

Goat Anti-Rabbit IgG (H+L)

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



Product Description

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

- 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)	Rabbit IgG
Species reactivity	Rabbit
Cross adsorption	Not cross-adsorbed
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

Conjugation	Ex/Em	Size	Catalog No.
CF@350	347/448 nm	50 uL (100 ug)	20141-1
		0.5 mL (1 mg)	20141
		1 mg (lyophilized)	20141-1mg
CF@405S	404/431 nm	50 uL (100 ug)	20082-1
		0.5 mL (1 mg)	20082
		1 mg (lyophilized)	20082-1mg
CF@405M	408/452 nm	50 uL (100 ug)	20181-1
		0.5 mL (1 mg)	20181
		1 mg (lyophilized)	20181-1mg
CF@405L	395/545 nm	50 uL (100 ug)	20409-50uL
		0.5 mL (1 mg)	20409-500uL
		1 mg (lyophilized)	20409-1mg
CF@488A	490/515 nm	50 uL (100 ug)	20012-1
		0.5 mL (1 mg)	20012
		1 mg (lyophilized)	20012-1mg
CF@514	516/548 nm	50 uL (100 ug)	20387-50uL
		0.5 mL (1 mg)	20387-500uL
		1 mg (lyophilized)	20387-1mg
CF@532	527/558 nm	50 uL (100 ug)	20366-50uL
		0.5 mL (1 mg)	20366-500uL
		1 mg (lyophilized)	20366-1mg
CF@543	541/560 nm	50 uL (100 ug)	20309-1
		0.5 mL (1 mg)	20309
		1 mg (lyophilized)	20309-1mg
CF@555	555/565 nm	50 uL (100 ug)	20033-1
		0.5 mL (1 mg)	20033
		1 mg (lyophilized)	20033-1mg
CF@568	562/583 nm	50 uL (100 ug)	20102-1
		0.5 mL (1 mg)	20102
		1 mg (lyophilized)	20102-1mg
CF@594	593/614 nm	50 uL (100 ug)	20112-1
		0.5 mL (1 mg)	20112
		1 mg (lyophilized)	20112-1mg
CF@633	630/650 nm	50 uL (100 ug)	20122-1
		0.5 mL (1 mg)	20122
		1 mg (lyophilized)	20122-1mg
CF@640R	642/662 nm	50 uL (100 ug)	20202-1
		0.5 mL (1 mg)	20202
		1 mg (lyophilized)	20202-1mg
CF@647	650/665 nm	50 uL (100 ug)	20043-1
		0.5 mL (1 mg)	20043
		1 mg (lyophilized)	20043-1mg
CF@660C	667/685 nm	50 uL (100 ug)	20053-1
		0.5 mL (1 mg)	20053
		1 mg (lyophilized)	20053-1mg
CF@660R	663/682 nm	50 uL (100 ug)	20055-1
		0.5 mL (1 mg)	20055
		1 mg (lyophilized)	20055-1mg
CF@750	755/777 nm	50 uL (100 ug)	20073-1
		0.5 mL (1 mg)	20073
		1 mg (lyophilized)	20073-1mg
CF@790	784/806 nm	50 uL (100 ug)	20379-50uL
R-PE	496, 546, 565/578 nm	1 mL (500 ug)	20353-1mL
APC	650/660 nm	200 uL (100 ug)	20353-200uL
		100 uL (50 ug)	20412-100uL
Biotin	N/A	0.5 mL (250 ug)	20412-500uL
		50 uL (100 ug)	20185
HRP	N/A	0.5 mL (1 mg)	20185-1
		1 mg (lyophilized)	20185-1mg
		100 uL (100 ug)	20402-100uL
AP	N/A	1 mL (1 mg)	20402-1mL
		1 mg (lyophilized)	20402-1mg
AP	N/A	100 uL (100 ug)	20465-100uL
		1 mL (1 mg)	20465-1mL