

Goat Anti-Rat IgG (H+L), Highly Cross-Adsorbed

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



Product Description

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

- Highly cross-adsorbed for specific staining with minimal background
- Available in 14 bright and photostable CF® Dyes
- APC, HRP, and R-PE conjugates also available
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

Note: Conjugates of blue fluorescent dyes like CF®350 and 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. 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](#).

Product attributes

Antibody type	Secondary
Clonality	Polyclonal
Host species	Goat
Antibody reactivity (target)	Rat IgG
Species reactivity	Rat
Cross adsorption	Bovine, Horse, Human, Rabbit
Concentration	2 mg/mL, 1 mg/mL (HRP, AP conjugates), 0.5 mg/mL (R-PE, APC 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, R-PE conjugates: PBS/2 mg/mL BSA/0.05% azide, 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

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF@350	347/448 nm	0.5 mL (1 mg) 1 mg 50 uL (100 ug)	20147 20147-1mg 20147-1	CF@350 Features
CF@405M	408/452 nm	0.5 mL (1 mg) 1 mg 50 uL (100 ug)	20374-50uL 20374-1mg 20374-50uL	CF@405M Features
CF@488A	490/515 nm	0.5 mL (1 mg) 1 mg 50 uL (100 ug)	20023 20023-1mg 20023-1	CF@488A Features
CF@543	541/560 nm	0.5 mL (1 mg) 1 mg 50 uL (100 ug)	20321 20321-1mg 20321-1	CF@543 Features
CF@555	555/565 nm	0.5 mL (1 mg) 1 mg 50 uL (100 ug)	20233 20233-1mg 20233-1	CF@555 Features
CF@568	562/583 nm	0.5 mL (1 mg) 1 mg 50 uL (100 ug)	20096 20096-1mg 20096-1	CF@568 Features
CF@583R	585/609 nm	50 uL (100 ug) 0.5 mL (1 mg)	20906-50uL 20906-500uL	CF@583R Features
CF@594	593/614 nm	0.5 mL (1 mg) 1 mg 50 uL (100 ug)	20155 20155-1mg 20155-1	CF@594 Features
CF@633	630/650 nm	0.5 mL (1 mg) 1 mg 50 uL (100 ug)	20133 20133-1mg 20133-1	CF@633 Features
CF@640R	642/662 nm	0.5 mL (1 mg) 1 mg 50 uL (100 ug)	20088 20088-1mg 20088-1	CF@640R Features
CF@647	650/665 nm	0.5 mL (1 mg) 1 mg 50 uL (100 ug)	20283 20283-1mg 20283-1	CF@647 Features
CF@660C	667/685 nm	0.5 mL (1 mg) 1 mg 50 uL (100 ug)	20370-500uL 20370-1mg 20370-50uL	CF@660C Features
CF@680	681/698 nm	0.25 mL 50 uL (100 ug)	20069 20069-1	CF@680 Features
CF@770	770/797 nm	0.25 mL 50 uL (100 ug)	20383-250uL 20383-50uL	CF@770 Features
R-PE	496, 546, 565/578 nm	200 uL (100 ug) 1 mL (500 ug)	20354-200uL 20354-1mL	
APC	650/660 nm	100 uL (50 ug) 0.5 mL (250 ug)	20413-100uL 20413-500uL	
HRP	N/A	100 uL (100 ug) 1 mL (1 mg) 1 mg	20406-100uL 20406-1mL 20406-1mg	

This datasheet was generated on May 20, 2024 at 05:32:42 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/goat-anti-rat-ighl-highly-cross-adsorbed/>