

Goat Anti-Human IgG (H+L), Fc Gamma Specific, Highly Cross-Adsorbed



Highly cross-adsorbed goat anti-human IgG (H+L), Fc gamma specific secondary antibody labeled with our superior CF® dyes.

Product Description

This is a highly cross-adsorbed goat anti-human IgG (H L), Fc gamma specific secondary antibody labeled with our superior CF® dyes. The antibody reacts specifically with the human IgG heavy chains and not with light chains or F(ab')₂ fragment of human immunoglobulins. To minimize cross-reactivity, the antibodies are cross-adsorbed against bovine, horse, mouse and rabbit serum IgG prior to conjugation.

- Highly cross-adsorbed for specific staining with minimal background
- Available in 5 bright and photostable CF® dyes
- Suitable for western, immunofluorescence, and immunohistology

Product attributes

| | |
|--|---|
| Clonality | Polyclonal |
| Antibody type | Secondary, Anti-Human Immunoglobulin |
| 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 |
| Species reactivity | Human |
| 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. |
| Host species | Goat |
| Antibody reactivity (target) | Human IgG Fc |
| Cross adsorption | Bovine, Horse, Mouse, Rabbit |

| Conjugation | Ex/Em | Size | Catalog No. | Dye Features |
|--|--|----------------|----------------------------|----------------------------------|
| CF@488A 0.5 mL (1 mg) 1 mg (lyophilized) | 490/515 nm 20444-500uL 20444-1mg | 50 uL (100 ug) | 20444-50uL | CF@488A Features |
| CF@543 0.5 mL (1 mg) 1 mg (lyophilized) | 541/560 nm 20445-500uL 20445-1mg | 50 uL (100 ug) | 20445-50uL | CF@543 Features |
| CF@594 0.5 mL (1 mg) 1 mg (lyophilized) | 593/614 nm 20446-500uL 20446-1mg | 50 uL (100 ug) | 20446-50uL | CF@594 Features |
| CF@640R 0.5 mL (1 mg) 1 mg (lyophilized) | 642/662 nm 20447-500uL 20447-1mg | 50 uL (100 ug) | 20447-50uL | CF@640R Features |
| CF@647 0.5 mL (1 mg) 1 mg (lyophilized) | 650/665 nm 20448-500uL 20448-1mg | 50 uL (100 ug) | 20448-50uL | CF@647 Features |

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 June 3, 2026 at 03:38:35 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/goat-anti-human-igg-hl-fc-gamma-specific-highly-cross-adsorbed/>