

Buccutite™ Rapid Protein Crosslinking Kit

Optimized for Crosslinking 100 ug Protein Per Reaction

Ordering Information

Cat#: 1315 (2 conjugations)

Storage Conditions

Multi-Storage Conditions

Introduction:

Protein-protein conjugations are commonly performed with a bifunctional linker SMCC. One end of the SMCC reacts (via NHS ester) with amines (-NH₂) found in the amino acid lysine and N-terminus, and the other end reacts (via maleimide) with the thiol groups (-SH) found in the amino acid cysteine. However, SMCC-modified protein is extremely unstable and often self-reactive since proteins often contain both amine and thiol groups that cause significant amount of homo-crosslinking. In addition it is quite difficult and tedious to quantify the number of maleimide groups on a protein.

A convenient and effective crosslinking method has been developed to link two biomolecules with a high conjugation yield. This method uses one pair of crosslinkers: Buccutite™ MTA and Buccutite™ FOL; MTA is added to one protein, while FOL is added to another protein. Protein-protein cross-linking reaction is initiated by mixing Protein-1- Buccutite™ MTA and Protein-2- Buccutite™ FOL. This crosslinking reaction occurs under extremely mild neutral conditions without any catalyst required, and it is robust and efficient.

Kit Components

Components	Amount	Storage
Component A: Buccutite™ MTA	2 Vial (lyophilized)	-20 °C
Component B: Buccutite™ FOL	2 Vial (lyophilized)	-20 °C
Component C: Reaction Buffer	1 Vial (50 µL)	4 °C (Do not freeze)
Component D: Spin Column	4 Columns	4 °C (Do not freeze)

Key Parameters to Achieve Best Performance:

1. Protein molecular weight: >25,000 Daltons
2. Protein concentration: >=1mg/ml
3. Protein sample volume: 60~120 µL

Operating Protocol (Labeling 100 µg Protein-1)

Upon receipt, store component A and B at -20°C. When stored properly, the kit should be stable for six months. Do not freeze 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 conjugate 100µg protein-1 with protein 2.

1. Prepare Protein Solution:

For labeling 100 µg protein-1 and protein-2 (assuming the concentrations are 1 mg/mL for both proteins), mix 5 µL (5% of the total reaction volume) of Reaction Buffer (Component C) with 100 µL of the each protein solution. The protein should be dissolved in 1X phosphate buffered saline (PBS), pH 7.2-7.4; If it 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 protein precipitation. Stabilizers like bovine serum albumin (BSA) or gelatin will affect the labeling reaction.

2. Run Protein 1-Buccutite™ MTA Reaction:

- 2.1 Add 105 µL Protein-1 solution directly into the vial of Buccutite™ MTA (Component A), and mix them well by repeatedly pipetting for a few times or vortex the vial for a few seconds.

- 2.2 Keep the Protein-1- Buccutite™ MTA reaction mixture at room temperature for 30 - 60 minutes. The Antibody-Buccutite™ MTA reaction mixture can be rotated or shaken for longer time if desired.
- 2.3 Purify Protein-1- Buccutite™ MTA through desalting column. Please refer to “Prepare spin column for Antibody purification” part for detailed operations.
- 2.4 Calculate the concentration of the Protein-1- Buccutite™ MTA with 75% yield after desalting. (For example: if starting with 100µg protein, after desalting column purification, the recovery protein amount is ~75µg.)

3. Run Protein 2-Buccutite™ FOL reaction:

- 3.1 Add the 105µL Protein-2 solution directly into the vial of Buccutite™ FOL (Component B), and mix them well by repeatedly pipetting for a few times or vortex the vial for a few seconds.
- 3.2 Keep the Protein-2- Buccutite™ FOL reaction mixture at room temperature for 30 - 60 minutes. The Antibody-Buccutite™ FOL reaction mixture can be rotated or shaken for longer time if desired.
- 3.3 Purify Protein-2- Buccutite™ FOL through desalting column. Please refer to “Prepare spin column for Antibody purification” part for detailed operations.
- 3.4 Calculate the concentration of the Protein-2- Buccutite™ FOL with 75% yield after desalting.

4. Cross-linking Protein-1 and Protein-2 Reaction:

- 4.1 Cross-linking reaction is initiated by mixing Protein-1- Buccutite™ MTA and Protein-2- Buccutite™ FOL at the desired molar ratio. Usually, Buccutite™ FOL-modified Protein-2 is mixed with Buccutite™ MTA-modified Protein-1 in 1.5-2.0 molar ratio to drive the crosslinking reaction to completion. The mixing ratio can be reversed depending on your downstream applications, and the cost of your proteins.
- 4.2 Rotate the mixture for 1~2 hour at room temperature or stay at 4°C overnight. The reaction mixture is now ready to use, or to be stored at 4 °C. Desalting is optional.

5. Optional: Conjugate Analysis and Purification:

- 5.1 A small amount of reaction mixture (for example: 2~4 µg) could be analyzed using 4-12% Bis-Tis Protein Gel in a SDS running buffer system to check the conjugation result.
- 5.2 The conjugation reaction mixture contains the desired conjugate along with small amount of unlinked Protein-2 (used in excess). If required, the reaction mixture can be purified by size exclusion chromatography (SEC), and the desired conjugate fractions are pooled and combined.

Prepare spin column for Antibody 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.
6. Place the column (from Step 3.5) in a **clean** Collecting Tube (1.5 mL, not provided). Carefully load the sample (~105 µL, from Step 2.2) directly to the center of the column.
7. After loading the sample, add 10 µL of 1X PBS (pH 7.2-7.4), centrifuge the column for 5-6 minutes at 1,000 x g, and collect the solution that contains the desired protein-1-Buccutite™ MTA solution or Protein-2- Buccutite™ FOL solution.

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*).

References

1. Duncan, R.J.S., *et al.* (1983). A new reagent which may be used to introduce sulfhydryl groups into proteins, and its use in the preparation of conjugates for immunoassay. *Anal Biochem* **132**:68-73.
2. Yoshitake, S., *et al.* (1979). Conjugation of glucose oxidase from *Aspergillus niger* and rabbit antibodies using *N*-hydroxysuccinimide ester of *N*-(4-carboxycyclohexyl-methyl)maleimide. *Eur J Biochem* **101**:395-9.
3. Hashida, S., *et al.* (1984). More useful maleimide compounds for the conjugation of Fab' to horseradish peroxidase through thiol groups in the hinge. *J Appl Biochem* **6**:56-63.
4. Imagawa, M., *et al.* (1982). Characteristics and evaluation of antibody- horseradish peroxidase conjugates prepared by using a maleimide compound, glutaraldehyde, and periodate. *J Appl Biochem* **4**:41-57.

Warning: This kit shall be only sold to our authorized distributors and end users. Neither resale nor transfer to a third party is allowed without written permission from AAT Bioquest. Chemical analysis of the kit components is strictly prohibited. Please call us at 408-733-1055 or e-mail us at info@aatbio.com if you have any questions.