

## Donkey Anti-Rabbit IgG (H+L), Highly Cross-Adsorbed

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



### Product Description

This is a highly cross-adsorbed donkey anti-rabbit 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, guinea pig, Syrian hamster, horse, human, mouse, rat, and sheep serum.

- Highly cross-adsorbed for specific staining with minimal background
- Available with 20 bright and photostable CF® Dyes
- Alkaline phosphatase, biotin, and HRP 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.

### Product attributes

<b>Antibody type</b>	Secondary
<b>Clonality</b>	Polyclonal
<b>Host species</b>	Donkey
<b>Antibody reactivity (target)</b>	Rabbit IgG
<b>Species reactivity</b>	Rabbit
<b>Cross adsorption</b>	Bovine, Chicken, Goat, Guinea pig, Horse, Human, Mouse, Rat, Sheep, Syrian hamster
<b>Concentration</b>	2 mg/mL, 1 mg/mL (HRP, AP conjugates)
<b>Antibody/conjugate formulation</b>	AP conjugates: 1 mg/mL in PBS/40% glycerol/5 mg/mL BSA/0.05% azide, Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide, Lyophilized: PBS/15 mg/mL BSA/20 mg/mL trehalose after reconstitution, HRP conjugates: PBS/50% glycerol/15 mg/mL BSA, HRP conjugates (lyophilized): PBS/10 mg/mL trehalose after reconstitution
<b>Secondary/tag antibody applications</b>	ELISA, Flow cytometry, IHC, IF (cells or tissue sections), Western blot
<b>Product origin</b>	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.

## Donkey Anti-Rabbit IgG (H+L), Highly Cross-Adsorbed

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF®350	347/448 nm	50 uL (100 ug)	<a href="#">20351-1</a>	<a href="#">CF®350 Features</a>
0.5 mL (1 mg)	<a href="#">20351</a>			
1 mg	<a href="#">20351-1mg</a>			
CF®405S	404/431 nm	50 uL (100 ug)	<a href="#">20420-50uL</a>	<a href="#">CF®405S Features</a>
0.5 mL (1 mg)	<a href="#">20420-500uL</a>			
1 mg	<a href="#">20420-1mg</a>			
CF®430	426/498 nm	50 uL (100 ug)	<a href="#">20462-50uL</a>	<a href="#">CF®430 Features</a>
0.5 mL (1 mg)	<a href="#">20462-500uL</a>			
1 mg	<a href="#">20462-1mg</a>			
CF®488A	490/515 nm	50 uL (100 ug)	<a href="#">20015-1</a>	<a href="#">CF®488A Features</a>
0.5 mL (1 mg)	<a href="#">20015</a>			
1 mg	<a href="#">20015-1mg</a>			
CF®543	541/560 nm	50 uL (100 ug)	<a href="#">20308-1</a>	<a href="#">CF®543 Features</a>
0.5 mL (1 mg)	<a href="#">20308</a>			
1 mg	<a href="#">20308-1mg</a>			
CF®555	555/565 nm	50 uL (100 ug)	<a href="#">20038-1</a>	<a href="#">CF®555 Features</a>
0.5 mL (1 mg)	<a href="#">20038</a>			
1 mg	<a href="#">20038-1mg</a>			
CF®568	562/583 nm	50 uL (100 ug)	<a href="#">20098-1</a>	<a href="#">CF®568 Features</a>
0.5 mL (1 mg)	<a href="#">20098</a>			
1 mg	<a href="#">20098-1mg</a>			
CF®583R	585/609 nm	50 uL (100 ug)	<a href="#">20895-50uL</a>	<a href="#">CF®583R Features</a>
0.5 mL (1 mg)	<a href="#">20895-500uL</a>			
CF®594	593/614 nm	50 uL (100 ug)	<a href="#">20152-1</a>	<a href="#">CF®594 Features</a>
0.5 mL (1 mg)	<a href="#">20152</a>			
1 mg	<a href="#">20152-1mg</a>			
CF®633	630/650 nm	50 uL (100 ug)	<a href="#">20125-1</a>	<a href="#">CF®633 Features</a>
0.5 mL (1 mg)	<a href="#">20125</a>			
1 mg	<a href="#">20125-1mg</a>			
CF®640R	642/662 nm	50 uL (100 ug)	<a href="#">20178-1</a>	<a href="#">CF®640R Features</a>
0.5 mL (1 mg)	<a href="#">20178</a>			
1 mg	<a href="#">20178-1mg</a>			
CF®647	650/665 nm	50 uL (100 ug)	<a href="#">20047-1</a>	<a href="#">CF®647 Features</a>
0.5 mL (1 mg)	<a href="#">20047</a>			
1 mg	<a href="#">20047-1mg</a>			
CF®660R	663/682 nm	50 uL (100 ug)	<a href="#">20389-50uL</a>	<a href="#">CF®660R Features</a>
0.5 mL (1 mg)	<a href="#">20389-500uL</a>			
1 mg	<a href="#">20389-1mg</a>			
CF®680	681/698 nm	50 uL (100 ug)	<a href="#">20418-50uL</a>	<a href="#">CF®680 Features</a>
0.25 mL (500 ug)	<a href="#">20418-250uL</a>			
CF®680R	680/701 nm	50 uL (100 ug)	<a href="#">20195-1</a>	<a href="#">CF®680R Features</a>

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
0.25 mL (500 ug)	<a href="#">20195</a>			
CF@725	729/750 nm	50 uL (100 ug)	<a href="#">20505-50uL</a>	<a href="#">CF@725 Features</a>
0.25 mL (500 ug)	<a href="#">20505-250uL</a>			
CF@740	742/767 nm	50 uL (100 ug)	<a href="#">20986-50uL</a>	<a href="#">CF@740 Features</a>
0.25 mL (500 ug)	<a href="#">20986-250uL</a>			
CF@750	755/777 nm	50 uL (100 ug)	<a href="#">20298-1</a>	<a href="#">CF@750 Features</a>
0.25 mL (500 ug)	<a href="#">20298</a>			
CF@770	770/797 nm	50 uL (100 ug)	<a href="#">20484-50uL</a>	<a href="#">CF@770 Features</a>
0.25 mL (500 ug)	<a href="#">20484-250uL</a>			
CF@790	784/806 nm	50 uL (100 ug)	<a href="#">20344-50uL</a>	<a href="#">CF@790 Features</a>
CF@800	797/816 nm	50 uL (100 ug)	<a href="#">20833-50uL</a>	<a href="#">CF@800 Features</a>
Biotin	N/A	50 uL (100 ug)	<a href="#">20191-1</a>	
0.5 mL (1 mg)	<a href="#">20191</a>			
1 mg	<a href="#">20191-1mg</a>			
HRP	N/A	100 uL (100 ug)	<a href="#">20405-100uL</a>	
1 mL (1 mg)	<a href="#">20405-1mL</a>			
1 mg	<a href="#">20405-1mg</a>			
AP	N/A	100 uL (100 ug)	<a href="#">20467-100uL</a>	
1 mL (1 mg)	<a href="#">20467-1mL</a>			

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 dH<sub>2</sub>O

HRP or DNP conjugates: add 1 mL dH<sub>2</sub>O

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](#).

This datasheet was generated on June 3, 2026 at 03:15:10 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/donkey-anti-rabbit-igg-hl-highly-cross-adsorbed/>