



Glowing products for science

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

Goat anti-guinea pig IgG (H L) secondary antibody labeled with our superior CF® Dyes.



### Product Description

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

- Available in 10 bright and photostable CF® Dyes
- 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.

Call us : [800-304-5357](tel:800-304-5357) Email: [btinfo@biotium.com](mailto:btinfo@biotium.com)

### Product attributes

|  |  |
|--|--|
| <b>Antibody type</b>                       | Secondary  |
| <b>Clonality</b>                           | Polyclonal   |
| <b>Host species</b>                        | Goat   |
| <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,<br>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 ( <i>Bos taurus</i> ), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot. |

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

| Conjugation | Ex/Em      | Size           | Catalog No.                 | Dye Features                     |
|-------------|------------|----------------|-----------------------------|----------------------------------|
| CF@350      | 347/448 nm | 50 uL (100 ug) | <a href="#">20198-1</a>     | <a href="#">CF@350 Features</a>  |
|             |            | 0.5 mL (1 mg)  | <a href="#">20198</a>       |                                  |
|             |            | 1 mg           | <a href="#">20198-1mg</a>   |                                  |
| CF@488A     | 490/515 nm | 50 uL (100 ug) | <a href="#">20017-1</a>     | <a href="#">CF@488A Features</a> |
|             |            | 0.5 mL (1 mg)  | <a href="#">20017</a>       |                                  |
|             |            | 1 mg           | <a href="#">20017-1mg</a>   |                                  |
| CF@543      | 541/560 nm | 50 uL (100 ug) | <a href="#">20317-1</a>     | <a href="#">CF@543 Features</a>  |
|             |            | 0.5 mL (1 mg)  | <a href="#">20317</a>       |                                  |
|             |            | 1 mg           | <a href="#">20317-1mg</a>   |                                  |
| CF@555      | 555/565 nm | 50 uL (100 ug) | <a href="#">20036-1</a>     | <a href="#">CF@555 Features</a>  |
|             |            | 0.5 mL (1 mg)  | <a href="#">20036</a>       |                                  |
|             |            | 1 mg           | <a href="#">20036-1mg</a>   |                                  |
| CF@568      | 562/583 nm | 50 uL (100 ug) | <a href="#">20108-1</a>     | <a href="#">CF@568 Features</a>  |
|             |            | 0.5 mL (1 mg)  | <a href="#">20108</a>       |                                  |
|             |            | 1 mg           | <a href="#">20108-1mg</a>   |                                  |
| CF@583R     | 585/609 nm | 50 uL (100 ug) | <a href="#">20899-50uL</a>  | <a href="#">CF@583R Features</a> |
|             |            | 0.5 mL (1 mg)  | <a href="#">20899-500uL</a> |                                  |
| CF@594      | 593/614 nm | 50 uL (100 ug) | <a href="#">20118-1</a>     | <a href="#">CF@594 Features</a>  |
|             |            | 0.5 mL (1 mg)  | <a href="#">20118</a>       |                                  |
|             |            | 1 mg           | <a href="#">20118-1mg</a>   |                                  |
| CF@633      | 630/650 nm | 50 uL (100 ug) | <a href="#">20129-1</a>     | <a href="#">CF@633 Features</a>  |
|             |            | 0.5 mL (1 mg)  | <a href="#">20129</a>       |                                  |
|             |            | 1 mg           | <a href="#">20129-1mg</a>   |                                  |
| CF@640R     | 642/662 nm | 50 uL (100 ug) | <a href="#">20085-1</a>     | <a href="#">CF@640R Features</a> |
|             |            | 0.5 mL (1 mg)  | <a href="#">20085</a>       |                                  |
|             |            | 1 mg           | <a href="#">20085-1mg</a>   |                                  |
| CF@647      | 650/665 nm | 50 uL (100 ug) | <a href="#">20041-1</a>     | <a href="#">CF@647 Features</a>  |
|             |            | 0.5 mL (1 mg)  | <a href="#">20041</a>       |                                  |
|             |            | 1 mg           | <a href="#">20041-1mg</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 curated [CF@ Dye references](#).

This datasheet was generated on November 21, 2024 at 07:51:46 AM. Visit product page to check for updated information before use.

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