



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357) Email: [btinfo@biotium.com](mailto:btinfo@biotium.com)

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

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



## Product Description

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

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](#).

### Product attributes

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

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
<a href="#">CF®488A</a>	490/515 nm	5 mL	<a href="#">20950</a>	<a href="#">CF®488A Features</a>
<a href="#">CF®543</a>	541/560 nm	5 mL	<a href="#">20966</a>	<a href="#">CF®543 Features</a>
<a href="#">CF®594</a>	593/614 nm	5 mL	<a href="#">20951</a>	<a href="#">CF®594 Features</a>
<a href="#">CF®640R</a>	642/662 nm	5 mL	<a href="#">20962</a>	<a href="#">CF®640R Features</a>

This datasheet was generated on April 28, 2024 at 03:44:08 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/drop-n-stain-cf-dye-donkey-anti-rabbit-igg-hl-highly-cross-adsorbed/>