Glowing products for science

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

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

Biotum

Product Description

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

- Highly cross-adsorbed for specific staining with minimal background
- Available in 11 bright and photostable CF® Dyes
- 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 Email: btinfo@biotium.com

Product attributes

Antibody type	Secondary		
Clonality	Polyclonal		
Host species	Donkey		
Antibody reactivity (target)	Sheep IgG		
Species reactivity	Sheep		
Cross adsorption	Chicken, Guinea pig, Horse, Human, Mouse, Rabbit, Rat, 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 (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.		

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

0.5 mL (1 mg)	Conjugation	Ex/Em	Size	Catalog No.	Dye Features
T mg	CF®350	347/448 nm	50 uL (100 ug)	<u>20148-1</u>	CF®350 Features
Fig. 488A 490/515 nm			0.5 mL (1 mg)	<u>20148</u>	
0.5 mL (1 mg) 20024 1 mg 20024-1 mg 20024-1 mg 20024-1 mg 20024-1 mg 20024-1 mg 20024-1 mg 20022-1 mg 20322-1 CF®543 Features 20.5 mL (100 ug) 20322-1 mg 20322-1 mg 20322-1 mg 20322-1 mg 20322-1 mg 20322-1 mg 20234-1 CF®555 Features 20.5 mL (1 mg) 20234 1 mg 20234-1 mg 20235-1			1 mg	20148-1mg	
Time	CF®488A	490/515 nm	50 uL (100 ug)	<u>20024-1</u>	CF®488A Features
Feb Section Feb Section Sould (100 ug) 20322-1 CF6543 Features CF6555 Section			0.5 mL (1 mg)	<u>20024</u>	
0.5 mL (1 mg) 20322 1 mg 20322-1mg 20322-1mg 20322-1mg 20322-1mg 20322-1mg 20322-1mg 20324-1 CF®555 Features 20.5 mL (1 mg) 20234 1 mg 20234-1mg 20235-1mg 20235-1mg			1 mg	20024-1mg	
1 mg 20322-1mg FB555 555/565 nm 50 uL (100 ug) 20234-1 CFB555 Features 0.5 mL (1 mg) 20234 1 mg 20234-1mg FB568 562/583 nm 50 uL (100 ug) 20095-1 CFB568 Features 0.5 mL (1 mg) 20095 1 mg 20095-1mg FB583R 585/609 nm 50 uL (100 ug) 20897-50uL CFB583R Features 0.5 mL (1 mg) 20897-50uL FB594 593/614 nm 50 uL (100 ug) 20156-1 CFB594 Features 0.5 mL (1 mg) 20156-1 CFB594 Features 0.5 mL (1 mg) 20156-1 CFB633 Features 0.5 mL (1 mg) 20156-1 CFB633 Features 0.5 mL (1 mg) 20134-1 CFB640R Features 0.5 mL (1 mg) 20083-1 CFB640R Features 0.5 mL (1 mg) 20083-1 CFB640R Features 0.5 mL (1 mg) 20083-1 CFB647 Features	CF®543	541/560 nm	50 uL (100 ug)	<u>20322-1</u>	CF®543 Features
Free			0.5 mL (1 mg)	<u>20322</u>	
D.5 mL (1 mg) 20234 1 mg 20234-1mg 20233-1mg 20233-1mg 20233-1mg 20234-1mg 20233-1mg 20234-1mg 20234-1mg			1 mg	20322-1mg	
