

ReadiLink Antibody Labeling Kit

Catalog #	Description
1351002	ReadiLink 492/516 Antibody Labeling Kit
1351004	ReadiLink 594/610 Antibody Labeling Kit
1351006	ReadiLink 647/674 Antibody Labeling Kit
1351011	ReadiLink 405/454 Antibody Labeling Kit

For Research Use Only. Not for use in diagnostic procedures.

Introduction

Bio-Rad's ReadiLink Antibody Labeling Kits offer easy fluorescence conjugations for microscale volumes (50–100 µg). Easily label antibodies in two simple steps without the need for a purification step (Figure 1).

Each ReadiLink Dye is coupled to a reactive moiety (a succinimidyl ester). The reactive dye selectively binds to primary amines of proteins to form a stable carboxamide bond, ensuring no dissociation between the fluorophore and antibody. After conjugation and the addition of the quencher buffer, any unbound ReadiLink Dye will bind to the quencher and become nonfluorescent.

The ReadiLink Antibody Labeling Kit provides all the essential components for performing two conjugation reactions (2 x 50 µg protein). Each kit can be used to label monoclonal or polyclonal antibodies, or other proteins (>10 kD).

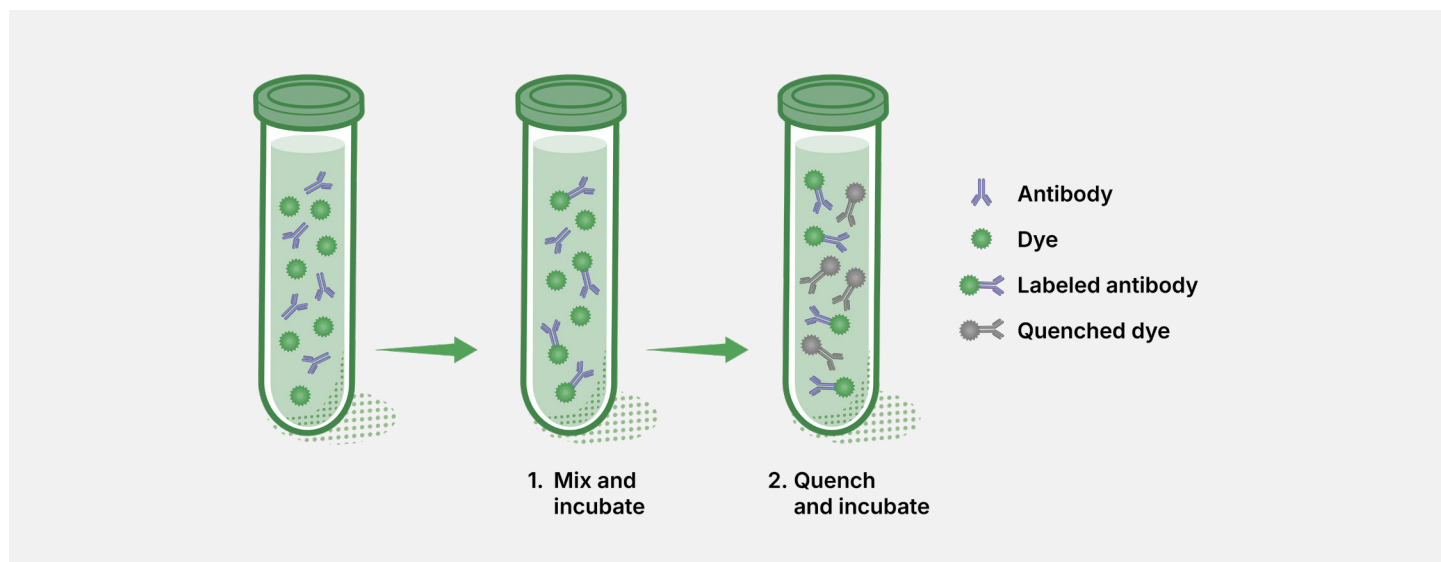


Fig. 1. Efficient labeling of antibodies in two easy steps.



Kit Contents and Storage

Follow the guidelines in Table 1 for storing kit components.

Table 1. Kit components and storage.

Kit Component	Quantity	Storage, °C
ReadiLink Labeling Dye (powder)	2 vials (Note: use one vial to label 50 µg of antibodies)	-20
Reaction buffer	1 vial (20 µL)	-20
Quench buffer	1 vial (20 µL)	-20

Directions for Labeling 50 µg of Antibodies

Important: Thaw all kit components prior to use.

Note: The antibody of interest must be suspended in phosphate buffered saline (PBS), pH 7.2–7.4 or dialyzed against PBS prior to conjugation to remove free amines or ammonium salts in the solution.

Note: The optimal antibody conjugation is 1 mg/mL. A conjugation performed at a different concentration may cause suboptimal labeling.

1. Suspend the antibody of interest in PBS to create a 1 mg/mL concentration.
2. Add 5 µL reaction buffer to 50 µL antibody solution from step 1.
3. Mix well by pipetting up and down a few times.
4. Add the entire volume (55 µL) to one vial of labeling dye and mix by pipetting.
5. Incubate for 60 min at room temperature.
6. Add 5 µL quench buffer to the reaction mixture.
7. Incubate for 10 min at room temperature.
8. Antibodies are now labeled and ready to use.

ReadiLink Antibody Labeling Kits with Excitation and Emission Wavelengths

Use Table 2 to select the appropriate ReadiLink Antibody Labeling Kit for your conjugation.

Table 2. Excitation and emission specifications.

Catalog Number	Labeling Dye Description	Maximum Excitation, nm	Maximum Emission, nm
1351002	ReadiLink 492/516	492	516
1351004	ReadiLink 594/610	594	610
1351006	ReadiLink 647/674	647	674
1351011	ReadiLink 405/454	405	454

Visit [bio-rad-antibodies.com/readilink](https://www.bio-rad-antibodies.com/readilink) for more information.

Bio-Rad is a trademark of Bio-Rad Laboratories, Inc. All trademarks used herein are the property of their respective owner. © 2026 Bio-Rad Laboratories, Inc.



Bio-Rad
Laboratories, Inc.

Life Science
Group

Website [bio-rad.com](https://www.bio-rad.com) USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 00 800 00 24 67 23 Belgium 00 800 00 24 67 23 Brazil 55 11 3065 7550 Canada 1 800 361 1808 China 86 21 6169 8500 Czech Republic 00 800 00 24 67 23 Denmark 00 800 00 24 67 23 Finland 00 800 00 24 67 23 France 00 800 00 24 67 23 Germany 00 800 00 24 67 23 Greece 30 210 7774396 Hong Kong 852 2789 3300 Hungary 00 800 00 24 67 23 India 91 124 4029300 Israel 000 800 00 24 67 23 Italy 00 800 00 24 67 23 Japan 81 3 6361 7000 Korea 82 080 007 7373 Luxembourg 00 800 00 24 67 23 Mexico 52 55 5488 7670 The Netherlands 00 800 00 24 67 23 New Zealand 64 9 415 2280 Norway 00 800 00 24 67 23 Poland 00 800 00 24 67 23 Portugal 00 800 00 24 67 23 Russian Federation 7 495 721 14 04 Singapore 65 6415-3170 South Africa 27 21 531 7504 Spain 00 800 00 24 67 23 Sweden 00 800 00 24 67 23 Switzerland 00 800 00 24 67 23 Taiwan 886 2 2578 7189 Thailand 662 651 8311 United Arab Emirates 971 4 818 7300 United Kingdom 00 800 00 24 67 23