

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 18 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. 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)	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	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.
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
		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@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
APC	650/660 nm	1 mL (500 ug)	20352-1mL
		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

This datasheet was generated on July 26, 2024 at 09:33:49 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/goat-anti-mouse-igg-hi/>