

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 attributes

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

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

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	20954	CF@488A Features
CF@543	541/560 nm	5 mL	20968	CF@543 Features
CF@594	593/614 nm	5 mL	20955	CF@594 Features
CF@640R	642/662 nm	5 mL	20964	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 11, 2026 at 04:46:06 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/drop-n-stain-cf-dye-goat-anti-rabbit-igg-hl-highly-cross-adsorbed/>