

## Rabbit Anti-Guinea Pig IgG (H+L)

Rabbit anti-guinea pig IgG (H+L) secondary antibody labeled with our superior CF® dyes.



### Product Description

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

- Available in 3 bright and photostable CF® dyes
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

### Product attributes

<b>Antibody type</b>	Secondary
<b>Clonality</b>	Polyclonal
<b>Host species</b>	Rabbit
<b>Antibody reactivity (target)</b>	Guinea Pig IgG
<b>Species reactivity</b>	Guinea pig
<b>Cross adsorption</b>	Not cross-adsorbed
<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>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 (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

Conjugation	Ex/Em	Size	Catalog No.
CF®543	541/560 nm	50 uL (100 ug)	<a href="#">20336-50uL</a>
0.5 mL (1 mg)	<a href="#">20336-500uL</a>		
1 mg	<a href="#">20336-1mg</a>		
CF®680	681/698 nm	50 uL (100 ug)	<a href="#">20243-1</a>
0.5 mL (1 mg)	<a href="#">20243</a>		
1 mg	<a href="#">20243-1mg</a>		
CF®770	770/797 nm	50 uL (100 ug)	<a href="#">20244-1</a>
0.5 mL (1 mg)	<a href="#">20244</a>		
1 mg	<a href="#">20244-1mg</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 06:06:14 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/rabbit-anti-guinea-pig-igg/>