Lightning-Link® Rapid Conjugation System

Applicable to: Product series 310 to 369

Release 1 © EXPEDEON, 07/03/2018

INTRODUCTION

Lightning-Link® Rapid conjugation technology works by targeting amine groups (e.g. lysines) and is widely used to label antibodies.

The Lightning-Link® Rapid conjugation kit allows conjugations to be set up in less than 30 seconds, simply by adding a solution of the antibody to a lyophilized mixture containing a proprietary activated ligand (Figure 1).

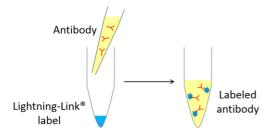


Figure 1. Lightning-Link® antibody conjugation

By circumventing the desalting or dialysis steps that commonly interrupt traditional antibody conjugation procedures, Lightning-Link® technology can be used to label both small (e.g. $10~\mu g$) and large quantities of primary antibodies with ease. Batch-to-batch variation upon scale up is minimal as the process is so simple, and recoveries are always 100%.

Directly labeled primary antibodies are advantageous as they eliminate the need for secondary reagents in immunoassay procedures, thus removing a tedious extra cycle of incubation and wash steps in applications, such as ELISA and Western blotting.

Lightning-Link® technology can also be used to label proteins, peptides and other biomolecules (See 'the What can I label using Lightning-Link®?' Section of the protocol for further information).

KIT CONTENTS

- 1 or 3 glass vial(s) of Lightning-Link® Rapid mix
- 1 vial of LL Rapid Modifier reagent
- 1 vial of LL Rapid Quencher reagent

SHIPPING CONDITIONS

The kit is shipped at ambient temperature in a tamper-evident polypropylene container. Store the kits at -20° C upon receipt.

Please note that the modifier and quencher FD after initial thawing can be stored at either 4° C or -20° C.

BUFFER CONSIDERATIONS

Please see the below table for recommended buffer conditions and components:

| Buffer Components | |
|-------------------------------------|---------|
| рН | 6.5-8.5 |
| Amine free buffer | ✓ |
| (e.g. MES, MOPS, HEPES, PBS) | |
| Non-buffering salts | ✓ |
| (e.g. sodium chloride) | |
| Chelating agents | ✓ |
| (e.g.EDTA) | |
| Sugars | ✓ |
| Glycerol | <50% |
| Thiomersal / Thimerosal | × |
| Merthiolate | × |
| Sodium Azide ¹ | <0.1% |
| BSA ^{1,2} | <0.1% |
| Gelatin ^{1,2} | <0.1% |
| Tris | <50mM |
| Glycine | × |
| Proclin | x |
| Borate buffer | ✓ |
| Nucleophilic components | × |
| (Primary amines e.g. amino acids or | |
| ethanolamine and thiols e.g. | |
| mercaptoethanol or DTT) | |

¹ Please note that individually the concentrations shown should not affect the reaction. However in combination with additional compounds that are not recommended above a certain concentration, the reaction may be affected.

AMOUNT AND VOLUME OF ANTIBODY

| Pack Size | Recommend amount of antibody | Maximum amount of antibody | Maximum Conjugation volume |
|-------------|------------------------------|----------------------------|----------------------------------|
| Up to 20µg | 10μg | 20μg | 10µl |
| Up to 200µg | 100μg | 200μg | 100µl |
| Up to 2mg | 1mg | 2mg | 1ml |

The recommended conjugation conditions are based on using a 1mg/ml antibody concentration. Antibodies greater than 1mg/ml should be diluted to 1mg/ml using either milli-Q water or PBS.

Superior conjugates are normally generated using the recommended amount of antibody. Using the maximum amount of antibody will still generate quality conjugates.

Antibodies less than 1mg/ml can still be used to generate good conjugates provided the maximum conjugation volume is not exceeded. Adding less than the recommended maximum amount of antibody may result in unbound label post conjugation. This excess

² If intending to use this kit for immunohistochemistry, it is recommended that there be no gelatin or BSA present.

label will be deactivated by the quencher and removed during the first wash step of any application. We would recommend that antibodies below 0.5mg/ml are concentrated prior to use.

