



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)

### Product attributes

<b>Clonality</b>	Polyclonal
<b>Antibody type</b>	Secondary
<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>Species reactivity</b>	Mouse
<b>Host species</b>	Goat
<b>Antibody reactivity (target)</b>	Mouse IgG3
<b>Cross adsorption</b>	Bovine, Human, Rabbit
<b>Conjugation</b>	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
<b>Secondary/tag antibody applications</b>	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 ( <i>Bos taurus</i> ), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

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

[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 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. 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 June 3, 2026 at 03:58:50 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cf750-goat-anti-mouse-igg3/>