



Glowing products for science

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

## 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.

### 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 (Bos taurus), 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)	<a href="#">20245-1</a>
		0.25 mL (500 ug)	<a href="#">20245</a>
CF®405S	404/431 nm	50 uL (100 ug)	<a href="#">20380-50uL</a>
		0.25 mL (500 ug)	<a href="#">20380-250uL</a>
CF®488A	490/515 nm	50 uL (100 ug)	<a href="#">20246-1</a>
		0.25 mL (500 ug)	<a href="#">20246</a>
CF®543	541/560 nm	50 uL (100 ug)	<a href="#">20325-1</a>
		0.25 mL (500 ug)	<a href="#">20325</a>
CF®555	555/565 nm	50 uL (100 ug)	<a href="#">20247-1</a>
		0.25 mL (500 ug)	<a href="#">20247</a>
CF®568	562/583 nm	50 uL (100 ug)	<a href="#">20248-1</a>
		0.25 mL (500 ug)	<a href="#">20248</a>
CF®594	593/614 nm	50 uL (100 ug)	<a href="#">20249-1</a>
		0.25 mL (500 ug)	<a href="#">20249</a>
CF®633	630/650 nm	50 uL (100 ug)	<a href="#">20250-1</a>
		0.25 mL (500 ug)	<a href="#">20250</a>
CF®640R	642/662 nm	50 uL (100 ug)	<a href="#">20251-1</a>
		0.25 mL (500 ug)	<a href="#">20251</a>
CF®647	650/665 nm	50 uL (100 ug)	<a href="#">20252-1</a>
		0.25 mL (500 ug)	<a href="#">20252</a>
CF®680	681/698 nm	50 uL (100 ug)	<a href="#">20253-1</a>
		0.25 mL (500 ug)	<a href="#">20253</a>
CF®770	770/797 nm	50 uL (100 ug)	<a href="#">20254-1</a>
		0.25 mL (500 ug)	<a href="#">20254</a>
Biotin	N/A	50 uL (100 ug)	<a href="#">20471-50uL</a>
		0.25 mL (500 ug)	<a href="#">20471-250uL</a>

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 dH<sub>2</sub>O

HRP or DNP conjugates: add 1 mL dH<sub>2</sub>O

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 August 29, 2025 at 01:47:13 PM. Visit product page to check for updated information before use.

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