

## Goat Anti-Llama IgG (H+L)

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



### Product Description

This is a goat anti-llama IgG (H L) secondary antibody labeled with our superior CF® Dyes, HRP, or DNP.

- Available in 14 bright and photostable CF® Dyes, HRP, or DNP
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

**Note:** Conjugates of blue fluorescent dyes like CF®405S 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. 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](#).

### Product attributes

Clonality	Polyclonal
Antibody type	Secondary
Host species	Goat
Antibody reactivity (target)	Llama IgG
Species reactivity	Llama
Cross adsorption	Not cross-adsorbed
Antibody/conjugate formulation	Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide, Lyophilized: PBS/15 mg/mL BSA/20 mg/mL trehalose after reconstitution, HRP conjugates: PBS/50% glycerol/15 mg/mL BSA, HRP conjugates (lyophilized): PBS/10 mg/mL trehalose after reconstitution
Concentration	1 mg/mL (HRP conjugate), 2 mg/mL
Secondary/tag antibody applications	ELISA, Flow cytometry, IHC, IF (cells or tissue sections), Western blot

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
<a href="#">CF@405S</a>	404/431 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	<a href="#">20844-50uL</a> <a href="#">20844-500uL</a> <a href="#">20844-1mg</a>	<a href="#">CF@405S Features</a>
<a href="#">CF@488A</a>	490/515 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	<a href="#">20845-50uL</a> <a href="#">20845-500uL</a> <a href="#">20845-1mg</a>	<a href="#">CF@488A Features</a>
<a href="#">CF@543</a>	541/560 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	<a href="#">20846-50uL</a> <a href="#">20846-500uL</a> <a href="#">20846-1mg</a>	<a href="#">CF@543 Features</a>
<a href="#">CF@555</a>	555/565 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	<a href="#">20847-50uL</a> <a href="#">20847-500uL</a> <a href="#">20847-1mg</a>	<a href="#">CF@555 Features</a>
<a href="#">CF@568</a>	562/583 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	<a href="#">20848-50uL</a> <a href="#">20848-500uL</a> <a href="#">20848-1mg</a>	<a href="#">CF@568 Features</a>
<a href="#">CF@583R</a>	585/609 nm	50 uL (100 ug) 0.5 mL (1 mg)	<a href="#">20902-50uL</a> <a href="#">20902-500uL</a>	<a href="#">CF@583R Features</a>
<a href="#">CF@594</a>	593/614 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	<a href="#">20849-50uL</a> <a href="#">20849-500uL</a> <a href="#">20849-1mg</a>	<a href="#">CF@594 Features</a>
<a href="#">CF@640R</a>	642/662 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	<a href="#">20850-50uL</a> <a href="#">20850-500uL</a> <a href="#">20850-1mg</a>	<a href="#">CF@640R Features</a>
<a href="#">CF@647</a>	650/665 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	<a href="#">20851-50uL</a> <a href="#">20851-500uL</a> <a href="#">20851-1mg</a>	<a href="#">CF@647 Features</a>
<a href="#">CF@660R</a>	663/682 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	<a href="#">20853-50uL</a> <a href="#">20853-500uL</a> <a href="#">20853-1mg</a>	<a href="#">CF@660R Features</a>
<a href="#">CF@660C</a>	667/685 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	<a href="#">20852-50uL</a> <a href="#">20852-500uL</a> <a href="#">20852-1mg</a>	<a href="#">CF@660C Features</a>
<a href="#">CF@680</a>	681/698 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	<a href="#">20854-50uL</a> <a href="#">20854-500uL</a> <a href="#">20854-1mg</a>	<a href="#">CF@680 Features</a>
<a href="#">CF@680R</a>	680/701 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	<a href="#">20855-50uL</a> <a href="#">20855-500uL</a> <a href="#">20855-1mg</a>	<a href="#">CF@680R Features</a>
<a href="#">CF@750</a>	755/777 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	<a href="#">20856-50uL</a> <a href="#">20856-500uL</a> <a href="#">20856-1mg</a>	<a href="#">CF@750 Features</a>
<a href="#">DNP</a>	N/A	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	<a href="#">20983-50uL</a> <a href="#">20983-500uL</a> <a href="#">20983-1mg</a>	
<a href="#">HRP</a>	N/A	100 uL (100 ug) 1 mL (1 mg) 1 mg (lyophilized)	<a href="#">20475-100uL</a> <a href="#">20475-1mL</a> <a href="#">20475-1mg</a>	

This datasheet was generated on July 17, 2024 at 09:35:32 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/goat-anti-llama-igg-hl/>