



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)

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	Flow cytometry, IHC, IF (cells or tissue sections), Western blot
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (<i>Bos taurus</i>), 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 0.25 mL (500 ug)	347/448 nm 20146	50 uL (100 ug)	20146-1	CF@350 Features
CF@488A 0.25 mL (500 ug)	490/515 nm 20013	50 uL (100 ug)	20013-1	CF@488A Features
CF@505 0.25 mL (500 ug)	505/519 nm 20915-250uL	50 uL (100 ug)	20915-50uL	
CF@543 0.25 mL (500 ug)	541/560 nm 20330	50 uL (100 ug)	20330-1	CF@543 Features
CF@555 0.25 mL (500 ug)	555/565 nm 20035	50 uL (100 ug)	20035-1	CF@555 Features
CF@568 0.25 mL (500 ug)	562/583 nm 20099	50 uL (100 ug)	20099-1	CF@568 Features
CF@583R 0.25 mL (500 ug)	585/609 nm 20913-250uL	50 uL (100 ug)	20913-50uL	CF@583R Features
CF@594 0.25 mL (500 ug)	593/614 nm 20153	50 uL (100 ug)	20153-1	CF@594 Features
CF@633 0.25 mL (500 ug)	630/650 nm 20131	50 uL (100 ug)	20131-1	CF@633 Features
CF@640R 0.25 mL (500 ug)	642/662 nm 20087	50 uL (100 ug)	20087-1	CF@640R Features
CF@647 0.25 mL (500 ug)	650/665 nm 20045	50 uL (100 ug)	20045-1	CF@647 Features
CF@660C 0.25 mL (500 ug)	667/685 nm 20917-250uL	50 uL (100 ug)	20917-50uL	CF@660C Features
CF@680 0.25 mL (500 ug)	681/698 nm 20064	50 uL (100 ug)	20064-1	CF@680 Features
CF@750 0.25 mL (500 ug)	755/779 nm 20919-250uL	50 uL (100 ug)	20919-50uL	CF@750 Features
R-PE 500 uL (250 ug)	496, 546, 565/578 nm 20358-500uL	200 uL (100 ug)	20358-200uL	
APC 500 uL (250 ug)	650/660 nm 20415-500uL	100 uL (50 ug)	20415-100uL	
Biotin 0.25 mL (500 ug)	N/A 20188	50 uL (100 ug)	20188-1	

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 June 3, 2026 at 04:06:52 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/goat-anti-rabbit-igg-hlfab2-fragment/>