

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

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](#).

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

Product attributes	
Clonality	Polyclonal
Cross adsorption	Not cross-adsorbed
Antibody type	Secondary
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
Species reactivity	Rabbit
Secondary/tag antibody applications	Flow cytometry, IHC, IF (cells or tissue sections), Western blot
Host species	Llama
Antibody reactivity (target)	Rabbit IgG, Rat IgG