

## Goat Anti-Mouse IgG (H+L), Highly Cross-Adsorbed

Highly cross-adsorbed goat anti-mouse IgG (H L) secondary antibody labeled with our superior CF® Dyes.



### Product Description

This is a highly cross-adsorbed goat anti-mouse IgG (H L) secondary antibody labeled with our bright and photostable CF® Dyes. To minimize cross-reactivity, the antibody has been adsorbed against bovine, horse, human, rabbit, and swine serum. This antibody is not cross-adsorbed against rat. For staining rat tissue, or co-staining with rat antibodies, we recommend using [Goat Anti-Mouse IgG \(H L\), Highly Cross-Adsorbed \(Min x Rat\)](#).

- Highly-cross adsorbed for specific staining with minimal background
- Available with 19 bright and photostable CF® Dyes
- Biotin and HRP conjugates also available
- 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](#).

### Product attributes

|                                     |   |
|-------------------------------------|---|
| Antibody type                       | Secondary   |
| Clonality                           | Polyclonal  |
| Host species                        | Goat  |
| Antibody reactivity (target)        | Mouse IgG   |
| Species reactivity                  | Mouse   |
| Cross adsorption                    | Bovine, Horse, Human, Rabbit, Swine   |
| Concentration                       | 2 mg/mL, 1 mg/mL (HRP, AP conjugates)   |
| 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, HRP conjugates: PBS/50% glycerol/15 mg/mL BSA, HRP conjugates (lyophilized): PBS/10 mg/mL trehalose after reconstitution |
| Secondary/tag antibody applications | ELISA, Flow cytometry, IHC, IF (cells or tissue sections), Western blot   |

| Conjugation | Ex/Em      | Size  | Catalog No.  | Dye Features                     |
|-------------|------------|---|--|----------------------------------|
| CF@350      | 347/448 nm | 50 uL (100 ug)<br>0.5 mL (1 mg)<br>1 mg (lyophilized) | <a href="#">20143-1</a><br><a href="#">20143</a><br><a href="#">20143-1mg</a>          | <a href="#">CF@350 Features</a>  |
| CF@405M     | 408/452 nm | 50 uL (100 ug)<br>0.5 mL (1 mg)<br>1 mg (lyophilized) | <a href="#">20182-1</a><br><a href="#">20182</a><br><a href="#">20182-1mg</a>          | <a href="#">CF@405M Features</a> |
| CF@430      | 426/498 nm | 50 uL (100 ug)<br>0.5 mL (1 mg)<br>1 mg (lyophilized) | <a href="#">20459-50uL</a><br><a href="#">20459-500uL</a><br><a href="#">20459-1mg</a> | <a href="#">CF@430 Features</a>  |
| CF@488A     | 490/515 nm | 50 uL (100 ug)<br>0.5 mL (1 mg)<br>1 mg (lyophilized) | <a href="#">20018-1</a><br><a href="#">20018</a><br><a href="#">20018-1mg</a>          | <a href="#">CF@488A Features</a> |
| CF@532      | 527/558 nm | 50 uL (100 ug)<br>0.5 mL (1 mg)<br>1 mg (lyophilized) | <a href="#">20468-50uL</a><br><a href="#">20468-500uL</a><br><a href="#">20468-1mg</a> | <a href="#">CF@532 Features</a>  |
| CF@543      | 541/560 nm | 50 uL (100 ug)<br>0.5 mL (1 mg)<br>1 mg (lyophilized) | <a href="#">20299-1</a><br><a href="#">20299</a><br><a href="#">20299-1mg</a>          | <a href="#">CF@543 Features</a>  |
| CF@555      | 555/565 nm | 50 uL (100 ug)<br>0.5 mL (1 mg)<br>1 mg (lyophilized) | <a href="#">20231-1</a><br><a href="#">20231</a><br><a href="#">20231-1mg</a>          | <a href="#">CF@555 Features</a>  |
| CF@568      | 562/583 nm | 50 uL (100 ug)<br>0.5 mL (1 mg)<br>1 mg (lyophilized) | <a href="#">20101-1</a><br><a href="#">20101</a><br><a href="#">20101-1mg</a>          | <a href="#">CF@568 Features</a>  |
| CF@583R     | 585/609 nm | 50 uL (100 ug)<br>0.5 mL (1 mg)                       | <a href="#">20904-50uL</a><br><a href="#">20904-500uL</a>                              | <a href="#">CF@583R Features</a> |
| CF@594      | 593/614 nm | 50 uL (100 ug)<br>0.5 mL (1 mg)<br>1 mg (lyophilized) | <a href="#">20111-1</a><br><a href="#">20111</a><br><a href="#">20111-1mg</a>          | <a href="#">CF@594 Features</a>  |
| CF@633      | 630/650 nm | 50 uL (100 ug)<br>0.5 mL (1 mg)<br>1 mg (lyophilized) | <a href="#">20121-1</a><br><a href="#">20121</a><br><a href="#">20121-1mg</a>          | <a href="#">CF@633 Features</a>  |
| CF@640R     | 642/662 nm | 50 uL (100 ug)<br>0.5 mL (1 mg)<br>1 mg (lyophilized) | <a href="#">20175-1</a><br><a href="#">20175</a><br><a href="#">20175-1mg</a>          | <a href="#">CF@640R Features</a> |
| CF@647      | 650/665 nm | 50 uL (100 ug)<br>0.5 mL (1 mg)<br>1 mg (lyophilized) | <a href="#">20281-1</a><br><a href="#">20281</a><br><a href="#">20281-1mg</a>          | <a href="#">CF@647 Features</a>  |
| CF@660C     | 667/685 nm | 50 uL (100 ug)<br>0.5 mL (1 mg)<br>1 mg (lyophilized) | <a href="#">20052-1</a><br><a href="#">20052</a><br><a href="#">20052-1mg</a>          | <a href="#">CF@660C Features</a> |
| CF@680R     | 680/701 nm | 50 uL (100 ug)<br>0.25 mL (500 ug)                    | <a href="#">20192-1</a><br><a href="#">20192</a>                                       | <a href="#">CF@680 Features</a>  |
| CF@680      | 681/698 nm | 50 uL (100 ug)<br>0.25 mL (500 ug)                    | <a href="#">20065-1</a><br><a href="#">20065</a>                                       | <a href="#">CF@680R Features</a> |
| CF@750      | 755/777 nm | 50 uL (100 ug)<br>250 uL (100 ug)                     | <a href="#">20463-50uL</a><br><a href="#">20463-250uL</a>                              | <a href="#">CF@750 Features</a>  |
| CF@770      | 770/797 nm | 50 uL (100 ug)<br>0.25 mL (500 ug)                    | <a href="#">20077-1</a><br><a href="#">20077</a>                                       | <a href="#">CF@770 Features</a>  |
| CF@790      | 784/806 nm | 50 uL (100 ug)  | <a href="#">20342-50uL</a>   | <a href="#">CF@790 Features</a>  |
| Biotin      | N/A        | 50 uL (100 ug)<br>0.5 mL (1 mg)<br>1 mg (lyophilized) | <a href="#">20184-1</a><br><a href="#">20184</a><br><a href="#">20184-1mg</a>          |                                  |
| HRP         | N/A        | 100 uL (100 ug)<br>1 mL (1 mg)<br>1 mg (lyophilized)  | <a href="#">20401-100uL</a><br><a href="#">20401-1mL</a><br><a href="#">20401-1mg</a>  |                                  |

This datasheet was generated on June 30, 2024 at 08:22:23 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/goat-anti-mouse-igg-hl-highly-cross-absorbed/>