

StarBright™ Blue 700 Fluorescent Secondary Antibodies

Catalog # Description

12004157	Goat Anti-Mouse IgG StarBright Blue 700, 5 x 400 µl
12004158	Goat Anti-Mouse IgG StarBright Blue 700, 400 µl
12004159	Goat Anti-Mouse IgG StarBright Blue 700, 80 µl
12004160	Goat Anti-Rabbit IgG StarBright Blue 700, 5 x 400 µl
12004161	Goat Anti-Rabbit IgG StarBright Blue 700, 400 µl
12004162	Goat Anti-Rabbit IgG StarBright Blue 700, 80 µl

Shelf life: 1 year at -20°C lyophilized; 6 months at 4°C after resuspension.

Instruction Manual

Visit bio-rad.com/web/StarBright for more detailed information about this product.

For technical support call your local Bio-Rad office. In the U.S. call 1-800-4BIORAD (1-800-424-6723).

BIO-RAD

Introduction

Bio-Rad's StarBright™ Blue 700 Secondary Antibodies are fluorescently tagged detection reagents for western blotting. These tagged antibodies are exceptionally bright and resistant to photobleaching, and they offer unparalleled sensitivity and reproducibility in western blotting.

Western blots visualized with StarBright Blue 700 Antibodies (excitation: 470 nm, emission: 700 nm) may be imaged using the fluorescence-enabled ChemiDoc™ MP Imaging System.

StarBright Blue 700 Secondary Antibodies may be used together with fluorescent antihousekeeping primary antibodies (hFAB™ Rhodamine Anti-HKP) for normalizing protein loading amounts.

Instructions for Use

Preparation

Resuspend contents of the tube in the indicated volume of distilled or deionized water and leave on ice for at least 30 min prior to use. The resuspended solution may be stored at 4°C in the dark for up to 6 months. **Do not freeze the solubilized material.**

Brief centrifugation (pulse spin for 2–3 sec at max speed in a tabletop microcentrifuge) may be used to collect the contents at the bottom of the tube. Do not centrifuge more than 10 sec.

General Guidelines

StarBright Blue 700 Secondary Antibodies may be used for detection on low fluorescence PVDF or nitrocellulose membranes. Choose LF PVDF when using any of the hFAB Rhodamine Anti-HKP Antibodies for protein loading normalization.

For blocking and washing, use 2 ml of solution per 10 cm² of membrane (15 ml for a mini gel). For primary and secondary antibody incubations, use 1 ml of solution per 10 cm² of membrane (10 ml for a mini gel). Use the smallest flat bottom tray that will accommodate the blot.

Protect the blot from light (for example, using aluminum foil) during incubations with fluorescent antibodies. Image immediately after step 5 for best results.

Do not allow the blot to dry out at any time prior to or during imaging.

We recommend using all blue protein standards (2–3 μ l) since red- or pink-colored standards may fluoresce brightly in the rhodamine imaging channel and can interfere with data acquisition. Shake or rock well (without spilling) during incubations with the StarBright 700 Secondary Antibodies.

Dilute StarBright Blue 700 Secondary Antibodies into incubation solution immediately prior to use.

StarBright Blue 700 Secondary Antibodies may be used successfully in many different immunodetection protocols. The following protocol is recommended for detection with high sensitivity, low background, and minimal nonspecific cross-reactivity.

Protocol

1. **Block:** 1 hr at room temperature (RT) with Tris buffered saline (TBS) + 1% casein.
2. **Incubate in primary antibody:** Dilute and incubate the primary antibody as specified in your protocol or by the vendor. If no protocol is provided with the primary antibody, it may be diluted in TBS + 1% casein buffer.
3. **Wash:** 5 x 5 min at RT with TBST (TBS + 0.05% Tween 20).

4. **Incubate with StarBright Blue 700 Secondary**

Antibody: Dilute secondary antibody into TBS + 1% casein. Recommended starting dilution is 1:2,500 (range: 1:2,500–1:5,000). Incubate for 1 hr at RT.

At this stage, hFAB Rhodamine Anti-HKP Primary Antibody (recommended dilution range: 1:1000–1:10,000) may be added as well for protein loading normalization.

5. **Wash:** 6 x 5 min at RT with TBST.

6. **Image** on the ChemiDoc™ MP Imaging System with Image Lab™ Touch Software. Select **StarBright B700** under **Application > Blots** (for Image Lab Touch Software, version 2.0 and later software releases).

If using an hFAB Rhodamine Anti-HKP Antibody for protein loading normalization, configure a multichannel imaging protocol using the **StarBright B700** and **Rhodamine** options on Image Lab Touch Software, version 2.0 and later software releases.

Ordering Information

Catalog # Description

StarBright Blue 700 Secondary Antibodies

12004157	Goat Anti-Mouse IgG StarBright Blue 700, 5 x 400 μ l
12004158	Goat Anti-Mouse IgG StarBright Blue 700, 400 μ l
12004159	Goat Anti-Mouse IgG StarBright Blue 700, 80 μ l
12004160	Goat Anti-Rabbit IgG StarBright Blue 700, 5 x 400 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, 500 μ l

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.

Tween is a trademark of ICI Americas Inc.

Bio-Rad Laboratories, Inc.

2000 Alfred Nobel Drive, Hercules, CA 94547 USA

510-741-1000