

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, IF (cells or tissue sections) |

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 December 21, 2024 at 05:06:38 PM. 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/>