



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

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



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

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, IHC, IF (cells or tissue sections), Western blot

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

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	20952	CF@488A Features
CF@543	541/560 nm	5 mL	20967	CF@543 Features
CF@594	593/614 nm	5 mL	20953	CF@594 Features
CF@640R	642/662 nm	5 mL	20963	CF@640R Features

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 June 15, 2026 at 08:27:50 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/drop-n-stain-cf-dye-donkey-anti-mouse-igg-hl-highly-cross-adsorbed/>