



Glowing products for science

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

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

Goat anti-mouse 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-mouse 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. 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](#).

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
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

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF@350	347/448 nm	50 uL (100 ug) 0.25 mL (500 ug)	20145-1 20145	CF@350 Features
CF@488A	490/515 nm	50 uL (100 ug) 0.25 mL (500 ug)	20011-1 20011	CF@488A Features
CF@505	505/519 nm	50 uL (100 ug) 0.25 mL (500 ug)	20916-50uL 20916-250uL	
CF@543	541/560 nm	50 uL (100 ug) 0.25 mL (500 ug)	20329-1 20329	CF@543 Features
CF@555	555/565 nm	50 uL (100 ug) 0.25 mL (500 ug)	20032-1 20032	CF@555 Features
CF@568	562/583 nm	50 uL (100 ug) 0.25 mL (500 ug)	20109-1 20109	CF@568 Features
CF@583R	585/609 nm	50 uL (100 ug) 0.25 mL (500 ug)	20914-50uL 20914-250uL	CF@583R Features
CF@594	593/614 nm	50 uL (100 ug) 0.25 mL (500 ug)	20119-1 20119	CF@594 Features
CF@633	630/650 nm	50 uL (100 ug) 0.25 mL (500 ug)	20130-1 20130	CF@633 Features
CF@640R	642/662 nm	50 uL (100 ug) 0.25 mL (500 ug)	20086-1 20086	CF@640R Features
CF@647	650/665 nm	50 uL (100 ug) 0.25 mL (500 ug)	20042-1 20042	CF@647 Features
CF@660C	667/685 nm	50 uL (100 ug) 0.25 mL (500 ug)	20918-50uL 20918-250uL	CF@660C Features
CF@680	681/698 nm	50 uL (100 ug) 0.25 mL (500 ug)	20063-1 20063	CF@680 Features
CF@750	755/779 nm	50 uL (100 ug) 0.25 mL (500 ug)	20920-50uL 20920-250uL	CF@750 Features
R-PE	496, 546, 565/578 nm	200 ul (100 ug) 500 ul (250 ug)	20357-200uL 20357-500uL	
APC	650/660 nm	100 ul (50 ug) 500 ul (250 ug)	20414-100uL 20414-500uL	
Biotin	N/A	50 uL (100 ug) 0.25 mL (500 ug)	20187-1 20187	

This datasheet was generated on July 26, 2024 at 04:47:38 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/goat-anti-mouse-igg-hlfab2-fragment/>