



Glowing products for science

## Donkey Anti-Guinea Pig IgG (H+L), Highly Cross-Adsorbed

Highly cross-adsorbed donkey anti-guinea pig IgG (H L) secondary antibody labeled with our superior CF® Dyes.



### Product Description

This is a highly cross-adsorbed donkey anti-guinea pig IgG (H L) secondary antibody labeled with our bright and photostable CF® Dyes. To minimize cross-reactivity, the antibody has been adsorbed against bovine, chicken, goat, Syrian hamster, horse, human, mouse, rabbit, and sheep serum.

- Highly cross-adsorbed for specific staining with minimal background
- Available in 12 bright and photostable CF® Dyes
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

**Note:** Conjugates of blue fluorescent dyes like CF®405M 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>	Donkey
<b>Antibody reactivity (target)</b>	Guinea Pig IgG
<b>Species reactivity</b>	Guinea pig
<b>Cross adsorption</b>	Bovine, Chicken, Goat, Horse, Human, Mouse, Rabbit, Rat, Sheep, Syrian hamster
<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 ( <i>Bos taurus</i> ), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

# Donkey Anti-Guinea Pig IgG (H+L), Highly Cross-Adsorbed

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF@405S	404/431 nm	50 uL (100 ug)	<a href="#">20356-50uL</a>	<a href="#">CF@405S Features</a>
		0.5 mL (1 mg)	<a href="#">20356-500uL</a>	
		1 mg	<a href="#">20356-1mg</a>	
CF@405M	408/452 nm	50 uL (100 ug)	<a href="#">20376-50uL</a>	<a href="#">CF@405M Features</a>
		0.5 mL (1 mg)	<a href="#">20376-500uL</a>	
		1 mg	<a href="#">20376-1mg</a>	
CF@488A	490/515 nm	50 uL (100 ug)	<a href="#">20169-1</a>	<a href="#">CF@488A Features</a>
		0.5 mL (1 mg)	<a href="#">20169</a>	
		1 mg	<a href="#">20169-1mg</a>	
CF@543	541/560 nm	50 uL (100 ug)	<a href="#">20316-1</a>	<a href="#">CF@543 Features</a>
		0.5 mL (1 mg)	<a href="#">20316</a>	
		1 mg	<a href="#">20316-1mg</a>	
CF@555	555/565 nm	50 uL (100 ug)	<a href="#">20276-1</a>	<a href="#">CF@555 Features</a>
		0.5 mL (1 mg)	<a href="#">20276</a>	
		1 mg	<a href="#">20276-1mg</a>	
CF@568	562/583 nm	50 uL (100 ug)	<a href="#">20377-50uL</a>	<a href="#">CF@568 Features</a>
		0.5 mL (1 mg)	<a href="#">20377-500uL</a>	
		1 mg	<a href="#">20377-1mg</a>	
CF@583R	585/609 nm	50 uL (100 ug)	<a href="#">20893-50uL</a>	<a href="#">CF@583R Features</a>
		0.5 mL (1 mg)	<a href="#">20893-500uL</a>	
CF@594	593/614 nm	50 uL (100 ug)	<a href="#">20170-1</a>	<a href="#">CF@594 Features</a>
		0.5 mL (1 mg)	<a href="#">20170</a>	
		1 mg	<a href="#">20170-1mg</a>	
CF@633	630/650 nm	50 uL (100 ug)	<a href="#">20171-1</a>	<a href="#">CF@633 Features</a>
		0.5 mL (1 mg)	<a href="#">20171</a>	
		1 mg	<a href="#">20171-1mg</a>	
CF@660C	667/685 nm	50 uL (100 ug)	<a href="#">20372-50uL</a>	<a href="#">CF@660C Features</a>
		0.5 mL (1 mg)	<a href="#">20372-500uL</a>	
		1 mg	<a href="#">20372-1mg</a>	
CF@680	681/698 nm	50 uL (100 ug)	<a href="#">20241-1</a>	<a href="#">CF@680 Features</a>
		0.25 mL (500 ug)	<a href="#">20241</a>	
CF@770	770/797 nm	50 uL (100 ug)	<a href="#">20242-1</a>	<a href="#">CF@770 Features</a>
		0.25 mL (500 ug)	<a href="#">20242</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 November 21, 2024 at 02:51:05 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/donkey-anti-guinea-pig-igg-hl-highly-cross-adsorbed/>