

Llama Anti-Mouse IgG (H+L)

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



Product Description

This is a llama anti-mouse 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.

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	Mouse
Secondary/tag antibody applications	Flow cytometry, IHC, IF (cells or tissue sections), Western blot
Host species	Llama
Antibody reactivity (target)	Mouse IgG

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)	20454-50uL	CF@488A Features
		0.5 mL (1 mg)	20454-500uL	
		1 mg (lyophilized)	20454-1mg	
CF@568	562/583 nm	50 uL (100 ug)	20455-50uL	CF@568 Features
		0.5 mL (1 mg)	20455-500uL	
		1 mg (lyophilized)	20455-1mg	
CF@583R	585/609 nm	50 uL (100 ug)	20908-50uL	CF@583R Features
		0.5 mL (1 mg)	20908-500uL	
		1 mg (lyophilized)	20456-1mg	
CF@594	593/614 nm	50 uL (100 ug)	20456-50uL	CF@594 Features
		0.5 mL (1 mg)	20456-500uL	
		1 mg (lyophilized)	20456-1mg	
CF@640R	642/662 nm	50 uL (100 ug)	20457-50uL	CF@640R Features
		0.5 mL (1 mg)	20457-500uL	
		1 mg (lyophilized)	20457-1mg	
CF@647	650/665 nm	50 uL (100 ug)	20458-50uL	CF@647 Features
		0.5 mL (1 mg)	20458-500uL	
		1 mg (lyophilized)	20458-1mg	