

## Rabbit Anti-Sheep IgG (H+L), Highly Cross-Adsorbed

Highly cross-adsorbed rabbit anti-sheep IgG (H+L) secondary antibody labeled with our superior CF® dyes.



### Product Description

This is a highly cross-adsorbed rabbit anti-sheep IgG (H L) secondary antibody labeled with our bright and photostable CF® dyes. To minimize cross-reactivity, the antibody has been adsorbed against human serum.

- Highly cross-adsorbed for specific staining with minimal background
- Available in 4 bright and photostable CF® dyes
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

### Product attributes

|                                     |   |
|-------------------------------------|---|
| Antibody type                       | Secondary   |
| Clonality                           | Polyclonal  |
| Host species                        | Rabbit  |
| Antibody reactivity (target)        | Sheep IgG   |
| Species reactivity                  | Sheep   |
| Cross adsorption                    | Human   |
| 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. |

## Rabbit Anti-Sheep IgG (H+L), Highly Cross-Adsorbed

| Conjugation | Ex/Em      | Size           | Catalog No.               |
|-------------|------------|----------------|---------------------------|
| CF®488A     | 490/515 nm | 50 uL (100 ug) | <a href="#">20172-1</a>   |
|             |            | 0.5 mL (1 mg)  | <a href="#">20172</a>     |
|             |            | 1 mg           | <a href="#">20172-1mg</a> |
| CF®543      | 541/560 nm | 50 uL (100 ug) | <a href="#">20323-1</a>   |
|             |            | 0.5 mL (1 mg)  | <a href="#">20323</a>     |
|             |            | 1 mg           | <a href="#">20323-1mg</a> |
| CF®594      | 593/614 nm | 50 uL (100 ug) | <a href="#">20173-1</a>   |
|             |            | 0.5 mL (1 mg)  | <a href="#">20173</a>     |
|             |            | 1 mg           | <a href="#">20173-1mg</a> |
| CF®633      | 630/650 nm | 50 uL (100 ug) | <a href="#">20174-1</a>   |
|             |            | 0.5 mL (1 mg)  | <a href="#">20174</a>     |
|             |            | 1 mg           | <a href="#">20174-1mg</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 [CF® dye references](#).

This datasheet was generated on August 2, 2025 at 07:00:59 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/rabbit-anti-sheep-igg-hl-highly-cross-adsorbed/>