

# Product Information

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

CF® Dye	Ex/Em	Catalog No.	Size
CF@350	347/448 nm	20275-1	50 uL
		20275	500 uL
		20275-1mg	1 mg (lyophilized)
CF@488A	490/515 nm	20166-1	50 uL
		20166	500 uL
		20166-1mg	1 mg (lyophilized)
CF@543	541/560 nm	20310-1	50 uL
		20310	500 uL
		20310-1mg	1 mg (lyophilized)
CF@594	593/614 nm	20167-1	50 uL
		20167	500 uL
		20167-1mg	1 mg (lyophilized)
CF@633	630/650 nm	20168-1	50 uL
		20168	500 uL
		20168-1mg	1 mg (lyophilized)

Bv: bovine; Gt: goat; GP: guinea pig; Hs: horse; Hu: human; Ms: mouse; Rb: rabbit; Rt: rat; Sh: Sheep; SHm: Syrian hamster

### Concentration

#### Liquid format:

CFdye: 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.

#### Lyophilized format (after reconstitution):

CFdye: 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.

### 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): add 0.5 mL dH<sub>2</sub>O and mix gently to dissolve. Store at -20°C, protected from light. Aliquot to avoid freeze-thaw cycles. Alternatively, glycerol can be added to the antibody so that it will not freeze at -20°C: add 0.25 mL dH<sub>2</sub>O to the lyophilized antibody and mix gently to dissolve, then add 0.25 mL glycerol and mix well. Optional: a preservative may be added, such as 0.05% (final concentration) sodium azide.

Note: storage of the antibody for more than a day at final working dilution is not recommended.

### Product Description

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

### Accessory products for immunofluorescence

Cat. No.	Product	Features
23001	EverBrite™ Mounting Medium	<ul style="list-style-type: none"> <li>• Excellent protection from photobleaching</li> <li>• Compatible with a wide variety of dyes including cyanine dyes</li> <li>• Available in wet-set or hardset formulations</li> <li>• With or without DAPI</li> </ul>
23002	EverBrite™ Mounting Medium with DAPI	
23003	EverBrite™ Hardset Mounting Medium	
23004	EverBrite™ Hardset Mounting Medium with DAPI	
23005	CoverGrip™ Coverslip Sealant	<ul style="list-style-type: none"> <li>• Replaces nail polish for coverslip sealing</li> <li>• Won't mix with aqueous mounting media</li> </ul>
23007	TrueBlack™ Lipofuscin Autofluorescence Quencher	<ul style="list-style-type: none"> <li>• Eliminates lipofuscin autofluorescence with less background than Sudan Black B</li> <li>• Reduces background from other sources</li> <li>• Can be used before or after IF staining</li> </ul>
40061	RedDot™2 Far Red Nuclear Counterstain	<ul style="list-style-type: none"> <li>• Far-red nuclear dye for the Cy@5 channel</li> <li>• More specific than Draq@7</li> </ul>
22005	Mini Super <sup>HT</sup> Pap Pen 2.5 mm tip, ~400 uses	<ul style="list-style-type: none"> <li>• Create hydrophobic barriers around tissue sections</li> <li>• Heat-stable to 120°C</li> <li>• Insoluble in aqueous buffers, detergents, alcohol and acetone; can be removed with xylene</li> </ul>
22006	Super <sup>HT</sup> Pap Pen 4 mm tip, ~800 uses	
22003	Mini-Cell Scrapers	<ul style="list-style-type: none"> <li>• Collect cells from 96-well to 24-well plates</li> <li>• Polypropylene, sterile, disposable</li> </ul>
23006	Flow Cytometry Fixation/Permeabilization Kit	<ul style="list-style-type: none"> <li>• Ready-to-use fix/perm buffers for intracellular staining</li> </ul>
22015	Fixation Buffer	<ul style="list-style-type: none"> <li>• Convenient buffers and blocking agents for immunofluorescence or western</li> </ul>
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	

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.