



Glowing products for science

Chicken Anti-Goat IgG (H+L)

Chicken anti-goat IgG (H L) secondary antibody labeled with our superior CF® Dyes or HRP.



Product Description

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

- Available in 7 bright and photostable CF® Dyes, and HRP
- 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

Antibody type	Secondary
Clonality	Polyclonal
Host species	Chicken
Antibody reactivity (target)	Goat IgG
Species reactivity	Goat
Cross adsorption	Not cross-adsorbed
Concentration	2 mg/mL (CF® dye conjugates), 1 mg/mL (HRP, AP conjugates)
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
Secondary/tag antibody applications	ELISA, 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.

Chicken Anti-Goat IgG (H+L)

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF@350	347/448 nm	50 uL (100 ug)	20364-50uL	CF@350 Features
0.5 mL (1 mg)	20364-500uL			
1 mg (lyophilized)	20364-1mg			
CF@488A	490/515 nm	50 uL (100 ug)	20225-1	CF@488A Features
0.5 mL (1 mg)	20225			
1 mg (lyophilized)	20225-1mg			
CF@543	541/560 nm	50 uL (100 ug)	20333-50uL	CF@543 Features
0.5 mL (1 mg)	20333-500uL			
1 mg (lyophilized)	20333-1mg			
CF@568	562/583 nm	50 uL (100 ug)	20337-50uL	CF@568 Features
0.5 mL (1 mg)	20337-500uL			
1 mg (lyophilized)	20337-1mg			
CF@583R	585/609 nm	50 uL (100 ug)	20889-50uL	CF@583R Features
0.5 mL (1 mg)	20889-500uL			
CF@594	593/614 nm	50 uL (100 ug)	20226-1	CF@594 Features
0.5 mL (1 mg)	20226			
1 mg (lyophilized)	20226-1mg			
CF@633	630/650 nm	50 uL (100 ug)	20227-1	CF@633 Features
0.5 mL (1 mg)	20227			
1 mg (lyophilized)	20227-1mg			
HRP	N/A	100 uL (100 ug)	20839-100uL	
1 mL (1 mg)	20839-1mL			
1 mg (lyophilized)	20839-1mg			

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 04:05:50 PM. Visit product page to check for updated information before use.

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