1 mg 20234-1mg F®568 562/583 nm 50 uL (100 ug) 20095-1 CF®568 Features 0.5 mL (1 mg) 20095 1 mg 20095-1mg F®583R 585/609 nm 50 uL (100 ug) 20897-50uL 0.5 mL (1 mg) 20897-50uL 0.5 mL (1 mg) 20897-500uL F®594 593/614 nm 50 uL (100 ug) 20156-1 CF®594 Features 0.5 mL (1 mg) 20156 1 mg 20156-1mg F®633 630/650 nm 50 uL (100 ug) 20134-1 CF®633 Features 0.5 mL (1 mg) 20134-1 1 mg 20134-1mg F®640R 642/662 nm 50 uL (100 ug) 20083-1 CF®640R Features 0.5 mL (1 mg) 20083-1mg F®647 650/665 nm 50 uL (100 ug) 20284-1 CF®647 Features	CF®555	555/565 nm	50 uL (100 ug)	<u>20234-1</u>	CF®555 Features
F8568 562/583 nm 50 uL (100 ug) 20095-1 CF8568 Features 50 uL (100 ug) 20095 1 mg 20095-1mg F8583R 585/609 nm 50 uL (100 ug) 20897-50uL 50 uL (100 ug) 20897-50uL 50 uL (100 ug) 20156-1 50 uL (100 ug) 20156-1 1 mg 20156-1mg F8633 630/650 nm 50 uL (100 ug) 20156-1mg F8663 630/650 nm 50 uL (100 ug) 20134-1 1 mg 20134-1mg F86640R 642/662 nm 50 uL (100 ug) 20083-1 1 mg 20083-1mg F86647 650/665 nm 50 uL (100 ug) 20284-1 CF86647 Features			0.5 mL (1 mg)	<u>20234</u>	
0.5 mL (1 mg) 20095 1 mg 20095-1mg 2F®583R 585/609 nm 50 uL (100 ug) 20897-50uL 0.5 mL (1 mg) 20897-50uL 0.5 mL (1 mg) 20156-1 0.5 mL (1 mg) 20156-1mg 0.5 mL (1 mg) 20134-1 0.5 mL (1 mg) 20134-1 0.5 mL (1 mg) 20134-1mg 0.5 mL (1 mg) 20134-1mg 0.5 mL (1 mg) 20083-1 0.5 mL (1 mg) 20083 1 mg 20083-1mg 0.5 mG647 0.5 mL (100 ug) 20083-1 005 0.5 mL (1 mg) 20083-1 0.5 mL (1 mg)			1 mg	20234-1mg	
1 mg 20095-1mg F®583R 585/609 nm 50 uL (100 ug) 20897-50uL CF®583R Features 0.5 mL (1 mg) 20897-50uL CF®594 593/614 nm 50 uL (100 ug) 20156-1 CF®594 Features 0.5 mL (1 mg) 20156 1 mg 20156-1mg CF®633 630/650 nm 50 uL (100 ug) 20134-1 CF®633 Features 0.5 mL (1 mg) 20134 1 mg 20134-1mg CF®640R 642/662 nm 50 uL (100 ug) 20083-1 CF®640R Features 0.5 mL (1 mg) 20083 1 mg 20083-1mg CF®647 650/665 nm 50 uL (100 ug) 20284-1 CF®647 Features	CF®568	562/583 nm	50 uL (100 ug)	<u>20095-1</u>	CF®568 Features
F®583R 585/609 nm 50 uL (100 ug) 20897-50uL CF®583R Features 0.5 mL (1 mg) 20897-500uL F®594 593/614 nm 50 uL (100 ug) 20156-1 CF®594 Features 0.5 mL (1 mg) 20156 1 mg 20156-1mg F®633 630/650 nm 50 uL (100 ug) 20134-1 CF®633 Features 0.5 mL (1 mg) 20134 1 mg 20134-1mg F®640R 642/662 nm 50 uL (100 ug) 20083-1 CF®640R Features 0.5 mL (1 mg) 20083 1 mg 20083-1mg F®647 650/665 nm 50 uL (100 ug) 20284-1 CF®647 Features			0.5 mL (1 mg)	<u>20095</u>	
0.5 mL (1 mg) 20897-500uL 50 uL (100 ug) 20156-1 CF®594 Features 0.5 mL (1 mg) 20156 1 mg 20156-1mg F®633 630/650 nm 50 uL (100 ug) 20134-1 CF®633 Features 0.5 mL (1 mg) 20134-1mg F®640R 642/662 nm 50 uL (100 ug) 20083-1 CF®640R Features 0.5 mL (1 mg) 20083 1 mg 20083-1mg F®647 650/665 nm 50 uL (100 ug) 20284-1 CF®647 Features			1 mg	20095-1mg	
F®594 593/614 nm 50 uL (100 ug) 20156-1 CF®594 Features 0.5 mL (1 mg) 20156 1 mg 20156-1mg F®633 630/650 nm 50 uL (100 ug) 20134-1 CF®633 Features 0.5 mL (1 mg) 20134 1 mg 20134-1mg F®640R 642/662 nm 50 uL (100 ug) 20083-1 CF®640R Features 0.5 mL (1 mg) 20083 1 mg 20083-1mg F®647 650/665 nm 50 uL (100 ug) 20284-1 CF®647 Features	CF®583R	585/609 nm	50 uL (100 ug)	20897-50uL	CF®583R Features
