



Glowing products for science

Goat Anti-Rabbit IgG (H+L), F(ab')₂ Fragment

Goat anti-rabbit IgG (H+L), F(ab')₂ fragment secondary antibody labeled with our superior CF® Dyes and other labels



Product Description

This is a goat anti-rabbit IgG (H L) F(ab')₂ fragment secondary antibody labeled with our superior CF® Dyes and other labels.

- Available in 14 bright and photostable CF® Dyes, and HRP
- Biotin, APC, and R-PE conjugates also available
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

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

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

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
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	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.

Goat Anti-Rabbit IgG (H+L), F(ab')₂ Fragment

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF@350	347/448 nm	50 uL (100 ug)	20146-1	CF@350 Features
		0.25 mL (500 ug)	20146	
CF@488A	490/515 nm	50 uL (100 ug)	20013-1	CF@488A Features
		0.25 mL (500 ug)	20013	
CF@505	505/519 nm	50 uL (100 ug)	20915-50uL	
		0.25 mL (500 ug)	20915-250uL	
CF@543	541/560 nm	50 uL (100 ug)	20330-1	CF@543 Features
		0.25 mL (500 ug)	20330	
CF@555	555/565 nm	50 uL (100 ug)	20035-1	CF@555 Features
		0.25 mL (500 ug)	20035	
CF@568	562/583 nm	50 uL (100 ug)	20099-1	CF@568 Features
		0.25 mL (500 ug)	20099	
CF@583R	585/609 nm	50 uL (100 ug)	20913-50uL	CF@583R Features
		0.25 mL (500 ug)	20913-250uL	
CF@594	593/614 nm	50 uL (100 ug)	20153-1	CF@594 Features
		0.25 mL (500 ug)	20153	
CF@633	630/650 nm	50 uL (100 ug)	20131-1	CF@633 Features
		0.25 mL (500 ug)	20131	
CF@640R	642/662 nm	50 uL (100 ug)	20087-1	CF@640R Features
		0.25 mL (500 ug)	20087	
CF@647	650/665 nm	50 uL (100 ug)	20045-1	CF@647 Features
		0.25 mL (500 ug)	20045	
CF@660C	667/685 nm	50 uL (100 ug)	20917-50uL	CF@660C Features
		0.25 mL (500 ug)	20917-250uL	
CF@680	681/698 nm	50 uL (100 ug)	20064-1	CF@680 Features
		0.25 mL (500 ug)	20064	
CF@750	755/779 nm	50 uL (100 ug)	20919-50uL	CF@750 Features
		0.25 mL (500 ug)	20919-250uL	
R-PE	496, 546, 565/578 nm	200 uL (100 ug)	20358-200uL	
		500 uL (250 ug)	20358-500uL	
APC	650/660 nm	100 uL (50 ug)	20415-100uL	
		500 uL (250 ug)	20415-500uL	
Biotin	N/A	50 uL (100 ug)	20188-1	
		0.25 mL (500 ug)	20188	

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 [CF® dye references](#).

This datasheet was generated on December 3, 2024 at 09:34:41 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/goat-anti-rabbit-igg-hlfab2-fragment/>