



Glowing products for science

Goat Anti-Mouse IgG1 (γ1)

Goat anti-mouse IgG1 isotype-specific secondary antibody labeled with our superior CF® dyes and biotin.



Product Description

This is a goat anti-mouse IgG1 isotype-specific secondary antibody labeled with our bright and photostable CF® Dyes and biotin. The conjugates are prepared from affinity-purified antibodies that react with Fc portion of the heavy chain of mouse IgG1. To minimize cross reactivity, the antibodies are cross-adsorbed against other mouse IgG subclasses (IgG2a, IgG2b, IgG3), and human, bovine and rabbit serum proteins.

- Cross-adsorbed for specific staining with minimal background
- Available in 12 bright and photostable CF® Dyes
- Biotin conjugate also available
- Suitable for western, immunofluorescence, and immunohistology

Note: Conjugates of blue fluorescent dyes like CF®350 and CF®405S are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors.

Call us : [800-304-5357](tel:800-304-5357) Email: btinfo@biotium.com

Product attributes

Antibody type	Secondary
Clonality	Polyclonal
Host species	Goat
Antibody reactivity (target)	Mouse IgG1 (γ1)
Species reactivity	Mouse
Cross adsorption	Bovine, Human, Rabbit
Concentration	2 mg/mL
Antibody/conjugate formulation	Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide, Lyophilized: PBS/15 mg/mL BSA/20 mg/mL trehalose after reconstitution
Secondary/tag antibody applications	ELISA, Flow cytometry, IHC, IF (cells or tissue sections), Western blot
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (<i>Bos taurus</i>), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

Goat Anti-Mouse IgG1

Conjugation	Ex/Em	Size	Catalog No.
CF@350	347/448 nm	50 uL (100 ug)	20245-1
		0.25 mL (500 ug)	20245
CF@405S	404/431 nm	50 uL (100 ug)	20380-50uL
		0.25 mL (500 ug)	20380-250uL
CF@488A	490/515 nm	50 uL (100 ug)	20246-1
		0.25 mL (500 ug)	20246
CF@543	541/560 nm	50 uL (100 ug)	20325-1
		0.25 mL (500 ug)	20325
CF@555	555/565 nm	50 uL (100 ug)	20247-1
		0.25 mL (500 ug)	20247
CF@568	562/583 nm	50 uL (100 ug)	20248-1
		0.25 mL (500 ug)	20248
CF@594	593/614 nm	50 uL (100 ug)	20249-1
		0.25 mL (500 ug)	20249
CF@633	630/650 nm	50 uL (100 ug)	20250-1
		0.25 mL (500 ug)	20250
CF@640R	642/662 nm	50 uL (100 ug)	20251-1
		0.25 mL (500 ug)	20251
CF@647	650/665 nm	50 uL (100 ug)	20252-1
		0.25 mL (500 ug)	20252
CF@680	681/698 nm	50 uL (100 ug)	20253-1
		0.25 mL (500 ug)	20253
CF@770	770/797 nm	50 uL (100 ug)	20254-1
		0.25 mL (500 ug)	20254
Biotin	N/A	50 uL (100 ug)	20471-50uL
		0.25 mL (500 ug)	20471-250uL

View our full selection of bright and specific [Secondary Antibodies](#), or search our catalog using our [Antibody Finder](#). Alternatively, you can view our [secondary antibody product listings](#) with catalog numbers.

CF® Dyes offer exceptional brightness and photostability. For more information see our [CF® Dye technology page](#).

Storage and Handling

Liquid format: Store at -20°C, protected from light. Product is stable for at least 6 months from date of receipt when stored as recommended. Liquid format antibodies contain 50% glycerol and will not freeze at -20°C.

Lyophilized format: Store at -20°C, protected from light. Product is stable for at least 6 months from date of receipt when stored as recommended. Reconstitute antibodies in water using the indicated volumes below:

CF® Dye and biotin conjugates: add 0.5 mL dH₂O

HRP or DNP conjugates: add 1 mL dH₂O

Add the indicated volume of water directly to the vial containing the lyophilized antibody and mix gently to dissolve. Store reconstituted antibody at -20°C and protect from light. Aliquot to avoid repeated freeze/thaw cycles. Alternatively, an equal volume of glycerol can be mixed with the reconstituted antibody so that it will remain liquid at -20°C.

Optional: A preservative such as 0.05% sodium azide (final concentration) can be added to CF® Dye and biotin conjugates. Do not add sodium azide to HRP conjugates.

Note: Storage of the antibody for more than a day at final working dilution is not recommended.

CF is a registered trademark of Biotium, Inc.

References

Download a list of curated [CF® Dye references](#).

This datasheet was generated on November 21, 2024 at 07:47:18 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/goat-anti-mouse-igg1/>