

## Goat Anti-Mouse IgG1 F(ab'), Fc Gamma Specific

Goat anti-mouse IgG1 F(ab'), Fc gamma fragment specific secondary antibody labeled with our superior CF® dyes.



### Product attributes

Clonality	Polyclonal
Cross adsorption	Not cross-adsorbed
Antibody type	Secondary
Concentration	2 mg/mL
Antibody/conjugate formulation	Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide
Species reactivity	Mouse
Host species	Goat
Antibody reactivity (target)	Mouse IgG1, Fc
Secondary/tag antibody applications	Flow cytometry, IF (cells or tissue sections), Western blot

### Product Description

This is a goat anti-mouse IgG1 F(ab'), Fc gamma fragment specific secondary antibody labeled with our superior CF® dyes.

- Available with 2 bright and photostable CF® dyes
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

View our full selection of [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.

### References

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

Conjugation	Ex/Em	Size	Catalog No.
CF®488A	490/515 nm	50 uL (100 ug)	<a href="#">20872-50uL</a>
CF®594	593/614 nm	50 uL (100 ug)	<a href="#">20873-50uL</a>

This datasheet was generated on June 30, 2024 at 09:30:01 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/goat-anti-mouse-igg1-fab-fragment-gamma-specific/>