

## 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 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

### Product attributes

<b>Clonality</b>	Polyclonal
<b>Cross adsorption</b>	Not cross-adsorbed
<b>Antibody type</b>	Secondary
<b>Concentration</b>	0.5 mg/mL
<b>Antibody/conjugate formulation</b>	Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide
<b>Species reactivity</b>	Mouse
<b>Host species</b>	Goat
<b>Antibody reactivity (target)</b>	Mouse IgG1, Fc
<b>Secondary/tag antibody applications</b>	Flow cytometry, 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.

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

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>

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](#).

This datasheet was generated on June 3, 2026 at 03:14:10 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/goat-anti-mouse-igg1-fab-fragment-gamma-specific/>