



Glowing products for science

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

Call us : [800-304-5357](tel:800-304-5357)

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
Product origin	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-Swine IgG (H+L)

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

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 dH₂O

HRP or DNP conjugates: add 1 mL dH₂O

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 June 3, 2026 at 03:58:59 PM. Visit product page to check for updated information before use.

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