

# Product Information

## Donkey Anti-Rabbit IgG (H+L), Whole Antibody, Highly Cross-Adsorbed (min x Bv, Ch, Gt, GP, Hs, Hu, Ms, Rt, Sh, SHm)

CF® Dye	Ex/Em	Catalog No.	Size
CF@350	347/448 nm	20351-1	50 uL
		20351	500 uL
		20351-1mg	1 mg (lyophilized)
CF@405S	404/431 nm	20420-50uL	50 uL
		20420-500uL	500 uL
		20420-1mg	1 mg (lyophilized)
CF@430	426/498 nm	20462-50uL	50 uL
		20462-500uL	500 uL
		20462-1mg	1 mg (lyophilized)
CF@488A	490/515 nm	20015-1	50 uL
		20015	500 uL
		20015-1mg	1 mg (lyophilized)
CF@543	541/560 nm	20308-1	50 uL
		20308	500 uL
		20308-1mg	1 mg (lyophilized)
CF@555	555/565 nm	20038-1	50 uL
		20038	500 uL
		20038-1mg	1 mg (lyophilized)
CF@568	562/583 nm	20098-1	50 uL
		20098	500 uL
		20098-1mg	1 mg (lyophilized)
CF@594	593/614 nm	20152-1	50 uL
		20152	500 uL
		20152-1mg	1 mg (lyophilized)
CF@633	630/650 nm	20125-1	50 uL
		20125	500 uL
		20125-1mg	1 mg (lyophilized)
CF@640R	642/662 nm	20178-1	50 uL
		20178	500 uL
		20178-1mg	1 mg (lyophilized)
CF@647	650/665 nm	20047-1	50 uL
		20047	500 uL
		20047-1mg	1 mg (lyophilized)

CF® Dye	Ex/Em	Catalog No.	Size
CF@660R	663/682 nm	20389-50uL	50 uL
		20389-500uL	500 uL
		20389-1mg	1 mg (lyophilized)
CF@680	681/698 nm	20418-50uL	50 uL
		20418-250uL	250 uL
CF@680R	680/701 nm	20195-1	50 uL
		20195	500 uL
CF@750	755/777 nm	20298-1	50 uL
		20298	250 uL
		20484-50uL	50 uL
CF@770	770/797 nm	20484-250uL	250 uL
		20344-50uL	50 uL
CF@790	784/806 nm	20344-50uL	50 uL
CF@800	797/816 nm	20833-50uL	50uL
Biotin	N/A	20191-1	50 uL
		20191	500 uL
		20191-1mg	1 mg (lyophilized)
AP	N/A	20467-100uL	100 uL
		20467-1mL	1 mL
HRP	N/A	20405-100uL	100 uL
		20405-1mL	1 mL
		20405-1mg	1 mg (lyophilized)

Bv: bovine; Ch: chicken; Gt: goat; GP: guinea pig; Hs: horse; Hu: human; Ms: mouse; Rt: rat; Sh: sheep; SHm: Syrian hamster

### Concentration

#### Liquid format:

CFdye and Biotin: 2 mg/mL in pH ~7.4 PBS containing 50% glycerol, 2 mg/ml bovine serum albumin (IgG-free and protease-free) and 0.05% sodium azide.

HRP: 1 mg/mL in pH ~7.4 PBS containing 50% glycerol and 15 mg/ml bovine serum albumin (IgG-free and protease-free).

#### Lyophilized format (after reconstitution):

CFdye and Biotin: 2 mg/mL in pH ~7.4 PBS containing 15 mg/mL bovine serum albumin (IgG-free and protease-free) and 20 mg/mL trehalose.

HRP: 1 mg/mL in pH ~7.4 PBS containing and 10 mg/mL trehalose.

## Storage and Handling

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.

Reconstitution (lyophilized format only):

CF® Dye and biotin conjugates: add 0.5 mL dH<sub>2</sub>O

HRP conjugate: 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 at -20°C and protect fluorescent conjugates 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.

## Product Description

Donkey anti-rabbit IgG (H+L), whole antibody, highly cross-adsorbed is an affinity purified donkey antibody that recognizes rabbit IgG. This antibody is highly-cross adsorbed for multiple labeling, and has minimal cross-reactivity with bovine, chicken, goat, guinea pig, Syrian hamster, horse, human, mouse, rat, and sheep serum proteins.

CF® dyes are superior to Alexa Fluor®, DyLight®, and Cy® dyes for antibody labeling by having combined advantages in brightness, photostability, specificity and novel features ideal for in vivo imaging. For more information about CF® dyes, download our CF® Dye Selection Guide at [www.biotium.com](http://www.biotium.com).

## Recommended Dilution Range

Fluorescence microscopy: 1-2 ug/mL

Flow cytometry: 1 ug/10<sup>6</sup> cells

Near-infrared western detection: 50-100 ng/mL

These concentrations are provided as a starting point for optimization; appropriate dilutions should be determined empirically. Generally IgG conjugates are used in the range of 1-10 ug/mL.

For more detailed protocols for immunofluorescence staining for fluorescence microscopy, flow cytometry, or fluorescence-based western detection, please visit our website.

## Related Products

Cat. No.	Product Name
23001	EverBrite™ Mounting Medium
23002	EverBrite™ Mounting Medium with DAPI
23003	EverBrite™ Hardset Mounting Medium
23004	EverBrite™ Hardset Mounting Medium with DAPI
23005	CoverGrip™ Coverslip Sealant
23007	TrueBlack™ Lipofuscin Autofluorescence Quencher
40061	RedDot™2 Far Red Nuclear Counterstain
22005	Mini Super <sup>HT</sup> Pap Pen 2.5 mm tip, ~400 uses
22006	Super <sup>HT</sup> Pap Pen 4 mm tip, ~800 uses
22003	Mini-Cell Scrapers
23006	Flow Cytometry Fixation/Permeabilization Kit
22015	Fixation Buffer
22016	Permeabilization Buffer
22017	Permeabilization and Blocking Buffer
22010	10X Fish Gelatin Blocking Agent
22011	Fish Gelatin Powder
22014	30% Bovine Serum Albumin Solution
22002	Tween®-20
22020	10X Phosphate-Buffered Saline (PBS)
22023	Paraformaldehyde, 4% in PBS, Ready-To-Use Fixative

Please visit [www.biotium.com](http://www.biotium.com) to view our full selection of products featuring bright and photostable fluorescent CF® dyes, including Mix-n-Stain™ antibody labeling kits, primary antibody conjugates, streptavidin, phalloidin, and other bioconjugates, as well as conjugates of biotin, HRP, AP, R-PE, APC and PerCP.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.

CF dye technology is covered by pending US and international patents. Alexa Fluor is a registered trademark of Molecular Probes; CYDYE is a registered trademark of GE Healthcare; DraQ7 is a registered trademark of Biostatus Ltd; DyLight is a registered trademark of Pierce Biotechnology; TWEEN is a registered trademark of Uniqema Americas LLC.