



Glowing products for science

Goat Anti-Swine IgG (H+L)

Goat anti-swine IgG (H L) secondary antibody labeled with our superior CF® Dyes and other labels.



Product Description

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

- Available in 9 bright and photostable CF® Dyes
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

Call us : [800-304-5357](tel:800-304-5357)

Product attributes

| | |
|--|---|
| Antibody type | Secondary |
| Clonality | Polyclonal |
| Host species | Goat |
| Antibody reactivity (target) | Swine IgG |
| Species reactivity | Swine |
| 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 origin | 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. |

Goat Anti-Swine IgG (H+L)

| Conjugation | Ex/Em | Size | Catalog No. | Dye Features |
|---------------|-----------------------------|----------------|----------------------------|----------------------------------|
| CF@488A | 490/515 nm | 50 uL (100 ug) | 20028-1 | CF@488A Features |
| 0.5 mL (1 mg) | 20028 | | | |
| 1 mg | 20028-1mg | | | |
| CF@543 | 541/560 nm | 50 uL (100 ug) | 20324-1 | CF@543 Features |
| 0.5 mL (1 mg) | 20324 | | | |
| 1 mg | 20324-1mg | | | |
| CF@555 | 555/565 nm | 50 uL (100 ug) | 20236-1 | CF@555 Features |
| 0.5 mL (1 mg) | 20236 | | | |
| 1 mg | 20236-1mg | | | |
| CF@568 | 562/583 nm | 50 uL (100 ug) | 20091-1 | CF@568 Features |
| 0.5 mL (1 mg) | 20091 | | | |
| 1 mg | 20091-1mg | | | |
| CF@583R | 585/609 nm | 50 uL (100 ug) | 20907-50uL | CF@583R Features |
| 0.5 mL (1 mg) | 20907-500uL | | | |
| CF@594 | 593/614 nm | 50 uL (100 ug) | 20160-1 | CF@594 Features |
| 0.5 mL (1 mg) | 20160 | | | |
| 1 mg | 20160-1mg | | | |
| CF@633 | 630/650 nm | 50 uL (100 ug) | 20138-1 | CF@633 Features |
| 0.5 mL (1 mg) | 20138 | | | |
| 1 mg | 20138-1mg | | | |
| CF@640R | 642/662 nm | 50 uL (100 ug) | 20089-1 | CF@640R Features |
| 0.5 mL (1 mg) | 20089 | | | |
| 1 mg | 20089-1mg | | | |
| CF@647 | 650/665 nm | 50 uL (100 ug) | 20286-1 | CF@647 Features |
| 0.5 mL (1 mg) | 20286 | | | |
| 1 mg | 20286-1mg | | | |

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 dH₂O

HRP or DNP conjugates: add 1 mL dH₂O

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 December 24, 2025 at 12:27:09 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/goat-anti-swine-igg-hl/>