

Goat Anti-Rabbit IgG (H+L)

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



Product Description

This is a goat anti-rabbit IgG (H L) secondary antibody labeled with our superior CF® dyes and other labels.

- Available in 18 bright and photostable CF® dyes
- Alkaline phosphatase, APC, HRP, R-PE and biotin conjugates also available
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

Note: Conjugates of blue fluorescent dyes like CF®350, CF®405S, and CF®405M 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 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 curated [CF® Dye references](#).

Product attributes

| | |
|-------------------------------------|---|
| Antibody type | Secondary |
| Clonality | Polyclonal |
| Host species | Goat |
| Antibody reactivity (target) | Rabbit IgG |
| Species reactivity | Rabbit |
| Cross adsorption | Not cross-adsorbed |
| 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. |
|-------------|----------------------|--------------------|-----------------------------|
| CF@350 | 347/448 nm | 50 uL (100 ug) | 20141-1 |
| | | 0.5 mL (1 mg) | 20141 |
| | | 1 mg (lyophilized) | 20141-1mg |
| CF@405S | 404/431 nm | 50 uL (100 ug) | 20082-1 |
| | | 0.5 mL (1 mg) | 20082 |
| | | 1 mg (lyophilized) | 20082-1mg |
| CF@405M | 408/452 nm | 50 uL (100 ug) | 20181-1 |
| | | 0.5 mL (1 mg) | 20181 |
| | | 1 mg (lyophilized) | 20181-1mg |
| CF@405L | 395/545 nm | 50 uL (100 ug) | 20409-50uL |
| | | 0.5 mL (1 mg) | 20409-500uL |
| | | 1 mg (lyophilized) | 20409-1mg |
| CF@488A | 490/515 nm | 50 uL (100 ug) | 20012-1 |
| | | 0.5 mL (1 mg) | 20012 |
| | | 1 mg (lyophilized) | 20012-1mg |
| CF@514 | 516/548 nm | 50 uL (100 ug) | 20387-50uL |
| | | 0.5 mL (1 mg) | 20387-500uL |
| | | 1 mg (lyophilized) | 20387-1mg |
| CF@532 | 527/558 nm | 50 uL (100 ug) | 20366-50uL |
| | | 0.5 mL (1 mg) | 20366-500uL |
| | | 1 mg (lyophilized) | 20366-1mg |
| CF@543 | 541/560 nm | 50 uL (100 ug) | 20309-1 |
| | | 0.5 mL (1 mg) | 20309 |
| | | 1 mg (lyophilized) | 20309-1mg |
| CF@555 | 555/565 nm | 50 uL (100 ug) | 20033-1 |
| | | 0.5 mL (1 mg) | 20033 |
| | | 1 mg (lyophilized) | 20033-1mg |
| CF@568 | 562/583 nm | 50 uL (100 ug) | 20102-1 |
| | | 0.5 mL (1 mg) | 20102 |
| | | 1 mg (lyophilized) | 20102-1mg |
| CF@594 | 593/614 nm | 50 uL (100 ug) | 20112-1 |
| | | 0.5 mL (1 mg) | 20112 |
| | | 1 mg (lyophilized) | 20112-1mg |
| CF@633 | 630/650 nm | 50 uL (100 ug) | 20122-1 |
| | | 0.5 mL (1 mg) | 20122 |
| | | 1 mg (lyophilized) | 20122-1mg |
| CF@640R | 642/662 nm | 50 uL (100 ug) | 20202-1 |
| | | 0.5 mL (1 mg) | 20202 |
| | | 1 mg (lyophilized) | 20202-1mg |
| CF@647 | 650/665 nm | 50 uL (100 ug) | 20043-1 |
| | | 0.5 mL (1 mg) | 20043 |
| | | 1 mg (lyophilized) | 20043-1mg |
| CF@660C | 667/685 nm | 50 uL (100 ug) | 20053-1 |
| | | 0.5 mL (1 mg) | 20053 |
| | | 1 mg (lyophilized) | 20053-1mg |
| CF@660R | 663/682 nm | 50 uL (100 ug) | 20055-1 |
| | | 0.5 mL (1 mg) | 20055 |
| | | 1 mg (lyophilized) | 20055-1mg |
| CF@750 | 755/777 nm | 50 uL (100 ug) | 20073-1 |
| | | 0.5 mL (1 mg) | 20073 |
| | | 1 mg (lyophilized) | 20073-1mg |
| CF@790 | 784/806 nm | 50 uL (100 ug) | 20379-50uL |
| | | 1 mL (500 ug) | 20353-1mL |
| R-PE | 496, 546, 565/578 nm | 200 uL (100 ug) | 20353-200uL |
| APC | 650/660 nm | 100 uL (50 ug) | 20412-100uL |
| | | 0.5 mL (250 ug) | 20412-500uL |
| Biotin | N/A | 50 uL (100 ug) | 20185 |
| | | 0.5 mL (1 mg) | 20185-1 |
| | | 1 mg (lyophilized) | 20185-1mg |
| HRP | N/A | 100 uL (100 ug) | 20402-100uL |
| | | 1 mL (1 mg) | 20402-1mL |
| | | 1 mg (lyophilized) | 20402-1mg |
| AP | N/A | 100 uL (100 ug) | 20465-100uL |
| | | 1 mL (1 mg) | 20465-1mL |