

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 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.

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)	20265-1
		0.25 mL (500 ug)	20265
CF®405S	404/431 nm	50 uL (100 ug)	20382-50uL
		0.25 mL (500 ug)	20382-250uL
CF®488A	490/515 nm	50 uL (100 ug)	20266-1
		0.25 mL (500 ug)	20266
CF®543	541/560 nm	50 uL (100 ug)	20327-1
		0.25 mL (500 ug)	20327
CF®555	555/565 nm	50 uL (100 ug)	20267-1
		0.25 mL (500 ug)	20267
CF®568	562/583 nm	50 uL (100 ug)	20268-1
		0.25 mL (500 ug)	20268
CF®594	593/614 nm	50 uL (100 ug)	20269-1
		0.25 mL (500 ug)	20269
CF®633	630/650 nm	50 uL (100 ug)	20270-1
		0.25 mL (500 ug)	20270
CF®640R	642/662 nm	50 uL (100 ug)	20271-1
		0.25 mL (500 ug)	20271
CF®647	650/665 nm	50 uL (100 ug)	20272-1
		0.25 mL (500 ug)	20272
CF®680	681/698 nm	50 uL (100 ug)	20273-1
		0.25 mL (500 ug)	20273
CF®750	755/777 nm	50 uL (100 ug)	20430-50uL
		250 uL (100 ug)	20430-250uL
CF®770	770/797 nm	50 uL (100 ug)	20274-1
		0.25 mL (500 ug)	20274
Biotin	N/A	50 uL (100 ug)	20473-50uL
		0.25 mL (500 ug) (500 ug)	20473-250uL

This datasheet was generated on July 17, 2024 at 09:37:03 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/goat-anti-mouse-igg2b-2b/>

