



Lightning-Link® PE-Cy5 Tandem Conjugation Kit

Applicable to:

760-0030 3 vials (to label 10µg Ab per vial)

760-0005 1 vial (to label 60µg Ab per vial)

Release 6

760-0010 3 vials (to label 60µg Ab per vial)

760-0015 1 vial (to label 0.6mg Ab per vial)

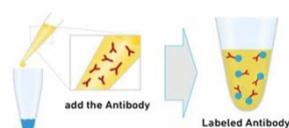
02/08/2016

Introduction

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

The Lightning-Link® PE-Cy5 Tandem conjugation kit allows PE-Cy5 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 PE-Cy5 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 µg) 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 this protocol for further information).

Kit contents

1 or 3 glass vial(s) of Lightning-Link® mix

1 vial of LL-modifier reagent

1 vial of LL-quencher FD 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	
pH	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	✗
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.

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

Amount and volume of antibody

Product code	Recommended maximum amount of antibody per vial	Recommended maximum antibody volume per vial
760-0030	10µg	10µl
760-0005 760-0010	60µg	60µl
760-0015	600µg	600µl

The recommended conjugation conditions are based on using a 1mg/ml antibody concentration and are designed to give a 1:1. Ab:PE-Cy5 conjugation molar ratio.

Antibodies greater than 1mg/ml should be diluted to 1mg/ml using either milli-Q water or PBS.

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 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® mix, add 1µl of LL-modifier reagent for each 10µl of antibody to be labeled. Mix gently.
2. Remove the screw cap from the vial of Lightning-Link® mix and pipette the antibody sample (with added LL-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 3 hours 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 in the dark, as the longer incubation time has no negative effect on the conjugate.
4. After incubating for 3 hours (or more), add 1µl of LL-quencher FD reagent for every 10µl of antibody used. The conjugate can be used after 30 minutes. No separation steps are necessary.

Storage of conjugates

Your PE-Cy5 tandem conjugate can be stored at 4°C for up to 3 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.

Limited use license

Innova Biosciences' 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 Innova Biosciences' 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 our Business Development Office, Babraham Hall, Babraham, Cambridge, UK, CB22 3AT. Tel +44(0)1223 661000; e-mail info@innovabiosciences.com

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, we would recommend when labeling other molecules you visit our dedicated alternative applications page, www.innovabiosciences.com/innova/advanced-users-guide.html

our Lightning-Link FAQ page for antibody related questions <https://www.innovabiosciences.com/faqs/antibody-labeling-faqs.html>

or contact our technical support team. technical.enquiries@innovabiosciences.com

Our custom conjugation service team will also be happy to generate your conjugate for you. Please contact our sales team for a quote. info@innovabiosciences.com

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. We have designed a handy flow chart on the AbSelect™ webpage to help you select the best kit; <https://www.innovabiosciences.com/images/stories/innova/pdfs/abselect%20purification%20flowchart.pdf>

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 glycine.

If your antibody contains AbSelect™ BSA, you can use our 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 (861-0010), which will work with any molecule greater than 10kDa.

I have a question regarding my conjugation

Please contact our technical support team at technical.enquiries@innovabiosciences.com if you require any further information or help about our products.