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

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)

Call us: 800-304-5357 Email: btinfo@biotium.com

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

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

See our other <u>Drop-n-Stain™ CF® Dye Secondary Antibodies</u>

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 November 17, 2025 at 07:47:20 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/