™ MTA Reaction

1. Add 2.5 µL of Buccutite™ MTA working solution into antibody working solution, 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.

#### Make Antibody-PE-iFluor® 710 Conjugation

1. Add 50 µL of Buccutite™ FOL-Activated PE-iFluor® 710 working solution with Antibody-Buccutite™ MTA solution, mix well by

repeatedly pipetting for a few times or vortex the vial for a few seconds.

2. Incubate for 1 to 2 hours.
3. The antibody-PE-iFluor® 710 conjugate is now ready to use.

**Note:** For immediate use, the antibody-PE-iFluor® 710 conjugate must be diluted with the buffer of your choice.

**Note:** For longer term storage, antibody-PE-iFluor® 710 conjugate solution must be concentrated or freeze dried.

#### 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](mailto:info@aatbio.com) if you have any questions.

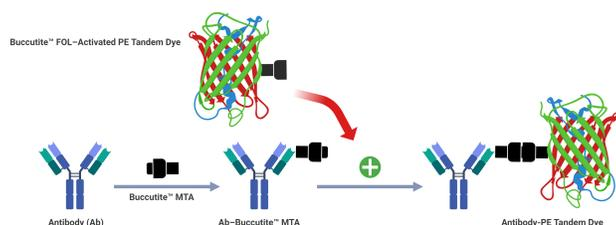
#### Storage of Antibody-PE-iFluor® 710 Conjugate

The antibody conjugate should be stored in the presence of a carrier protein (e.g., 0.1% bovine serum albumin) and 0.02-0.05% sodium azide. The Ab-PE-iFluor® 710 conjugate solution could be stored at 4 °C for two months without significant change and kept from light.

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

| Cat# | Labels          | Ex (nm) | Em (nm) |
|------|-----------------|---------|---------|
| 1325 | PerCP           | 482     | 677     |
| 1310 | PE              | 565     | 575     |
| 1318 | PE-Texas Red    | 565     | 600     |
| 1356 | PE-iFluor® 594  | 565     | 606     |
| 1322 | PE-Cy5          | 565     | 674     |
| 1316 | PE-Cy5.5        | 565     | 700     |
| 1358 | PE-iFluor® 710  | 565     | 710     |
| 1317 | PE-Cy7          | 565     | 780     |
| 1311 | APC             | 651     | 662     |
| 1320 | APC-Cy5.5       | 651     | 700     |
| 1319 | APC-iFluor® 700 | 651     | 713     |
| 1321 | APC-Cy7         | 651     | 780     |

#### EXAMPLE DATA ANALYSIS AND FIGURES



**Figure 1.** AAT Bioquest offers the Buccutite™ rapid labeling kit to streamline PE tandem dye conjugation for antibodies and other proteins, including streptavidin and secondary reagents. This kit utilizes preactivated PE modified with Buccutite™ FOL, while your antibody or protein is modified with Buccutite™ MTA to produce MTA-modified proteins. The MTA-modified proteins react efficiently with FOL-modified PE, yielding the desired PE-antibody conjugate with significantly higher efficiency compared to traditional SMCC chemistry. Additionally, the reaction requires much lower biopolymer concentrations, enhancing efficiency and reducing material usage compared to SMCC-based methods.