



Glowing products for science

## Goat Anti-Guinea Pig IgG (H+L)

Goat anti-guinea pig IgG (H L) secondary antibody labeled with our superior CF® Dyes.



### Product Description

This is a goat anti-guinea pig IgG (H L) secondary antibody labeled with our superior CF® Dyes.

- Available in 10 bright and photostable CF® Dyes
- 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)

### Product attributes

<b>Antibody type</b>	Secondary
<b>Clonality</b>	Polyclonal
<b>Host species</b>	Goat
<b>Antibody reactivity (target)</b>	Guinea Pig IgG
<b>Species reactivity</b>	Guinea pig
<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>	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 (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

# Goat Anti-Guinea Pig IgG (H+L)

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF@350	347/448 nm	50 uL (100 ug)	<a href="#">20198-1</a>	<a href="#">CF@350 Features</a>
0.5 mL (1 mg)	<a href="#">20198</a>			
1 mg	<a href="#">20198-1mg</a>			
CF@488A	490/515 nm	50 uL (100 ug)	<a href="#">20017-1</a>	<a href="#">CF@488A Features</a>
0.5 mL (1 mg)	<a href="#">20017</a>			
1 mg	<a href="#">20017-1mg</a>			
CF@543	541/560 nm	50 uL (100 ug)	<a href="#">20317-1</a>	<a href="#">CF@543 Features</a>
0.5 mL (1 mg)	<a href="#">20317</a>			
1 mg	<a href="#">20317-1mg</a>			
CF@555	555/565 nm	50 uL (100 ug)	<a href="#">20036-1</a>	<a href="#">CF@555 Features</a>
0.5 mL (1 mg)	<a href="#">20036</a>			
1 mg	<a href="#">20036-1mg</a>			
CF@568	562/583 nm	50 uL (100 ug)	<a href="#">20108-1</a>	<a href="#">CF@568 Features</a>
0.5 mL (1 mg)	<a href="#">20108</a>			
1 mg	<a href="#">20108-1mg</a>			
CF@583R	585/609 nm	50 uL (100 ug)	<a href="#">20899-50uL</a>	<a href="#">CF@583R Features</a>
0.5 mL (1 mg)	<a href="#">20899-500uL</a>			
CF@594	593/614 nm	50 uL (100 ug)	<a href="#">20118-1</a>	<a href="#">CF@594 Features</a>
0.5 mL (1 mg)	<a href="#">20118</a>			
1 mg	<a href="#">20118-1mg</a>			
CF@633	630/650 nm	50 uL (100 ug)	<a href="#">20129-1</a>	<a href="#">CF@633 Features</a>
0.5 mL (1 mg)	<a href="#">20129</a>			
1 mg	<a href="#">20129-1mg</a>			
CF@640R	642/662 nm	50 uL (100 ug)	<a href="#">20085-1</a>	<a href="#">CF@640R Features</a>
0.5 mL (1 mg)	<a href="#">20085</a>			
1 mg	<a href="#">20085-1mg</a>			
CF@647	650/665 nm	50 uL (100 ug)	<a href="#">20041-1</a>	<a href="#">CF@647 Features</a>
0.5 mL (1 mg)	<a href="#">20041</a>			
1 mg	<a href="#">20041-1mg</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 curated [CF® Dye references](#).

This datasheet was generated on June 3, 2026 at 04:39:22 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/goat-anti-guinea-pig-igg-hl/>