



Glowing products for science

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

## Drop-n-Stain™ CF® Dye Goat Anti-Mouse IgG (H+L), Highly Cross-Adsorbed



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

### Product attributes

|                                     |  |
|-------------------------------------|--|
| Clonality                           | Polyclonal   |
| Antibody reactivity (target)        | Mouse IgG  |
| Host species                        | Goat   |
| Cross adsorption                    | Bovine, Horse, Human, Rabbit, Swine                          |
| Antibody type                       | Secondary  |
| 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)                |

### Product Description

This is a CF® dye labeled, highly cross-adsorbed, goat anti-mouse IgG (H L) secondary antibody provided in a convenient dropper bottle for easy dispensing. To minimize cross-reactivity, the antibody has been adsorbed against human, mouse, and rat 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

## Drop-n-Stain™ CF® Dye Goat Anti-Mouse IgG (H+L), Highly Cross-Adsorbed

| Conjugation             | Ex/Em      | Size | Catalog No.           | Dye Features                     |
|-------------------------|------------|------|-----------------------|----------------------------------|
| <a href="#">CF@488A</a> | 490/515 nm | 5 mL | <a href="#">20956</a> | <a href="#">CF@488A Features</a> |
| <a href="#">CF@543</a>  | 541/560 nm | 5 mL | <a href="#">20969</a> | <a href="#">CF@543 Features</a>  |
| <a href="#">CF@594</a>  | 593/614 nm | 5 mL | <a href="#">20957</a> | <a href="#">CF@594 Features</a>  |
| <a href="#">CF@640R</a> | 642/662 nm | 5 mL | <a href="#">20965</a> | <a href="#">CF@640R Features</a> |

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](#).

This datasheet was generated on December 21, 2024 at 05:17:27 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/drop-n-stain-cf-dye-goat-anti-mouse-igg-hl-highly-cross-adsorbed/>