

Goat Anti-Mouse IgG1 (γ1)

Goat anti-mouse IgG1 isotype-specific secondary antibody labeled with our superior CF® dyes and biotin.



Product Description

This is a goat anti-mouse IgG1 isotype-specific secondary antibody labeled with our bright and photostable CF® Dyes and biotin. The conjugates are prepared from affinity-purified antibodies that react with Fc portion of the heavy chain of mouse IgG1. To minimize cross reactivity, the antibodies are cross-adsorbed against other mouse IgG subclasses (IgG2a, IgG2b, IgG3), and human, bovine and rabbit serum proteins.

- Cross-adsorbed for specific staining with minimal background
- Available in 12 bright and photostable CF® Dyes
- Biotin conjugate also available
- Suitable for western, immunofluorescence, and immunohistology

Note: Conjugates of blue fluorescent dyes like CF®350 and CF®405S 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 bright and specific [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.

Product attributes

| | |
|-------------------------------------|---|
| Antibody type | Secondary |
| Clonality | Polyclonal |
| Host species | Goat |
| Antibody reactivity (target) | Mouse IgG1 (γ1) |
| Species reactivity | Mouse |
| Cross adsorption | Bovine, Human, Rabbit |
| 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 | ELISA, Flow cytometry, IHC, IF (cells or tissue sections), Western blot |

References

Download a list of [CF® dye references](#).

| Conjugation | Ex/Em | Size | Catalog No. |
|-------------|------------|------------------------------------|---|
| CF®350 | 347/448 nm | 50 uL (100 ug) 0.25 mL (500 ug) | 20245-1 20245 |
| CF®405S | 404/431 nm | 50 uL (100 ug) 0.25 mL (500 ug) | 20380-50uL 20380-250uL |
| CF®488A | 490/515 nm | 50 uL (100 ug) 0.25 mL (500 ug) | 20246-1 20246 |
| CF®543 | 541/560 nm | 50 uL (100 ug) 0.25 mL (500 ug) | 20325-1 20325 |
| CF®555 | 555/565 nm | 50 uL (100 ug) 0.25 mL (500 ug) | 20247-1 20247 |
| CF®568 | 562/583 nm | 50 uL (100 ug) 0.25 mL (500 ug) | 20248-1 20248 |
| CF®594 | 593/614 nm | 50 uL (100 ug) 0.25 mL (500 ug) | 20249-1 20249 |
| CF®633 | 630/650 nm | 50 uL (100 ug) 0.25 mL (500 ug) | 20250-1 20250 |
| CF®640R | 642/662 nm | 50 uL (100 ug) 0.25 mL (500 ug) | 20251-1 20251 |
| CF®647 | 650/665 nm | 50 uL (100 ug) 0.25 mL (500 ug) | 20252-1 20252 |
| CF®680 | 681/698 nm | 50 uL (100 ug) 0.25 mL (500 ug) | 20253-1 20253 |
| CF®770 | 770/797 nm | 50 uL (100 ug) 0.25 mL (500 ug) | 20254-1 20254 |
| Biotin | N/A | 50 uL (100 ug) 0.25 mL (500 ug) | 20471-50uL 20471-250uL |