

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

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

### **Product Description**

This is a CF® dye labeled, highly cross-adsorbed, goat 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 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

Call us: 800-304-5357

Clonality	Polyclonal	
Antibody type	Secondary	
Host species	Goat	
Antibody reactivity (target)	Rabbit IgG	
Cross adsorption	Human, Mouse, Rat	
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)	

Email: techsupport@biotium.com

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

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF®488A	490/515 nm	5 mL	<u>20954</u>	CF®488A Features
<u>CF®543</u>	541/560 nm	5 mL	<u>20968</u>	CF®543 Features
CF®594	593/614 nm	5 mL	<u>20955</u>	CF®594 Features
CF®640R	642/662 nm	5 mL	<u>20964</u>	CF®640R Features

See our other Drop-n-Stain™ CF® Dye Secondary Antibodies

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 October 18, 2025 at 02:38:47 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/drop-n-stain-cf-dye-goat-anti-rabbit-igg-hl-highly-cross-adsorbed/">https://biotium.com/product/drop-n-stain-cf-dye-goat-anti-rabbit-igg-hl-highly-cross-adsorbed/</a>