

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) | 20172-1 |
| | | 0.5 mL (1 mg) | 20172 |
| | | 1 mg | 20172-1mg |
| CF®543 | 541/560 nm | 50 uL (100 ug) | 20323-1 |
| | | 0.5 mL (1 mg) | 20323 |
| | | 1 mg | 20323-1mg |
| CF®594 | 593/614 nm | 50 uL (100 ug) | 20173-1 |
| | | 0.5 mL (1 mg) | 20173 |
| | | 1 mg | 20173-1mg |
| CF®633 | 630/650 nm | 50 uL (100 ug) | 20174-1 |
| | | 0.5 mL (1 mg) | 20174 |
| | | 1 mg | 20174-1mg |

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 November 21, 2024 at 02:32:36 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/>