

# Donkey Anti-Guinea Pig IgG (H+L), Highly Cross-Adsorbed, CF® Dye Conjugates, Single Label for STORM



Highly cross-adsorbed donkey anti-guinea pig IgG (H+L) secondary antibody with single CF® Dye label for STORM super-resolution microscopy.

## Product Description

This is a highly cross-adsorbed donkey anti-guinea pig IgG (H L) secondary antibody that has a low degree of labeling (DOL) with one of our bright and photostable CF® Dyes.

- CF® Dye single label secondary antibody ideal for STORM imaging
- Highly cross-adsorbed for specific staining with minimal background
- Available in 3 bright and photostable CF® Dyes

Secondary antibodies with a low DOL, or number of dye molecules per antibody molecule, have been reported to be optimal for STORM ([Bittel et al. \(2015\) Proc. SPIE 9331](#)). This product is prepared by single labeling (DOL=1) of highly cross-adsorbed goat anti-mouse IgG (H L) with a selection of compatible CF® Dyes for **(d)-STORM super-resolution microscopy**. To minimize cross-reactivity, the antibody has been adsorbed against human, bovine, horse, rabbit, and swine serum. Learn more about [CF® Dyes for super-resolution microscopy](#). View our full selection of bright and specific [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](#). **Storage and Handling Liquid format:** Store at -20°C, protected from light. Product is stable for at least 6 months from date of receipt when stored as recommended. Liquid format antibodies contain 50% glycerol and will not freeze at -20°C. **Lyophilized format:** Store at -20°C, protected from light. Product is stable for at least 6 months from date of receipt when stored as recommended. Reconstitute antibodies in water using the indicated volumes below: CF® Dye and biotin conjugates: add 0.5 mL dH<sub>2</sub>O HRP or DNP conjugates: add 1 mL dH<sub>2</sub>O Add the indicated volume of water directly to the vial containing the lyophilized antibody and mix gently to dissolve. Store reconstituted antibody at -20°C and protect from light. Aliquot to avoid repeated freeze/thaw cycles. Alternatively, an equal volume of glycerol can be mixed with the reconstituted antibody so that it will remain liquid at -20°C. Optional: A preservative such as 0.05% sodium azide (final concentration) can be added to CF® Dye and biotin conjugates. Do not add sodium azide to HRP conjugates. **Note:** Storage of the antibody for more than a day at final working dilution is not recommended. CF is a registered trademark of Biotium, Inc.

## References

Download a list of curated [CF® Dye references](#).

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
<a href="#">CF®505</a>	505/519 nm	50 uL (50 ug) 0.5 mL (500 ug)	<a href="#">20881-50uL</a> <a href="#">20881-500uL</a>	
<a href="#">CF®568</a>	562/583 nm	50 uL (50 ug) 0.5 mL (500 ug)	<a href="#">20838-50uL</a> <a href="#">20838-500uL</a>	<a href="#">CF®568 Features</a>
<a href="#">CF®647</a>	650/665 nm	50 uL (50 ug) 0.5 mL (500 ug)	<a href="#">20837-50uL</a> <a href="#">20837-500uL</a>	<a href="#">CF®647 Features</a>

## Product attributes

Antibody type	Secondary
Cross adsorption	Bovine, Chicken, Goat, Horse, Human, Mouse, Rabbit, Rat, Sheep, Syrian hamster
Concentration	1 mg/mL
Clonality	Polyclonal
Host species	Donkey
Species reactivity	Guinea pig
Antibody/conjugate formulation	Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide
Secondary/tag antibody applications	IF (cells or tissue sections), STORM