



ReadiLink Antibody Labeling Kits

Bio-Rad's ReadiLink Antibody Labeling Kits offer rapid fluorescent labeling of up to 100 μg of antibody or other proteins for flow cytometry and cell imaging studies. Choose to label at the bench when a commercially available conjugate is not available, an antibody quantity is limited, or secondary antibodies don't work.

With ReadiLink Antibody Labeling Kits you can:

- Save time and expense with no purification column required
- Use as little as 50 μg /reaction, ideal for limited antibodies
- Have choice in multicolor experiments with 10 different fluorophores
- Label monoclonal and polyclonal antibodies and other proteins >10 kD
- Get equivalent or better brightness than a pre-conjugated alternative
- Simplify experiments by eliminating the need for secondary antibodies

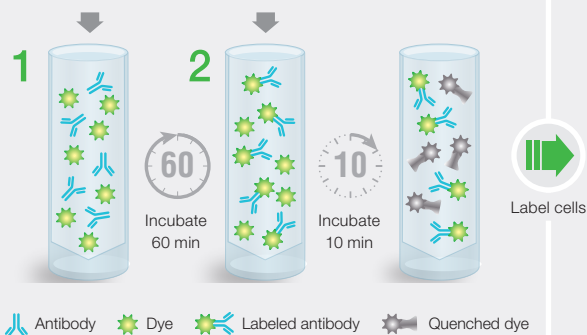
Visit bio-rad.com/ReadiLink1 for more information.



Conjugate Antibodies in Two Easy Steps

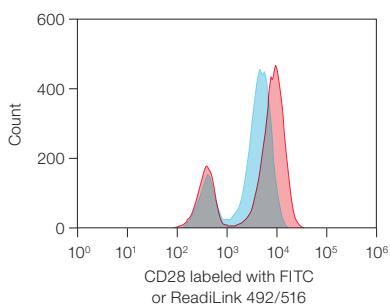
50 µl 1 mg/ml antibody
5 µl reaction buffer
1 vial labeling dye

5 µl quench buffer

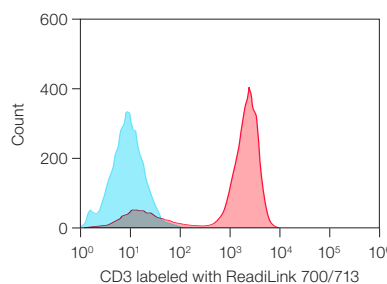


Sort, Analyze, or Image Samples

Use any of these instruments.



ReadiLink 492/516 shows excellent separation and a brighter signal between positive and negative populations on the S3e Cell Sorter. A commercially labeled CD28 FITC (clone CD28.2) (■) was compared to the same antibody labeled with ReadiLink 492/516 (■). FITC, fluorescein isothiocyanate.



ReadiLink 700/713 provides a strong signal to differentiate between positive and negative populations on the ZE5 Cell Analyzer. CD3 labeled with ReadiLink 700/713 at 0.25 µg/ml on peripheral blood mononuclear cells (■) was compared to unstained cells (■).

Ordering Information

Catalog Number	Description	Fluorophore Equivalent	Maximum Excitation, nm	Maximum Emission, nm	Optimal Excitation Laser, nm	S3e Cell Sorter	ZE5 Cell Analyzer	ZOE Fluorescent Cell Imager
1351011	ReadiLink 405/454	Pacific Blue	403	454	405	•	•	
1351002	ReadiLink 492/516	FITC, Alexa Fluor 488, DyLight 488	492	516	488	•	•	•
1351003	ReadiLink 555/570	Cy3, Alexa Fluor 555, DyLight 550	555	570	561	•	•	•
1351004	ReadiLink 594/610	Texas Red, Alexa Fluor 594, DyLight 594	594	610	561	•	•	•
1351005	ReadiLink 633/655	Cy5, Alexa Fluor 633, DyLight 633	633	655	640	•	•	
1351006	ReadiLink 647/674	Cy5, Alexa Fluor 647, DyLight 650	647	674	640	•	•	
1351007	ReadiLink 680/701	Cy5.5, IRDye 700, Alexa Fluor 680, DyLight 680	680	701	640	•	•	
1351008	ReadiLink 700/713	Alexa Fluor 700	700	713	640	•	•	
1351009	ReadiLink 750/780	Alexa Fluor 750, DyLight 755	750	780	640		•	
1351010	ReadiLink 790/811	IRDye 800, Alexa Fluor 790, DyLight 800	790	811				

Each ReadiLink Kit includes 2 vials of labeling dye, reaction buffer, and dye quench for labeling 2 x 50 µg of antibodies. FITC, fluorescein isothiocyanate.

Alexa Fluor, DyLight, Pacific Blue, and Texas Red are trademarks of Thermo Fisher Scientific Inc. Cy is a trademark of GE Healthcare. IRDye is a trademark of LI-COR, Inc. ReadiLink is a trademark of AAT Bioquest, Inc.

BIO-RAD

Bio-Rad Laboratories, Inc.

Life Science Group

Web site bio-rad.com USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 43 01 877 89019 Belgium 32 03 710 53 00 Brazil 55 11 3065 7550 Canada 1 905 364 3435 China 86 21 6169 8500 Czech Republic 36 01 459 6192 Denmark 45 04 452 10 00 Finland 35 08 980 422 00 France 33 01 479 593 00 Germany 49 089 3188 4393 Hong Kong 852 2789 3300 Hungary 36 01 459 6190 India 91 124 4029300 Israel 972 03 963 6050 Italy 39 02 49486600 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 31 0 318 540 666 New Zealand 64 9 415 2280 Norway 47 0 233 841 30 Poland 36 01 459 6191 Portugal 351 21 4727717 Russia 7 495 721 14 04 Singapore 65 6415 3188 South Africa 36 01 459 6193 Spain 34 091 49 06 580 Sweden 46 08 555 127 00 Switzerland 41 0617 17 9555 Taiwan 886 2 2578 7189 Thailand 66 2 651 8311 United Arab Emirates 971 4 8187300 United Kingdom 44 01923 47 1301

