

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](#)

[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 November 22, 2024 at 11:59:47 PM. Visit product page to check for updated information before use.

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

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 (<i>Bos taurus</i>), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.