

Goat Anti-Llama IgG (H+L)

Goat anti-llama IgG (H L) secondary antibody labeled with our superior CF® Dyes.



Product Description

This is a goat anti-llama IgG (H L) secondary antibody labeled with our superior CF® Dyes, HRP, or DNP.

- Available in 14 bright and photostable CF® Dyes, HRP, or DNP
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

Note: Conjugates of blue fluorescent dyes like CF@405S 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

Clonality	Polyclonal
Antibody type	Secondary
Host species	Goat
Antibody reactivity (target)	Llama IgG
Species reactivity	Llama
Cross adsorption	Not cross-adsorbed
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, HRP conjugates: PBS/50% glycerol/15 mg/mL BSA, HRP conjugates (lyophilized): PBS/10 mg/mL trehalose after reconstitution
Concentration	1 mg/mL (HRP conjugate), 2 mg/mL
Secondary/tag antibody applications	ELISA, 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.

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF@405S 0.5 mL (1 mg) 1 mg (lyophilized)	404/431 nm 20844-500uL 20844-1mg	50 uL (100 ug)	20844-50uL	CF@405S Features
CF@488A 0.5 mL (1 mg) 1 mg (lyophilized)	490/515 nm 20845-500uL 20845-1mg	50 uL (100 ug)	20845-50uL	CF@488A Features
CF@543 0.5 mL (1 mg) 1 mg (lyophilized)	541/560 nm 20846-500uL 20846-1mg	50 uL (100 ug)	20846-50uL	CF@543 Features
CF@555 0.5 mL (1 mg) 1 mg (lyophilized)	555/565 nm 20847-500uL 20847-1mg	50 uL (100 ug)	20847-50uL	CF@555 Features
CF@568 0.5 mL (1 mg) 1 mg (lyophilized)	562/583 nm 20848-500uL 20848-1mg	50 uL (100 ug)	20848-50uL	CF@568 Features
CF@583R 0.5 mL (1 mg)	585/609 nm 20902-500uL	50 uL (100 ug)	20902-50uL	CF@583R Features
CF@594 0.5 mL (1 mg) 1 mg (lyophilized)	593/614 nm 20849-500uL 20849-1mg	50 uL (100 ug)	20849-50uL	CF@594 Features
CF@640R 0.5 mL (1 mg) 1 mg (lyophilized)	642/662 nm 20850-500uL 20850-1mg	50 uL (100 ug)	20850-50uL	CF@640R Features
CF@647 0.5 mL (1 mg) 1 mg (lyophilized)	650/665 nm 20851-500uL 20851-1mg	50 uL (100 ug)	20851-50uL	CF@647 Features
CF@660R 0.5 mL (1 mg) 1 mg (lyophilized)	663/682 nm 20853-500uL 20853-1mg	50 uL (100 ug)	20853-50uL	CF@660R Features
CF@660C 0.5 mL (1 mg) 1 mg (lyophilized)	667/685 nm 20852-500uL 20852-1mg	50 uL (100 ug)	20852-50uL	CF@660C Features
CF@680 0.5 mL (1 mg) 1 mg (lyophilized)	681/698 nm 20854-500uL 20854-1mg	50 uL (100 ug)	20854-50uL	CF@680 Features
CF@680R 0.5 mL (1 mg) 1 mg (lyophilized)	680/701 nm 20855-500uL 20855-1mg	50 uL (100 ug)	20855-50uL	CF@680R Features
CF@750 0.5 mL (1 mg) 1 mg (lyophilized)	755/777 nm 20856-500uL 20856-1mg	50 uL (100 ug)	20856-50uL	CF@750 Features
DNP 0.5 mL (1 mg) 1 mg (lyophilized)	N/A 20983-500uL 20983-1mg	50 uL (100 ug)	20983-50uL	
HRP 1 mL (1 mg)	N/A 20475-1mL	100 uL (100 ug)	20475-100uL	

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
1 mg (lyophilized)	20475-1mg			

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 June 3, 2026 at 03:29:29 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/goat-anti-lama-igg-hl/>