

Goat Anti-Mouse IgG (H+L)

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



Product Description

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

- Available in 19 bright and photostable CF® dyes
- Alkaline phosphatase, APC, HRP, R-PE and biotin conjugates also available
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

Note: Conjugates of blue fluorescent dyes like CF®350, CF®405S, 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.

Goat Anti-Mouse IgG (H+L)

Conjugation	Ex/Em	Size	Catalog No.
CF®350	347/448 nm	50 uL (100 ug)	20140-1
0.5 mL (1 mg)	20140		
1 mg (lyophilized)	20140-1mg		
CF®405S	404/431 nm	50 uL (100 ug)	20080-1
0.5 mL (1 mg)	20080		
1 mg (lyophilized)	20080-1mg		
CF®405M	408/452 nm	50 uL (100 ug)	20180-1
0.5 mL (1 mg)	20180		
1 mg (lyophilized)	20180-1mg		
CF®405L	395/545 nm	50 uL (100 ug)	20408-50uL
0.5 mL (1 mg)	20408-500uL		
1 mg (lyophilized)	20408-1mg		
CF®488A	490/515 nm	50 uL (100 ug)	20010-1
0.5 mL (1 mg)	20010		
1 mg (lyophilized)	20010-1mg		
CF®514	516/548 nm	50 uL (100 ug)	20386-50uL
0.5 mL (1 mg)	20386-500uL		
1 mg (lyophilized)	20386-1mg		
CF®532	527/558 nm	50 uL (100 ug)	20365-50uL
0.5 mL (1 mg)	20365-500uL		
1 mg (lyophilized)	20365-1mg		
CF®543	541/560 nm	50 uL (100 ug)	20306-1
0.5 mL (1 mg)	20306		
1 mg (lyophilized)	20306-1mg		
CF®555	555/565 nm	50 uL (100 ug)	20030-1
0.5 mL (1 mg)	20030		
1 mg (lyophilized)	20030-1mg		
CF®568	562/583 nm	50 uL (100 ug)	20100-1
0.5 mL (1 mg)	20100		
1 mg (lyophilized)	20100-1mg		
CF®594	593/614 nm	50 uL (100 ug)	20110-1
0.5 mL (1 mg)	20110		
1 mg (lyophilized)	20110-1mg		
CF®633	630/650 nm	50 uL (100 ug)	20120-1
0.5 mL (1 mg)	20120		
1 mg (lyophilized)	20120-1mg		
CF®640R	642/662 nm	50 uL (100 ug)	20197-1
0.5 mL (1 mg)	20197		
1 mg (lyophilized)	20197-1mg		
CF®647	650/665 nm	50 uL (100 ug)	20040-1
0.5 mL (1 mg)	20040		
1 mg (lyophilized)	20040-1mg		
CF®660R	663/682 nm	50 uL (100 ug)	20054-1
0.5 mL (1 mg)	20054		

Product attributes

Antibody type	Secondary
Clonality	Polyclonal
Host species	Goat
Antibody reactivity (target)	Mouse IgG
Species reactivity	Mouse
Cross adsorption	Not cross-adsorbed
Concentration	2 mg/mL, 1 mg/mL (HRP, AP conjugates), 0.5 mg/mL (R-PE, APC conjugates)
Antibody/conjugate formulation	AP conjugates: 1 mg/mL in PBS/40% glycerol/5 mg/mL BSA/0.05% azide, 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
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

Conjugation	Ex/Em	Size	Catalog No.
1 mg (lyophilized)	20054-1mg		
CF@660C	667/685 nm	50 uL (100 ug)	20050-1
0.5 mL (1 mg)	20050		
1 mg (lyophilized)	20050-1mg		
CF@740	742/767 nm	50 uL (100 ug)	20984-50uL
0.5 mL (1 mg)	20984-500uL		
CF@750	755/777 nm	50 uL (100 ug)	20070-1
0.5 mL (1 mg)	20070		
1 mg (lyophilized)	20070-1mg		
CF@790	784/806 nm	50 uL (100 ug)	20378-50uL
R-PE	496, 546, 565/578 nm	200 uL (100 ug)	20352-200uL
1 mL (500 ug)	20352-1mL		
APC	650/660 nm	100 uL (50 ug)	20411-100uL
0.5 mL (250 ug)	20411-0.5mL		
Biotin	N/A	50 uL (100 ug)	20183-1
0.5 mL (1 mg)	20183		
1 mg (lyophilized)	20183-1mg		
HRP	N/A	100 uL (100 ug)	20400-100uL
1 mL (1 mg)	20400-1mL		
1 mg (lyophilized)	20400-1mg		
AP	N/A	100 uL (100 ug)	20464-100uL
1 mL (1 mg)	20464-1mL		

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 03:58:51 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/goat-anti-mouse-igg-h/>