

# 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. 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.

## Product attributes

Antibody type	Secondary
Clonality	Polyclonal
Host species	Goat
Antibody reactivity (target)	Mouse IgG2b (γ2b)
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

## References

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

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>



