



## 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 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](#).

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