

Chicken Anti-Rabbit IgG (H+L)

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



Product attributes

Antibody type	Secondary
Clonality	Polyclonal
Host species	Chicken
Antibody reactivity (target)	Rabbit IgG
Species reactivity	Rabbit
Cross adsorption	Not cross-adsorbed
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 Description

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

- Available in 6 bright and photostable CF® Dyes, and HRP
- 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. 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 curated [CF® Dye references](#).

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF®350	347/448 nm	50 uL (100 ug)	20332-50uL	CF®350 Features
		0.5 mL (1 mg)	20332-500uL	
		1 mg (lyophilized)	20332-1mg	
CF®488A	490/515 nm	50 uL (100 ug)	20209-1	CF®488A Features
		0.5 mL (1 mg)	20209	
		1 mg (lyophilized)	20209-1mg	
CF®543	541/560 nm	50 uL (100 ug)	20335-50uL	CF®543 Features
		0.5 mL (1 mg)	20335-500uL	
		1 mg (lyophilized)	20335-1mg	
CF®568	562/583 nm	50 uL (100 ug)	20339-50uL	CF®568 Features
		0.5 mL (1 mg)	20339-500uL	
		1 mg (lyophilized)	20339-1mg	
CF®583R	585/609 nm	50 uL (100 ug)	20891-50uL	CF®583R Features
		0.5 mL (1 mg)	20891-500uL	
CF®594	593/614 nm	50 uL (100 ug)	20223-1	CF®594 Features
		0.5 mL (1 mg)	20223	
		1 mg (lyophilized)	20223-1mg	
CF®633	630/650 nm	50 uL (100 ug)	20224-1	CF®633 Features
		0.5 mL (1 mg)	20224	
		1 mg (lyophilized)	20224-1mg	