

Fluorescent Western Blotting with StarBright™ Blue Secondary Antibodies

Unmatched Fluorescence Sensitivity



Bio-Rad's new StarBright Blue Secondary Antibodies are labeled with novel, exceptionally bright fluorophores. These antibodies offer unparalleled sensitivity for western blotting, with exposure times in seconds instead of 5–10 minutes, compared to traditional dye-labeled antibodies.

Easy Multiplexing

The fluorophores (excitation [Ex]/emission [Em]: 440/520 and 470/700 nm) are fully compatible with Bio-Rad's new ChemiDoc™ MP Imaging System.* StarBright Blue Secondary Antibodies may be used together with:

- Stain-free gels for total protein normalization
- hFAB™ Rhodamine Antibodies (Ex/Em: 530/580 nm), which are pre-labeled housekeeping protein (HKP) primary antibodies for normalizing protein loading amounts
- Other common dyes, for example DyLight 800 (Ex/Em: 775/795 nm)

Recommendations for Multiplex Fluorescent Western Blotting

Target Description	Fluorophore**	Protein Normalization
1-plex: target of interest	StarBright Blue 700***	—
2-plex: target with protein loading normalization	StarBright Blue 700***	Total protein (stain-free) and/or HKP (hFAB Rhodamine Antibody†)
3-plex: two targets with protein loading normalization	Target 1 (low abundance): StarBright Blue 700*** Target 2 (higher abundance): StarBright Blue 520***	Total protein (stain-free) and/or HKP (hFAB Rhodamine Antibody†)

* StarBright Antibodies are also compatible with other imagers enabled with the appropriate excitation/emission capabilities.

** These fluorophores have been validated for use with Bio-Rad's ChemiDoc MP Imaging System and offer highly sensitive detection with minimal cross talk. Other fluorophores with similar excitation and emission properties may be used but optimization may be needed for best results.

*** Use StarBright B520 or B700 option under Application > Blots in Image Lab™ Touch Software, version 2.3 and later software releases.

† Use Rhodamine option under Application > Blots in Image Lab Touch Software, version 2.3 and later software releases.

Fluorescent Western Blotting with Other Fluorophores

Alternative Multiplexing Options

Number of Targets	Fluorophores		
1	DyLight 800		
2	DyLight 800	DyLight 650	
3	DyLight 800	DyLight 650	hFAB HKP

Ordering Information

Catalog # Description

StarBright Blue 520 Secondary Antibodies

12005866 Goat Anti-Mouse IgG StarBright Blue 520, 400 μ l
12005867 Goat Anti-Mouse IgG StarBright Blue 520, 80 μ l
12005869 Goat Anti-Rabbit IgG StarBright Blue 520, 400 μ l
12005870 Goat Anti-Rabbit IgG StarBright Blue 520, 80 μ l

StarBright Blue 700 Secondary Antibodies

12004158 Goat Anti-Mouse IgG StarBright Blue 700, 400 μ l
12004159 Goat Anti-Mouse IgG StarBright Blue 700, 80 μ l
12004161 Goat Anti-Rabbit IgG StarBright Blue 700, 400 μ l
12004162 Goat Anti-Rabbit IgG StarBright Blue 700, 80 μ l

Related Products

12004163 Anti-Actin hFAB Rhodamine Antibody, 200 μ l
12004164 Anti-Actin hFAB Rhodamine Antibody, 40 μ l
12004165 Anti-Tubulin hFAB Rhodamine Antibody, 200 μ l
12004166 Anti-Tubulin hFAB Rhodamine Antibody, 40 μ l
12004167 Anti-GAPDH hFAB Rhodamine Antibody, 200 μ l
12004168 Anti-GAPDH hFAB Rhodamine Antibody, 40 μ l
1610782 1x Tris Buffered Saline (TBS) with 1% Casein, 1 L
1706435 10x Tris Buffered Saline (TBS), 1 L
1610781 10% Tween 20, 1 L
1610373 Precision Plus Protein™ All Blue Standards

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

DyLight is a trademark of Thermo Fisher Scientific Inc. Tween is a trademark of ICI Americas Inc.

Precision Plus Protein Standards are sold under license from Life Technologies Corporation, Carlsbad, CA for use only by the buyer of the product. The buyer is not authorized to sell or resell this product or its components.

The BIO-RAD logo consists of the words "BIO-RAD" in a bold, white, sans-serif font, centered within a green rounded rectangular background.