



Glowing products for science

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

### 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)

## 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 July 31, 2025 at 02:51:28 AM. 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/>