0.5 mL (1 mg) 20156 1 mg 20156-1mg 50 uL (100 ug) 20134-1 CF®633 Features 0.5 mL (1 mg) 20134 1 mg 20134-1mg FF®640R 642/662 nm 50 uL (100 ug) 20083-1 CF®640R Features 0.5 mL (1 mg) 20083 1 mg 20083-1mg FF®647 650/665 nm 50 uL (100 ug) 20284-1 CF®647 Features			0.5 mL (1 mg)	20897-500uL	
1 mg 20156-1mg F®633 630/650 nm 50 uL (100 ug) 20134-1 CF®633 Features 0.5 mL (1 mg) 20134-1mg 1 mg 20134-1mg F®640R 642/662 nm 50 uL (100 ug) 20083-1 CF®640R Features 0.5 mL (1 mg) 20083 1 mg 20083-1mg F®647 650/665 nm 50 uL (100 ug) 20284-1 CF®647 Features	CF®594	593/614 nm	50 uL (100 ug)	<u>20156-1</u>	CF®594 Features
F®633 630/650 nm 50 uL (100 ug) 20134-1 CF®633 Features 0.5 mL (1 mg) 20134 1 mg 20134-1mg F®640R 642/662 nm 50 uL (100 ug) 20083-1 CF®640R Features 0.5 mL (1 mg) 20083 1 mg 20083-1mg F®647 650/665 nm 50 uL (100 ug) 20284-1 CF®647 Features			0.5 mL (1 mg)	<u>20156</u>	
0.5 mL (1 mg) 20134 1 mg 20134-1mg 5 mL (100 ug) 20083-1 CF®640R Features 0.5 mL (1 mg) 20083 1 mg 20083-1mg F®647 650/665 nm 50 uL (100 ug) 20284-1 CF®647 Features			1 mg	20156-1mg	
1 mg 20134-1mg F®640R 642/662 nm 50 uL (100 ug) 20083-1 CF®640R Features 0.5 mL (1 mg) 20083 1 mg 20083-1mg F®647 650/665 nm 50 uL (100 ug) 20284-1 CF®647 Features	CF®633	630/650 nm	50 uL (100 ug)	<u>20134-1</u>	CF®633 Features
F®640R 642/662 nm 50 uL (100 ug) 20083-1 CF®640R Features 0.5 mL (1 mg) 20083 1 mg 20083-1 mg F®647 650/665 nm 50 uL (100 ug) 20284-1 CF®647 Features			0.5 mL (1 mg)	<u>20134</u>	
0.5 mL (1 mg) 20083 1 mg 20083-1mg CF®647 650/665 nm 50 uL (100 ug) 20284-1 CF®647 Features			1 mg	20134-1mg	
1 mg 20083-1mg CF®647 650/665 nm 50 uL (100 ug) 20284-1 CF®647 Features	CF®640R	642/662 nm	50 uL (100 ug)	<u>20083-1</u>	CF®640R Features
F®647 650/665 nm 50 uL (100 ug) <u>20284-1</u> <u>CF®647 Features</u>			0.5 mL (1 mg)	<u>20083</u>	
			1 mg	20083-1mg	
	CF®647	650/665 nm	50 uL (100 ug)	20284-1	CF®647 Features
0.5 mL (1 mg) <u>20284</u>			0.5 mL (1 mg)	<u>20284</u>	
1 mg <u>20284-1mg</u>			1 mg	20284-1mg	
F®680 681/698 nm 50 uL (100 ug) <u>20062-1</u> <u>CF®680 Features</u>	CF®680	681/698 nm	50 uL (100 ug)	<u>20062-1</u>	CF®680 Features
0.25 mL (500 ug) <u>20062</u>			0.25 mL (500 ug)	20062	

View our full selection of <u>Secondary Antibodies</u>, or search our catalog using our <u>Antibody Finder</u>. Alternatively, you can view our <u>secondary antibody product listings</u> 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 curated CF® Dye references.

This datasheet was generated on November 6, 2025 at 08:57:05 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/donkey-anti-sheep-igghl-highly-cross-adsorbed/