



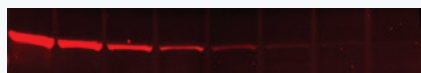
Fluorescent Western Blotting Antibodies

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 nanoparticles, excited by blue light (430–470 nm), that give a signal that is considerably brighter than conventional dyes. 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.

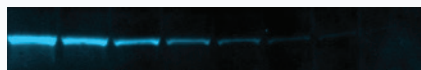
StarBright Blue Antibodies

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



▲
32 pg

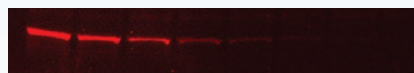
StarBright Blue 520
(blue-excited fluorescence, 20 sec)



▲
32 pg

Traditional Detection Methods

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



▲
63 pg

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



▲
500 pg

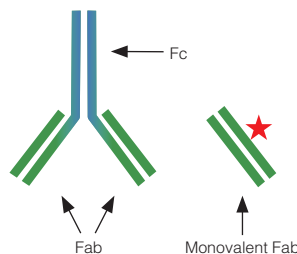
SuperSignal West Femto ECL
(horseradish peroxidase chemiluminescence, 60 sec)



▲
32 pg

hFAB Rhodamine Housekeeping Antibodies

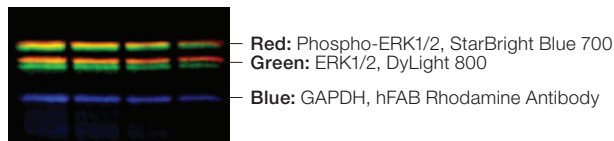
One-step detection of common housekeeping proteins like actin, tubulin, and GAPDH is possible using hFAB Housekeeping Antibodies. These fluorescent antibodies are directly labeled with a rhodamine derivative to make secondary antibodies unnecessary. hFAB Antibodies are generated with human combinatorial 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 Fluor fluorophores. Confidently use them to normalize your data with either Bio-Rad 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

Blocking Buffer

12010020 EveryBlot Blocking Buffer, 500 ml

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

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc. HuCAL is a trademark of MorphoSys AG in certain jurisdictions.

All trademarks used herein are the property of their respective owner.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website 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 4003 0399
Canada 1 905 364 3435 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 Hong Kong 852 2789 3300 Hungary 00 800 00 24 67 23 India 91 124 4029300 Israel 0 3 9636050
Italy 00 800 00 24 67 23 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Luxembourg 00 800 00 24 67 23 Mexico 52 555 488 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 00 800 00 24 67 23 Singapore 65 6415 3188 South Africa 00 800 00 24 67 23 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 66 2 651 8311 United Arab Emirates 36 1 459 6150 United Kingdom 00 800 00 24 67 23

