



Glowing products for science

## Goat Anti-Mouse IgG2b (γ2b)

Goat anti-mouse IgG2b isotype-specific secondary antibody labeled with our superior CF® Dyes and biotin.



### Product Description

This is a goat anti-mouse IgG2b 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 IgG2b. To minimize cross reactivity, the antibodies are cross-adsorbed against other mouse IgG subclasses (IgG1, IgG2a, IgG3), and human, bovine and rabbit serum proteins.

- Cross-adsorbed for specific staining with minimal background
- Available in 13 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](mailto:btinfo@biotium.com)

### Product attributes

<b>Antibody type</b>	Secondary
<b>Clonality</b>	Polyclonal
<b>Host species</b>	Goat
<b>Antibody reactivity (target)</b>	Mouse IgG2b (γ2b)
<b>Species reactivity</b>	Mouse
<b>Cross adsorption</b>	Bovine, Human, Rabbit
<b>Concentration</b>	2 mg/mL
<b>Antibody/conjugate formulation</b>	Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide, Lyophilized: PBS/15 mg/mL BSA/20 mg/mL trehalose after reconstitution
<b>Secondary/tag antibody applications</b>	ELISA, Flow cytometry, IHC, IF (cells or tissue sections), Western blot
<b>Product origin</b>	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 IgG2b

Conjugation	Ex/Em	Size	Catalog No.
CF®350	347/448 nm	50 uL (100 ug)	<a href="#">20265-1</a>
		0.25 mL (500 ug)	<a href="#">20265</a>
CF®405S	404/431 nm	50 uL (100 ug)	<a href="#">20382-50uL</a>
		0.25 mL (500 ug)	<a href="#">20382-250uL</a>
CF®488A	490/515 nm	50 uL (100 ug)	<a href="#">20266-1</a>
		0.25 mL (500 ug)	<a href="#">20266</a>
CF®543	541/560 nm	50 uL (100 ug)	<a href="#">20327-1</a>
		0.25 mL (500 ug)	<a href="#">20327</a>
CF®555	555/565 nm	50 uL (100 ug)	<a href="#">20267-1</a>
		0.25 mL (500 ug)	<a href="#">20267</a>
CF®568	562/583 nm	50 uL (100 ug)	<a href="#">20268-1</a>
		0.25 mL (500 ug)	<a href="#">20268</a>
CF®594	593/614 nm	50 uL (100 ug)	<a href="#">20269-1</a>
		0.25 mL (500 ug)	<a href="#">20269</a>
CF®633	630/650 nm	50 uL (100 ug)	<a href="#">20270-1</a>
		0.25 mL (500 ug)	<a href="#">20270</a>
CF®640R	642/662 nm	50 uL (100 ug)	<a href="#">20271-1</a>
		0.25 mL (500 ug)	<a href="#">20271</a>
CF®647	650/665 nm	50 uL (100 ug)	<a href="#">20272-1</a>
		0.25 mL (500 ug)	<a href="#">20272</a>
CF®680	681/698 nm	50 uL (100 ug)	<a href="#">20273-1</a>
		0.25 mL (500 ug)	<a href="#">20273</a>
CF®750	755/777 nm	50 uL (100 ug)	<a href="#">20430-50uL</a>
		250 uL (100 ug)	<a href="#">20430-250uL</a>
CF®770	770/797 nm	50 uL (100 ug)	<a href="#">20274-1</a>
		0.25 mL (500 ug)	<a href="#">20274</a>
Biotin	N/A	50 uL (100 ug)	<a href="#">20473-50uL</a>
		0.25 mL (500 ug) (500 ug)	<a href="#">20473-250uL</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 [CF® dye references](#).

This datasheet was generated on December 3, 2024 at 09:39:33 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/goat-anti-mouse-igg2b-2b/>