

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

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



Product Description

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

- Highly cross-adsorbed for specific staining with minimal background
- Available in 6 bright and photostable CF® Dyes
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

Product attributes

Antibody type	Secondary, Anti-Human Immunoglobulin
Clonality	Polyclonal
Host species	Donkey
Antibody reactivity (target)	Human IgG
Species reactivity	Human
Cross adsorption	Bovine, Chicken, Goat, Guinea pig, Horse, 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
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.

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

Conjugation	Ex/Em	Size	Catalog No.
CF®488A	490/515 nm	50 uL (100 ug)	20074-1
		0.5 mL (1 mg)	20074
		1 mg	20074-1mg
CF®543	541/560 nm	50 uL (100 ug)	20318-1
		0.5 mL (1 mg)	20318
		1 mg	20318-1mg
CF®594	593/614 nm	50 uL (100 ug)	20075-1
		0.5 mL (1 mg)	20075
		1 mg	20075-1mg
CF®633	630/650 nm	50 uL (100 ug)	20076-1
		0.5 mL (1 mg)	20076
		1 mg	20076-1mg
CF®680	681/698 nm	50 uL (100 ug)	20278-1
		0.25 mL (500 ug)	20278
CF®770	770/797 nm	50 uL (100 ug)	20279-1
		0.25 mL (500 ug)	20279

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 April 16, 2025 at 09:29:31 AM. Visit product page to check for updated information before use.

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