

Rabbit Anti-Sheep IgG (H+L), Highly Cross-Adsorbed

Highly cross-adsorbed rabbit anti-sheep IgG (H+L) secondary antibody labeled with our superior CF® dyes.

Product Description

This is a highly cross-adsorbed rabbit anti-sheep IgG (H L) secondary antibody labeled with our bright and photostable CF® dyes. To minimize cross-reactivity, the antibody has been adsorbed against human serum.

- Highly cross-adsorbed for specific staining with minimal background
- Available in 4 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

Antibody type	Secondary
Clonality	Polyclonal
Host species	Rabbit
Antibody reactivity (target)	Sheep IgG
Species reactivity	Sheep
Cross adsorption	Human
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

References

Download a list of [CF® dye references](#).

Conjugation	Ex/Em	Size	Catalog No.
CF®488A	490/515 nm	50 uL (100 ug)	20172-1
		0.5 mL (1 mg)	20172
		1 mg	20172-1mg
CF®543	541/560 nm	50 uL (100 ug)	20323-1
		0.5 mL (1 mg)	20323
		1 mg	20323-1mg
CF®594	593/614 nm	50 uL (100 ug)	20173-1
		0.5 mL (1 mg)	20173
		1 mg	20173-1mg
CF®633	630/650 nm	50 uL (100 ug)	20174-1
		0.5 mL (1 mg)	20174
		1 mg	20174-1mg