

Donkey Anti-Guinea Pig IgG (H+L), Highly Cross-Adsorbed

Highly cross-adsorbed donkey anti-guinea pig IgG (H L) secondary antibody labeled with our superior CF® Dyes.



Product Description

This is a highly cross-adsorbed donkey anti-guinea pig IgG (H L) secondary antibody labeled with our bright and photostable CF® Dyes. To minimize cross-reactivity, the antibody has been adsorbed against bovine, chicken, goat, Syrian hamster, horse, human, mouse, rabbit, and sheep serum.

- Highly cross-adsorbed for specific staining with minimal background
- Available in 12 bright and photostable CF® Dyes
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

Note: Conjugates of blue fluorescent dyes like 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 [CF® dye references](#).

Product attributes

Antibody type	Secondary
Clonality	Polyclonal
Host species	Donkey
Antibody reactivity (target)	Guinea Pig IgG
Species reactivity	Guinea pig
Cross adsorption	Bovine, Chicken, Goat, Horse, Human, Mouse, Rabbit, Rat, Sheep, Syrian hamster
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

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF®405S	404/431 nm	50 uL (100 ug)	20356-50uL	CF®405S Features
		0.5 mL (1 mg)	20356-500uL	
		1 mg	20356-1mg	
CF®405M	408/452 nm	50 uL (100 ug)	20376-50uL	CF®405M Features
		0.5 mL (1 mg)	20376-500uL	
		1 mg	20376-1mg	
CF®488A	490/515 nm	50 uL (100 ug)	20169-1	CF®488A Features
		0.5 mL (1 mg)	20169	
		1 mg	20169-1mg	
CF®543	541/560 nm	50 uL (100 ug)	20316-1	CF®543 Features
		0.5 mL (1 mg)	20316	
		1 mg	20316-1mg	
CF®555	555/565 nm	50 uL (100 ug)	20276-1	CF®555 Features
		0.5 mL (1 mg)	20276	
		1 mg	20276-1mg	
CF®568	562/583 nm	50 uL (100 ug)	20377-50uL	CF®568 Features
		0.5 mL (1 mg)	20377-500uL	
		1 mg	20377-1mg	
CF®583R	585/609 nm	50 uL (100 ug)	20893-50uL	CF®583R Features
		0.5 mL (1 mg)	20893-500uL	
		1 mg	20893-1mg	
CF®594	593/614 nm	50 uL (100 ug)	20170-1	CF®594 Features
		0.5 mL (1 mg)	20170	
		1 mg	20170-1mg	
CF®633	630/650 nm	50 uL (100 ug)	20171-1	CF®633 Features
		0.5 mL (1 mg)	20171	
		1 mg	20171-1mg	
CF®660C	667/685 nm	50 uL (100 ug)	20372-50uL	CF®660C Features
		0.5 mL (1 mg)	20372-500uL	
		1 mg	20372-1mg	
CF®680	681/698 nm	50 uL (100 ug)	20241-1	CF®680 Features
		0.25 mL (500 ug)	20241	
		0.25 mL (500 ug)	20241-1	
CF®770	770/797 nm	50 uL (100 ug)	20242-1	CF®770 Features
		0.25 mL (500 ug)	20242	
		0.25 mL (500 ug)	20242	

This datasheet was generated on April 28, 2024 at 12:01:32 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/donkey-anti-guinea-pig-igg-hl-highly-cross-adsorbed/>