Please contact our technical support team for more advice.

INSTRUCTIONS

Setting up your conjugation reaction

- 1. Before you add antibody to the Lightning-Link® Rapid mix, add 1µl of LL-Rapid modifier reagent for each 10µl of antibody to be labeled. Mix gently.
- Remove the screw cap from the vial of Lightning-Link® Rapid
 mix and pipette the antibody sample (with added LL-Rapid
 modifier) directly onto the lyophilized material. Resuspend
 gently by withdrawing and re-dispensing the liquid once or
 twice using a pipette.
- 3. Place the cap back on the vial and leave the vial standing for 15 minutes at room temperature (20-25° C) in the dark. Alternatively, and sometimes more conveniently, conjugations can be set up and left at room temperature overnight, as the longer incubation time has no negative effect on the conjugate.
- **4.** After incubating for 15 minutes (or more), add 1µl of LL-Rapid quencher reagent for every 10µl of antibody used. The conjugate can be used after 5 minutes. The conjugates do not require purification.

STORAGE OF CONJUGATES

Your Rapid conjugate can be stored at 4° C for up to 18 months. For longer storage the conjugate can be stored at -20° C with a cryoprotectant such as 50% glycerol.

The best storage conditions for any particular conjugate must be determined by experimentation.

The conjugate should always be stored in the dark.

WHAT CAN I LABEL USING LIGHTNING-LINK®?

Lightning-Link® technology works by targeting free amine groups. It can be used to label antibodies, peptides, proteins and other molecules with free amine groups.

The protocol provided here is optimized for labeling IgGs, when labeling other molecules please contact our technical support team at: www.expedeon.com/contact/

Our custom conjugation service team will also be happy to generate your conjugate for you. Please contact our sales team for a quote. www.expedeon.com/contact/

WHAT IF MY BUFFER DOESN'T FIT THE REQUIREMENTS?

The AbSelect™ purification kit range allows you to quickly and simply purify your antibody using our conjugation friendly buffers.

The appropriate kit to use depends on your particular sample. Please view our AbSelect ™ range of products on our website to select the best kit.

If your antibody is already purified but its concentration is too low, you can concentrate it by using our AbSelect™ Antibody Concentration and Clean Up Kit (product code 861-0010). This kit can also be used

to remove low molecular weight contaminants such as azide, Tris or alvoine.

If your antibody contains BSA, you can use our AbSelect™ BSA removal kit (product code 820-0100) to purify your antibody in one simple step.

Our AbSelect™ kits are fully compatible with Lightning-Link® and are designed to work with IgG antibodies. The only exception is the Concentration and Clean Up kit (product code 861-0010) which will work with any molecule greater than 10kDa.

TECHNICAL SUPPORT

For technical enquiries get in touch with our technical support team at: www.expedeon.com/contact/

Limited use license

Expedeon's products are offered for research purposes alone, and are not intended for human, therapeutic or diagnostic use. The purchase of this product conveys to the buyer (whether the buyer is a not-for-profit, academic or for-profit entity) the non-transferable right to use the amount of product purchased and the components of the product for in-house research. The buyer shall not sell or otherwise transfer this product, its components, or materials prepared therefrom to any third party. The buyer shall not use this product or its components for commercial purposes. For the avoidance of doubt, 'commercial purposes' means any activity by a party for consideration and includes, without limitation, use of the product or its components (i) in the manufacturing of other materials (ii) to provide a service, information or data, (iii) for therapeutic, diagnostic or prophylactic purposes, or (iv) for repackaging/resale, whether or not such product or its components are resold for use in research. The use of this product by the buyer constitutes agreement with the terms of this limited use label

Many of Expedeon's products and technologies are protected by patents or are the subject of patent filings. For information on purchasing a license for commercial use please contact us at +44 (0) 1223 873364

