

Donkey Anti-Human IgG (H+L), Highly Cross-Adsorbed

Highly cross-adsorbed donkey anti-human IgG (H+L) secondary antibody labeled with our superior CF® Dyes.



Product Description

This is a highly cross-adsorbed donkey anti-human IgG (H L) secondary antibody labeled with our bright and photostable CF® Dyes. To minimize cross-reactivity, the antibody has been adsorbed against bovine, chicken, goat, guinea pig, horse, mouse, rabbit, rat, sheep, and Syrian hamster serum.

- Highly cross-adsorbed for specific staining with minimal background
- Available in 6 bright and photostable CF® Dyes
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

Product attributes

| | |
|--|--|
| Antibody type | Secondary, Anti-Human Immunoglobulin |
| Clonality | Polyclonal |
| Host species | Donkey |
| Antibody reactivity (target) | Human IgG |
| Species reactivity | Human |
| Cross adsorption | Bovine, Chicken, Goat, Guinea pig, Horse, Mouse, Rabbit, Rat, Sheep, Syrian hamster |
| 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 (<i>Bos taurus</i>), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot. |

Donkey Anti-Human IgG (H+L), Highly Cross-Adsorbed

| Conjugation | Ex/Em | Size | Catalog No. |
|-------------|------------|------------------|---------------------------|
| CF®488A | 490/515 nm | 50 uL (100 ug) | 20074-1 |
| | | 0.5 mL (1 mg) | 20074 |
| | | 1 mg | 20074-1mg |
| CF®543 | 541/560 nm | 50 uL (100 ug) | 20318-1 |
| | | 0.5 mL (1 mg) | 20318 |
| | | 1 mg | 20318-1mg |
| CF®594 | 593/614 nm | 50 uL (100 ug) | 20075-1 |
| | | 0.5 mL (1 mg) | 20075 |
| | | 1 mg | 20075-1mg |
| CF®633 | 630/650 nm | 50 uL (100 ug) | 20076-1 |
| | | 0.5 mL (1 mg) | 20076 |
| | | 1 mg | 20076-1mg |
| CF®680 | 681/698 nm | 50 uL (100 ug) | 20278-1 |
| | | 0.25 mL (500 ug) | 20278 |
| CF®770 | 770/797 nm | 50 uL (100 ug) | 20279-1 |
| | | 0.25 mL (500 ug) | 20279 |

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 22, 2024 at 11:59:44 PM. Visit product page to check for updated information before use.

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