



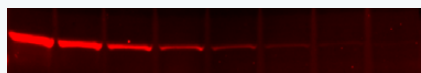
Fluorescent Western Blotting Antibodies

Experience Unmatched Sensitivity and Simpler Multiplexing

StarBright Blue Fluorescent Secondary Antibodies deliver lower limits of detection than traditional methods and even match high sensitivity chemiluminescence. This is possible because they utilize novel fluorescent polymer-based molecules excited by blue light (430–470 nm). StarBright Antibodies are available in two colors to provide multiplexing flexibility and are conjugated to highly cross-absorbed goat anti-mouse and goat anti-rabbit IgG to eliminate cross-reactivity.

StarBright Blue Antibodies

StarBright Blue 700
(blue-excited fluorescence, 20 s)



▲
32 pg

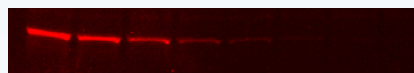
StarBright Blue 520
(blue-excited fluorescence, 20 s)



▲
32 pg

Traditional Detection Methods

IRDye 800CW
(IR-excited fluorescence, Odyssey 9120 Imaging System)



▲
63 pg

Alexa Fluor 488
(blue-excited fluorescence, 20 s)



▲
500 pg

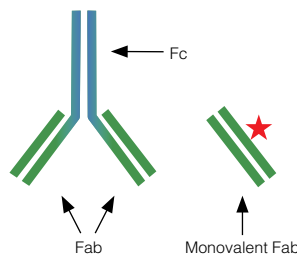
SuperSignal West Femto ECL
(HRP chemiluminescence, 60 s)



▲
32 pg

hFAB Rhodamine Housekeeping Antibodies

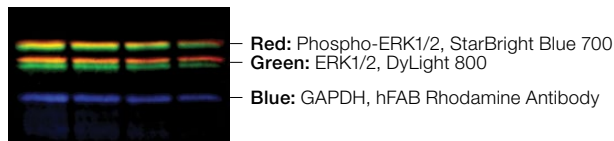
One-step detection of common housekeeping proteins like actin, tubulin, and GAPDH is possible using hFAB Housekeeping Antibodies. These fluorescent antibodies are directly labeled with a rhodamine derivative to make secondary antibodies unnecessary. hFAB Antibodies are generated with human combinatorial antibody library (HuCAL[®]) technology to provide exceptional specificity and eliminate cross-reactivity.



Comparison of typical IgG and an hFAB Rhodamine Antibody.

Detecting Posttranslationally Modified Targets

Fluorescent multiplexing with StarBright Antibodies lets you detect targets with posttranslational modifications. Their flexibility and high sensitivity let you visualize unmodified and modified targets simultaneously — even if only a fraction of the protein is modified.



Example of detecting posttranslationally modified targets.

Easy Multiplexing and Versatile Normalization

StarBright Antibodies are easily adapted into your current protocols. They are the ideal multiplexing partners to traditional secondary antibodies conjugated to DyLight or Alexa fluorophores. Confidently use them to normalize your data with either Bio-Rad's hFAB Rhodamine Housekeeping Antibodies or a total protein approach using stain-free technology.



Example of quadruplex western blotting targets.

Ordering information

Catalog #	Description
StarBright Blue 700 Secondary Antibodies	
12004158	Goat Anti-Mouse IgG StarBright Blue 700, 400 µl
12004159	Goat Anti-Mouse IgG StarBright Blue 700, 80 µl
12004161	Goat Anti-Rabbit IgG StarBright Blue 700, 400 µl
12004162	Goat Anti-Rabbit IgG StarBright Blue 700, 80 µl
StarBright Blue 520 Secondary Antibodies	
12005866	Goat Anti-Mouse IgG StarBright Blue 520, 400 µl
12005867	Goat Anti-Mouse IgG StarBright Blue 520, 80 µl
12005869	Goat Anti-Rabbit IgG StarBright Blue 520, 400 µl
12005870	Goat Anti-Rabbit IgG StarBright Blue 520, 80 µl

Catalog #	Description
hFAB Rhodamine Antibodies	
12004163	Anti-Actin hFAB Rhodamine Antibody, 200 µl
12004164	Anti-Actin hFAB Rhodamine Antibody, 40 µl
12004165	Anti-Tubulin hFAB Rhodamine Antibody, 200 µl
12004166	Anti-Tubulin hFAB Rhodamine Antibody, 40 µl
12004167	Anti-GAPDH hFAB Rhodamine Antibody, 200 µl
12004168	Anti-GAPDH hFAB Rhodamine Antibody, 40 µl

Visit bio-rad.com/FluorescentAbs for more information on StarBright Antibodies

Bio-Rad is a trademark of Bio-Rad Laboratories, Inc. in certain jurisdictions. HuCAL is a trademark of MorphoSys AG. All trademarks used herein are the property of their respective owner.



**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

