

Buccutite™ ALP (Alkaline Phosphatase) Antibody Conjugation Kit *Optimized for Labeling 25 ug Protein*

 Catalog number: 5512
 Unit size: 2 Labelings

Component	Storage	Amount
Component A: Buccutite™ FOL-Activated ALP	Refrigerated (2-8 °C), Minimize light exposure	2 vials
Component B: Buccutite™ MTA	Refrigerated (2-8 °C), Minimize light exposure	2 vials (Lyophilized)
Component C: Reaction Buffer	Refrigerated (2-8 °C), Minimize light exposure	1 vial (20 µL)

OVERVIEW

Protein-protein conjugations are commonly performed with a bifunctional linker (such as SMCC) having different reactivity at each end for linking two different proteins. One end of the crosslinker reacts (via NHS ester) with amines (-NH₂) of lysine and N-terminus, and the other end reacts (via maleimide) with the thiol groups (-SH) of cysteine. However, SMCC-modified protein is extremely unstable and often self-reactive since proteins contain multiple amine and thiol groups that cause significant amount of crosslinkings. In addition, it is quite difficult and tedious to quantify the number of maleimide groups in a protein. Buccutite™ ALP (Alkaline Phosphatase) Antibody Conjugation Kit is designed to rapidly prepare ALP conjugates with pre-activated ALP. The kit is optimized for labeling 25 ug protein. The Buccutite™ FOL-activated ALP readily reacts with Buccutite™ MTA-modified antibody under extremely mild neutral conditions without any catalyst required. Compared with commonly used SMCC and other similar technologies, Buccutite™ bioconjugation system is much more robust and easier to use. It enables faster and quantitative conjugation of biomolecules with higher efficiency and yields.

AT A GLANCE

Protocol summary

1. Add 1.25 µL of Reaction Buffer (Component C) into antibody (25 µL)
2. Add 2.5 µL of reconstituted Buccutite™ MTA (Component B) to the antibody working solution
3. Incubate the mixture at room temperature for 30 minutes
4. Mix the antibody- Buccutite™ MTA reaction mixture with Buccutite™ FOL-Activated ALP (Component A)
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, Component B can be stored at -20 °C. Warm up all the components at room temperature and centrifuge the vials briefly before opening. Immediately prepare the required solutions before starting your conjugation. The following SOP is an example for labeling 25 µg goat anti-mouse IgG antibody.

PREPARATION OF WORKING SOLUTION

Antibody working solution

For labeling 25 µg 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 you have a different concentration, adjust the antibody volume accordingly to make ~25 µg antibody /25 µL 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 10 kDa Filter (Cat. # 60502 from AAT Bioquest) to remove free amines or ammonium salts (such as ammonium sulfate and ammonium acetate).

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 10 µL of DMSO (not provided in the kit) into the vial of Buccutite™ MTA to reconstitute Buccutite™ MTA.
2. Add 2.5 µL of Buccutite™ MTA (Component B) to Antibody working solution, and mix them well by repeatedly pipetting for a few times or vortex the vial for a few seconds.
3. Keep the antibody- Buccutite™ MTA reaction mixture at room temperature for 30 minutes.

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

Make antibody-ALP conjugation

1. Add antibody-Buccutite™ MTA reaction mixture to the vial of Buccutite™ FOL-Activated ALP (Component A), mix well and incubate the mixture at room temperature for 1 hour.

Note The antibody-ALP reaction mixture can be incubated for longer time if desired.

2. The antibody-ALP conjugate is now ready to use.

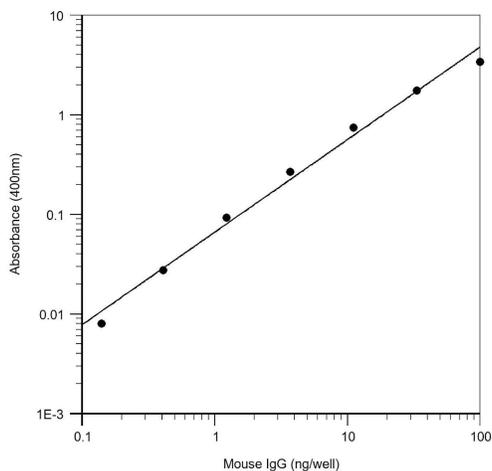
Note For immediate use, the antibody-ALP conjugate needs to be diluted with the buffer of your choice.

3. The concentration of the conjugate can be calculated as follows:
Antibody Concentration (µg/µL) = 25 µg (total amount of antibody)/(25 µL + 1.25 µL+2.5 µL) = 0.86 µg/µL

Storage of Antibody-ALP Conjugate

The antibody conjugate-ALP conjugate should be stored in the presence of a carrier protein (e.g., 0.1% bovine serum albumin) at 4 °C and kept from light for two months.

EXAMPLE DATA ANALYSIS AND FIGURES



Mouse IgG dose curve was detected using the ALP- goat-anti-mouse IgG conjugate prepared with Buccutite™ ALP (Alkaline Phosphatase) Antibody Conjugation Kit. 3-fold serial diluted mouse IgG was coated on a 96-well plate, and 100uL GAM IgG-ALP conjugate (1ug/ml) was added to each well followed the standard ELISA method. pNPP substrate solution was used to detect the immobilized mouse IgG with 30 min incubation and read at 400 nm.

Image generated with Quest™ Graph, ©2021 AAT Bioquest

Figure 1. Mouse IgG dose curve was detected using the ALP- goat-anti-mouse IgG conjugate prepared with Buccutite™ ALP (Alkaline Phosphatase) Antibody Conjugation Kit. 3-fold serial diluted mouse IgG was coated on a 96-well plate, and 100uL GAM IgG-ALP conjugate (1ug/ml) was added to each well followed the standard ELISA method. pNPP substrate solution was used to detect the immobilized mouse IgG with 30 min incubation and read at 400 nm.

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.