

## Llama Anti-Rabbit IgG (H+L)

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



### Product Description

This is a llama anti-rabbit IgG (H L) secondary antibody labeled with our superior CF® Dyes.

- Available in 6 bright and photostable CF® Dyes
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

### Product attributes

<b>Clonality</b>	Polyclonal
<b>Cross adsorption</b>	Not cross-adsorbed
<b>Antibody type</b>	Secondary
<b>Concentration</b>	2 mg/mL
<b>Antibody/conjugate formulation</b>	Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide, Lyophilized: PBS/15 mg/mL BSA/20 mg/mL trehalose after reconstitution
<b>Species reactivity</b>	Rabbit
<b>Secondary/tag antibody applications</b>	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.
<b>Host species</b>	Llama
<b>Antibody reactivity (target)</b>	Rabbit IgG, Rat IgG

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
<a href="#">CF@488A</a>	490/515 nm	50 uL (100 ug)	<a href="#">20449-50uL</a>	<a href="#">CF@488A Features</a>
0.5 mL (1 mg)	<a href="#">20449-500uL</a>			
1 mg (lyophilized)	<a href="#">20449-1mg</a>			
<a href="#">CF@568</a>	562/583 nm	50 uL (100 ug)	<a href="#">20450-50uL</a>	<a href="#">CF@568 Features</a>
0.5 mL (1 mg)	<a href="#">20450-500uL</a>			
1 mg (lyophilized)	<a href="#">20450-1mg</a>			
<a href="#">CF@583R</a>	585/609 nm	50 uL (100 ug)	<a href="#">20909-50uL</a>	<a href="#">CF@583R Features</a>
0.5 mL (1 mg)	<a href="#">20909-500uL</a>			
<a href="#">CF@594</a>	593/614 nm	50 uL (100 ug)	<a href="#">20451-50uL</a>	<a href="#">CF@594 Features</a>
0.5 mL (1 mg)	<a href="#">20451-500uL</a>			
1 mg (lyophilized)	<a href="#">20451-1mg</a>			
<a href="#">CF@640R</a>	642/662 nm	50 uL (100 ug)	<a href="#">20452-50uL</a>	<a href="#">CF@640R Features</a>
0.5 mL (1 mg)	<a href="#">20452-500uL</a>			
1 mg (lyophilized)	<a href="#">20452-1mg</a>			
<a href="#">CF@647</a>	650/665 nm	50 uL (100 ug)	<a href="#">20453-50uL</a>	<a href="#">CF@647 Features</a>
0.5 mL (1 mg)	<a href="#">20453-500uL</a>			
1 mg (lyophilized)	<a href="#">20453-1mg</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 [CF® dye references](#).

This datasheet was generated on June 3, 2026 at 04:38:19 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/llama-anti-rabbit-igg-hl/>