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

The same of the sa

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

## **Product Description**

This is a CF® dye labeled, highly cross-adsorbed, donkey 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 bovine, chicken, goat, guinea pig, Syrian hamster, horse, human, rabbit, 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

#### **Product attributes**

Clonality	Polyclonal		
Antibody type	Secondary		
Host species	Donkey		
Antibody reactivity (target)	Mouse IgG		
Cross adsorption	Bovine, Chicken, Goat, Guinea pig, Horse, Human, Rabbit, 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)		

Call us: 800-304-5357 Email: btinfo@biotium.com

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

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF®488A	490/515 nm	5 mL	<u>20952</u>	CF®488A Features
<u>CF®543</u>	541/560 nm	5 mL	<u>20967</u>	CF®543 Features
CF®594	593/614 nm	5 mL	<u>20953</u>	CF®594 Features
CF®640R	642/662 nm	5 mL	<u>20963</u>	CF®640R Features

See our other <u>Drop-n-Stain™ CF® Dye Secondary Antibodies</u>

View our full selection of <u>Secondary Antibodies</u>, or search our catalog using our <u>Antibody Finder</u>. Alternatively, you can view our <u>secondary antibody product listings</u> 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 November 17, 2025 at 07:44:28 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/drop-n-stain-cf-dye-donkey-anti-mouse-igg-hl-highly-cross-adsorbed/">https://biotium.com/product/drop-n-stain-cf-dye-donkey-anti-mouse-igg-hl-highly-cross-adsorbed/</a>