



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357) Email: btinfo@biotium.com

Drop-n-Stain™ CF® Dye Donkey Anti-Rabbit IgG (H+L), Highly Cross-Adsorbed



CF® dye labeled, highly cross-adsorbed, donkey anti-rabbit IgG (H+L) secondary antibody provided in a convenient dropper bottle.

Product Description

This is a CF® dye labeled, highly cross-adsorbed, donkey anti-rabbit IgG (H L) secondary antibody provided in a convenient dropper bottle for easy dispensing. To minimize cross-reactivity, the antibody has been adsorbed against bovine, chicken, goat, guinea pig, Syrian hamster, horse, human, mouse, rat, and sheep serum.

- Highly cross-adsorbed for specific staining with minimal background
- Convenient dropper bottle for easy dispensing without pipetting
- Available with 4 bright and photostable CF® dyes
- Suitable for western, immunofluorescence, and immunohistology

See our other [Drop-n-Stain™ CF® Dye Secondary Antibodies](#) 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](#). CF is a registered trademark of Biotium, Inc.

References

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

Product attributes

| | |
|-------------------------------------|---|
| Clonality | Polyclonal |
| Antibody type | Secondary |
| Host species | Donkey |
| Antibody reactivity (target) | Rabbit IgG |
| Cross adsorption | Bovine, Chicken, Goat, Guinea pig, Horse, Human, Mouse, Sheep, Syrian hamster |
| Concentration | Use 2 drops/mL for IF, or 1 drop per tube for flow cytometry |
| Secondary/tag antibody applications | Flow cytometry, IF (cells or tissue sections) |

| Conjugation | Ex/Em | Size | Catalog No. | Dye Features |
|-------------------------|------------|------|-----------------------|----------------------------------|
| CF@488A | 490/515 nm | 5 mL | 20950 | CF@488A Features |
| CF@543 | 541/560 nm | 5 mL | 20966 | CF@543 Features |
| CF@594 | 593/614 nm | 5 mL | 20951 | CF@594 Features |
| CF@640R | 642/662 nm | 5 mL | 20962 | CF@640R Features |

This datasheet was generated on July 24, 2024 at 05:10:27 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/drop-n-stain-cf-dye-donkey-anti-rabbit-igg-hl-highly-cross-adsorbed/>