



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357) Email: btinfo@biotium.com

Rabbit Anti-Goat IgG (H+L)

Rabbit anti-goat IgG (H L) secondary antibody labeled with our superior CF® Dyes.



Product attributes

Antibody type	Secondary
Clonality	Polyclonal
Host species	Rabbit
Antibody reactivity (target)	Goat IgG
Species reactivity	Goat
Cross adsorption	Not cross-adsorbed
Concentration	2 mg/mL
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
Secondary/tag antibody applications	Flow cytometry, IHC, IF (cells or tissue sections), Western blot

Product Description

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

- Available in 10 bright and photostable CF® Dyes
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

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](#).

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF@488A	490/515 nm	50 uL (100 ug)	20021-1	CF@488A Features
		0.5 mL (1 mg)	20021	
		1 mg	20021-1mg	
CF@543	541/560 nm	50 uL (100 ug)	20315-1	CF@543 Features
		0.5 mL (1 mg)	20315	
		1 mg	20315-1mg	
CF@555	555/565 nm	50 uL (100 ug)	20031-1	CF@555 Features
		0.5 mL (1 mg)	20031	
		1 mg	20031-1mg	
CF@568	562/583 nm	50 uL (100 ug)	20107-1	CF@568 Features
		0.5 mL (1 mg)	20107	
		1 mg	20107-1mg	
CF@583R	585/609 nm	50 uL (100 ug)	20910-50uL	CF@583R Features
		0.5 mL (1 mg)	20910-500uL	
		1 mg	20910-1mg	
CF@594	593/614 nm	50 uL (100 ug)	20117-1	CF@594 Features
		0.5 mL (1 mg)	20117	
		1 mg	20117-1mg	
CF@633	630/650 nm	50 uL (100 ug)	20128-1	CF@633 Features
		0.5 mL (1 mg)	20128	
		1 mg	20128-1mg	
CF@640R	642/662 nm	50 uL (100 ug)	20090-1	CF@640R Features
		0.5 mL (1 mg)	20090	
		1 mg	20090-1mg	
CF@647	650/665 nm	50 uL (100 ug)	20049-1	CF@647 Features
		0.5 mL (1 mg)	20049	
		1 mg	20049-1mg	
CF@680	681/698 nm	50 uL (100 ug)	20068-1	CF@680 Features
		0.5 mL (1 mg)	20068	
		1 mg	20068-1mg	