



Glowing products for science

## Goat Anti-Mouse IgG (H+L), F(ab')<sub>2</sub> Fragment

Goat anti-mouse IgG (H+L), F(ab')<sub>2</sub> fragment secondary antibody labeled with our superior CF® Dyes and other labels.



### Product Description

This is a goat anti-mouse IgG (H L), F(ab')<sub>2</sub> fragment secondary antibody labeled with our superior CF® Dyes and other labels.

- Available in 14 bright and photostable CF® Dyes, and HRP
- Biotin, APC, and R-PE conjugates also available
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

**Note:** Conjugates of blue fluorescent dyes like CF®350 are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors.

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

### Product attributes

<b>Antibody type</b>	Secondary
<b>Clonality</b>	Polyclonal
<b>Host species</b>	Goat
<b>Antibody reactivity (target)</b>	Mouse IgG
<b>Species reactivity</b>	Mouse
<b>Cross adsorption</b>	Not cross-adsorbed
<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>Secondary/tag antibody applications</b>	ELISA, 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.

# Goat Anti-Mouse IgG (H+L), F(ab')<sub>2</sub> Fragment

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
<a href="#">CF@350</a>	347/448 nm	50 uL (100 ug)	<a href="#">20145-1</a>	<a href="#">CF@350 Features</a>
		0.25 mL (500 ug)	<a href="#">20145</a>	
<a href="#">CF@488A</a>	490/515 nm	50 uL (100 ug)	<a href="#">20011-1</a>	<a href="#">CF@488A Features</a>
		0.25 mL (500 ug)	<a href="#">20011</a>	
<a href="#">CF@505</a>	505/519 nm	50 uL (100 ug)	<a href="#">20916-50uL</a>	
		0.25 mL (500 ug)	<a href="#">20916-250uL</a>	
<a href="#">CF@543</a>	541/560 nm	50 uL (100 ug)	<a href="#">20329-1</a>	<a href="#">CF@543 Features</a>
		0.25 mL (500 ug)	<a href="#">20329</a>	
<a href="#">CF@555</a>	555/565 nm	50 uL (100 ug)	<a href="#">20032-1</a>	<a href="#">CF@555 Features</a>
		0.25 mL (500 ug)	<a href="#">20032</a>	
<a href="#">CF@568</a>	562/583 nm	50 uL (100 ug)	<a href="#">20109-1</a>	<a href="#">CF@568 Features</a>
		0.25 mL (500 ug)	<a href="#">20109</a>	
<a href="#">CF@583R</a>	585/609 nm	50 uL (100 ug)	<a href="#">20914-50uL</a>	<a href="#">CF@583R Features</a>
		0.25 mL (500 ug)	<a href="#">20914-250uL</a>	
<a href="#">CF@594</a>	593/614 nm	50 uL (100 ug)	<a href="#">20119-1</a>	<a href="#">CF@594 Features</a>
		0.25 mL (500 ug)	<a href="#">20119</a>	
<a href="#">CF@633</a>	630/650 nm	50 uL (100 ug)	<a href="#">20130-1</a>	<a href="#">CF@633 Features</a>
		0.25 mL (500 ug)	<a href="#">20130</a>	
<a href="#">CF@640R</a>	642/662 nm	50 uL (100 ug)	<a href="#">20086-1</a>	<a href="#">CF@640R Features</a>
		0.25 mL (500 ug)	<a href="#">20086</a>	
<a href="#">CF@647</a>	650/665 nm	50 uL (100 ug)	<a href="#">20042-1</a>	<a href="#">CF@647 Features</a>
		0.25 mL (500 ug)	<a href="#">20042</a>	
<a href="#">CF@660C</a>	667/685 nm	50 uL (100 ug)	<a href="#">20918-50uL</a>	<a href="#">CF@660C Features</a>
		0.25 mL (500 ug)	<a href="#">20918-250uL</a>	
<a href="#">CF@680</a>	681/698 nm	50 uL (100 ug)	<a href="#">20063-1</a>	<a href="#">CF@680 Features</a>
		0.25 mL (500 ug)	<a href="#">20063</a>	
<a href="#">CF@750</a>	755/779 nm	50 uL (100 ug)	<a href="#">20920-50uL</a>	<a href="#">CF@750 Features</a>
		0.25 mL (500 ug)	<a href="#">20920-250uL</a>	
R-PE	496, 546, 565/578 nm	200 ul (100 ug)	<a href="#">20357-200uL</a>	
		500 ul (250 ug)	<a href="#">20357-500uL</a>	
APC	650/660 nm	100 ul (50 ug)	<a href="#">20414-100uL</a>	
		500 ul (250 ug)	<a href="#">20414-500uL</a>	
Biotin	N/A	50 uL (100 ug)	<a href="#">20187-1</a>	
		0.25 mL (500 ug)	<a href="#">20187</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 November 21, 2024 at 02:54:42 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/goat-anti-mouse-igg-hlfab2-fragment/>