

ReadiLink Antibody Labeling Kits



Bio-Rad's ReadiLink Antibody Labeling Kits offer easy fluorescence conjugations for microscale volumes (50–100 μ g).

Key Benefits of ReadiLink Antibody Labeling Kits

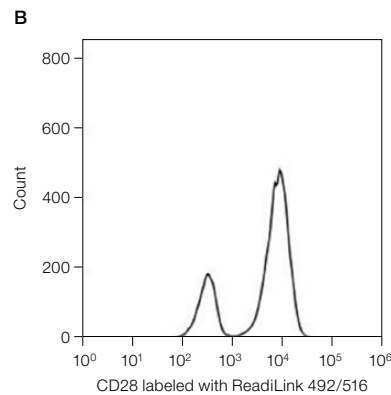
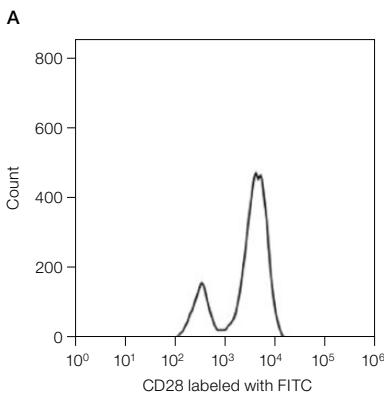
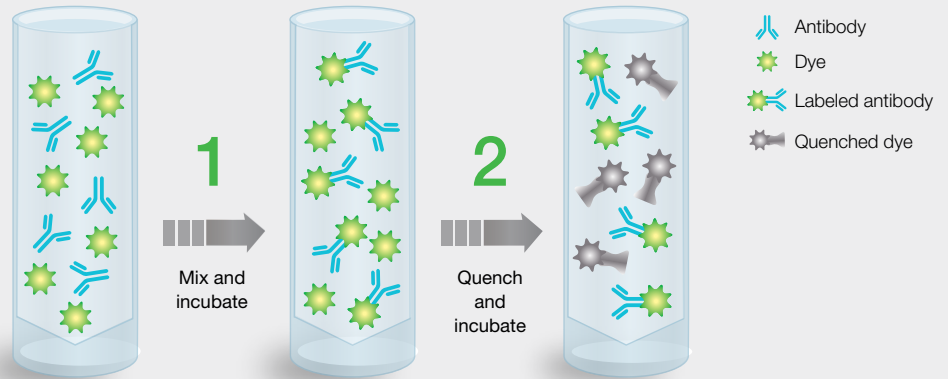
- Save time with no purification column
- Start your experiment sooner with a 2-step antibody conjugation
- Easily conjugate 1 of 13 different fluorophores to an antibody to fit into your multicolor fluorometric experiment

Visit bio-rad.com/web/readilink1 for more information.

BIO-RAD

Efficiently Label Antibodies in Two Easy Steps

Quickly label any antibody of interest in two easy steps with the unique chemistry of ReadILink Dyes. Each ReadILink Dye includes a succinimidyl ester moiety that reacts with primary amines of proteins to form a stable carboxamide bond. After conjugation and the addition of the quencher buffer, any unbound ReadILink Dye will bind to the quencher and become nonfluorescent.



A ReadILink Antibody Labeling Kit shows excellent separation between positive and negative populations. A commercially labeled CD28 FITC (clone CD28.2) (A) compared to the same antibody labeled with ReadILink 492/516 (B). FITC, fluorescein isothiocyanate.

Ordering Information

Catalog Number	ReadILink Antibody Labeling Kit	Maximum Excitation, nm	Maximum Emission, nm	S3™ Cell Sorter Laser, nm	488/561 nm Filter Channel
135-1001	ReadILink 350/440	350	440	–	–
135-1002	ReadILink 492/516	492	516	488	FL1 (525/30)
135-1003	ReadILink 555/570	555	570	561	FL2 (586/25)
135-1004	ReadILink 594/610	594	610	561	FL3 (615/25)
135-1005	ReadILink 633/655	633	655	–	–
135-1006	ReadILink 647/674	647	674	–	–
135-1007	ReadILink 680/701	680	701	–	–
135-1008	ReadILink 700/713	700	713	–	–
135-1009	ReadILink 750/780	750	780	–	–
135-1010	ReadILink 790/811	790	811	–	–
135-1011	ReadILink 405/454	403	454	–	–
135-1012	ReadILink 405/508	414	508	–	–
135-1013	ReadILink 405/537	405	537	–	–

ReadILink is a trademark of AAT Bioquest, Inc.



BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA 800 424 6723 Australia 61 2 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 11 3065 7550
Canada 905 364 3435 China 86 21 6169 8500 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00
France 01 47 95 69 65 Germany 089 31 884 0 Greece 30 210 9532 220 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300
Israel 03 963 6050 Italy 39 02 216091 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666
New Zealand 64 9 415 2280 Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04
Singapore 65 6415 3188 South Africa 27 861 246 723 Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 026 674 55 05
Taiwan 886 2 2578 7189 Thailand 1800 88 22 88 United Kingdom 020 8328 2000