



Glowing products for science

Goat Anti-Chicken IgY (H+L), Highly Cross-Adsorbed

Highly cross-adsorbed goat anti-chicken IgY (H L) secondary antibody labeled with our superior CF® Dyes.



Product Description

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

- Highly cross-adsorbed for specific staining with minimal background
- Available in HRP and 11 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)

Product attributes

Antibody type	Secondary
Clonality	Polyclonal
Host species	Goat
Antibody reactivity (target)	Chicken IgY
Species reactivity	Chicken
Cross adsorption	Bovine, Goat, Guinea pig, Horse, Human, Mouse, Rabbit, Rat, Sheep, Syrian hamster
Concentration	2 mg/mL, 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.

Goat Anti-Chicken IgY (H+L), Highly Cross-Adsorbed

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF@405M	408/452 nm	50 uL (100 ug)	20375-50uL	CF@405M Features
0.5 mL (1 mg)	20375-500uL			
1 mg	20375-1mg			
CF@488A	490/515 nm	50 uL (100 ug)	20020-1	CF@488A Features
0.5 mL (1 mg)	20020			
1 mg	20020-1mg			
CF@543	541/560 nm	50 uL (100 ug)	20311-1	CF@543 Features
0.5 mL (1 mg)	20311			
1 mg	20311-1mg			
CF@555	555/565 nm	50 uL (100 ug)	20034-1	CF@555 Features
0.5 mL (1 mg)	20034			
1 mg	20034-1mg			
CF@568	562/583 nm	50 uL (100 ug)	20104-1	CF@568 Features
0.5 mL (1 mg)	20104			
1 mg	20104-1mg			
CF@583R	585/609 nm	50 uL (100 ug)	20898-50uL	CF@583R Features
0.5 mL (1 mg)	20898-500uL			
CF@594	593/614 nm	50 uL (100 ug)	20114-1	CF@594 Features
0.5 mL (1 mg)	20114			
1 mg	20114-1mg			
CF@633	630/650 nm	50 uL (100 ug)	20126-1	CF@633 Features
0.5 mL (1 mg)	20126			
1 mg	20126-1mg			
CF@640R	642/662 nm	50 uL (100 ug)	20084-1	CF@640R Features
0.5 mL (1 mg)	20084			
1 mg	20084-1mg			
CF@647	650/665 nm	50 uL (100 ug)	20044-1	CF@647 Features
0.5 mL (1 mg)	20044			
1 mg	20044-1mg			
CF@660C	667/685 nm	50 uL (100 ug)	20371-50uL	CF@660C Features
0.5 mL (1 mg)	20371-500uL			
1 mg	20371-1mg			
HRP	N/A	100 uL (100 ug)	20474-100uL	
1 mL (1 mg)	20474-1mL			
1 mg	20474-1mg			

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 curated [CF® Dye references](#).

This datasheet was generated on June 3, 2026 at 03:41:27 PM. Visit product page to check for updated information before use.

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