



Glowing products for science

Goat Anti-Mouse IgG3

Goat anti-mouse IgG3 isotype-specific secondary antibody labeled with CF® Dyes.



Product Description

This is a goat anti-mouse IgG3 isotype-specific secondary antibody labeled with a selection of CF® Dyes and biotin. The conjugate is prepared from affinity-purified antibody that reacts with the Fc portion of the heavy chain of mouse IgG3. To minimize cross reactivity, the antibody is cross-adsorbed against other mouse IgG subclasses (IgG1, IgG2a, IgG2b, IgG2c), and human, bovine and rabbit serum proteins.

- Cross-adsorbed for specific staining with minimal background
- Available with 12 CF® Dyes and biotin
- Suitable for western, immunofluorescence, and immunohistology

Advantages of CF® Dyes

- Exceptionally bright and stable
- Highly water soluble
- Superior signal-to-noise
- Ideal for flow cytometry and microscopy
- CF®680 and CF®770 are ideal for near-IR Western/In Cell Western®
- Options validated in [super-resolution STORM and photoacoustic imaging](#)

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

Product attributes

Clonality	Polyclonal
Antibody type	Secondary
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
Species reactivity	Mouse
Host species	Goat
Antibody reactivity (target)	Mouse IgG3
Cross adsorption	Bovine, Human, Rabbit
Conjugation	CF®350, CF®405S, CF®488A, CF®543, CF®555, CF®594, CF®633, CF®640R, CF®647, CF®680, CF®750, CF®770, Biotin
Secondary/tag antibody applications	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.

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF®350	347/448 nm	50 uL (100 ug)	20922-50uL	CF®350 Features
		0.25 mL (500 ug)	20922-250uL	
CF®405S	404/431 nm	50 uL (100 ug)	20923-50uL	CF®405S Features
		0.25 mL (500 ug)	20923-250uL	
CF®488A	490/515 nm	50 uL (100 ug)	20924-50uL	CF®488A Features
		0.25 mL (500 ug)	20924-250uL	
CF®543	541/560 nm	50 uL (100 ug)	20925-50uL	CF®543 Features
		0.25 mL (500 ug)	20925-250uL	
CF®555	555/565 nm	50 uL (100 ug)	20926-50uL	CF®555 Features
		0.25 mL (500 ug)	20926-250uL	
CF®594	593/614 nm	50 uL (100 ug)	20927-50uL	CF®594 Features
		0.25 mL (500 ug)	20927-250uL	
CF®633	630/650 nm	50 uL (100 ug)	20928-50uL	CF®633 Features
		0.25 mL (500 ug)	20928-250uL	
CF®640R	642/662 nm	50 uL (100 ug)	20929-50uL	CF®640R Features
		0.25 mL (500 ug)	20929-250uL	
CF®647	650/665 nm	50 uL (100 ug)	20930-50uL	CF®647 Features
		0.25 mL (500 ug)	20930-250uL	
CF®680	681/698 nm	50 uL (100 ug)	20931-50uL	CF®680 Features
		0.25 mL (500 ug)	20931-250uL	
CF®750	755/777 nm	50 uL (100 ug)	20431-50uL	CF®750 Features
		250 uL (100 ug)	20431-250uL	
CF®770	770/797 nm	50 uL (100 ug)	20932-50uL	CF®770 Features
		0.25 mL (500 ug)	20932-250uL	
Biotin	N/A	50 uL (100 ug)	20921-50uL	
		0.25 mL (500 ug) (500 ug)	20921-250uL	

[Browse Secondary Antibodies](#), search using our [Antibody Finder](#), or see all [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. Alexa Fluor and DyLight are registered trademarks of Thermo Fisher Scientific. Cy Dye is a registered trademark of Cytiva.

References

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

This datasheet was generated on July 9, 2025 at 11:35:54 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/cf750-goat-anti-mouse-igg3/>