

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

Goat anti-swine IgG (H L) secondary antibody labeled with our superior CF® Dyes and other labels.



### Product Description

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

- Available in 9 bright and photostable CF® Dyes
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

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.

### Product attributes

Antibody type	Secondary
Clonality	Polyclonal
Host species	Goat
Antibody reactivity (target)	Swine IgG
Species reactivity	Swine
Cross adsorption	Not cross-adsorbed
Concentration	2 mg/mL
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
Secondary/tag antibody applications	Flow cytometry, IHC, IF (cells or tissue sections), Western blot

### References

Download a list of [CF® dye references](#).

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF®488A	490/515 nm	50 uL (100 ug)	<a href="#">20028-1</a>	<a href="#">CF®488A Features</a>
		0.5 mL (1 mg)	<a href="#">20028</a>	
		1 mg	<a href="#">20028-1mg</a>	
CF®543	541/560 nm	50 uL (100 ug)	<a href="#">20324-1</a>	<a href="#">CF®543 Features</a>
		0.5 mL (1 mg)	<a href="#">20324</a>	
		1 mg	<a href="#">20324-1mg</a>	
CF®555	555/565 nm	50 uL (100 ug)	<a href="#">20236-1</a>	<a href="#">CF®555 Features</a>
		0.5 mL (1 mg)	<a href="#">20236</a>	
		1 mg	<a href="#">20236-1mg</a>	
CF®568	562/583 nm	50 uL (100 ug)	<a href="#">20091-1</a>	<a href="#">CF®568 Features</a>
		0.5 mL (1 mg)	<a href="#">20091</a>	
		1 mg	<a href="#">20091-1mg</a>	
CF®583R	585/609 nm	50 uL (100 ug)	<a href="#">20907-50uL</a>	<a href="#">CF®583R Features</a>
		0.5 mL (1 mg)	<a href="#">20907-500uL</a>	
		1 mg	<a href="#">20907-500uL</a>	
CF®594	593/614 nm	50 uL (100 ug)	<a href="#">20160-1</a>	<a href="#">CF®594 Features</a>
		0.5 mL (1 mg)	<a href="#">20160</a>	
		1 mg	<a href="#">20160-1mg</a>	
CF®633	630/650 nm	50 uL (100 ug)	<a href="#">20138-1</a>	<a href="#">CF®633 Features</a>
		0.5 mL (1 mg)	<a href="#">20138</a>	
		1 mg	<a href="#">20138-1mg</a>	
CF®640R	642/662 nm	50 uL (100 ug)	<a href="#">20089-1</a>	<a href="#">CF®640R Features</a>
		0.5 mL (1 mg)	<a href="#">20089</a>	
		1 mg	<a href="#">20089-1mg</a>	
CF®647	650/665 nm	50 uL (100 ug)	<a href="#">20286-1</a>	<a href="#">CF®647 Features</a>
		0.5 mL (1 mg)	<a href="#">20286</a>	
		1 mg	<a href="#">20286-1mg</a>	