

## Rabbit Anti-Chicken IgY (H+L)

Rabbit anti-chicken IgY (H+L) secondary antibody labeled with our superior CF® dyes.



### Product Description

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

- 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 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.

### Product attributes

Antibody type	Secondary
Clonality	Polyclonal
Host species	Rabbit
Antibody reactivity (target)	Chicken IgY
Species reactivity	Chicken
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

### References

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

Conjugation	Ex/Em	Size	Catalog No.
CF®488A	490/515 nm	50 uL (100 ug)	<a href="#">20079-1</a>
		0.5 mL (1 mg)	<a href="#">20079</a>
		1 mg	<a href="#">20079-1mg</a>
CF®543	541/560 nm	50 uL (100 ug)	<a href="#">20312-1</a>
		0.5 mL (1 mg)	<a href="#">20312</a>
		1 mg	<a href="#">20312-1mg</a>
CF®594	593/614 nm	50 uL (100 ug)	<a href="#">20164-1</a>
		0.5 mL (1 mg)	<a href="#">20164</a>
		1 mg	<a href="#">20164-1mg</a>
CF®633	630/650 nm	50 uL (100 ug)	<a href="#">20165-1</a>
		0.5 mL (1 mg)	<a href="#">20165</a>
		1 mg	<a href="#">20165-1mg</a>

This datasheet was generated on July 17, 2024 at 09:18:59 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/rabbit-anti-chicken-igy-hl/>