

## XFD350 Goat Anti-human IgG (H+L) Antibody \*Cross Adsorbed\*

Catalog Number: 50154, 50155

Unit Size: 200 ug, 1 mg

### Product Details

---

Storage Conditions	2-6°C and kept from light. To extend the shelf-life of this product, add an equal volume of glycerol to make a final concentration of approximately 50% glycerol and store at -20°C.
Expiration Date	12 months upon receiving
Concentration	1 mg/mL
Formulation	PBS, 2 mg/mL BSA

### Antibody Properties

---

Species Reactivity	Human
Class	Secondary
Clonality	Polyclonal
Host	Goat

### Biological Properties

---

Stabilizer	None
Appearance	Off-white solid
Preparation	'Goat anti-human IgG (H+L) is produced in goat with pooled total human IgG, and affinity purified with human IgG coupled beads. The purified IgG has a minimal cross-reaction to human, horse, mouse, rabbit and bovine IgG. The antibody is conjugated with XFD350 under optimal condition.
Soluble In	Water
Application	Flow Cytometry (FACS), ELISA, HC, Western Blot
Recommended Dilutions	For IF, the suggested staining concentration is at 75-750 ng/ml. For FACS, the suggested concentration is at 500 ng-5 µg/million cells in 1 mL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined. <i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i>

### Spectral Properties

---

Conjugate	XFD350
Excitation	343 nm
Wavelength	
Emission Wavelength	441 nm

## Applications

---

AAT Bioquest's anti-human secondary antibodies have well-characterized specificity for human immunoglobulins and are useful in the detection, sorting or purification of its specified target. This Alexa Fluor® 350-labeled secondary antibody was prepared using AAT Bioquest's proprietary labeling technology. It demonstrated much brighter signal compared to the similar Alexa Fluor® 350 goat anti-human IgG antibodies from other commercial sources, and thus can significantly increase assay sensitivities. Secondary antibodies offer increased versatility enabling users to use many detection systems (e.g. HRP, AP, fluorescence). They can also provide greater sensitivity through signal amplification as multiple secondary antibodies can bind to a single primary antibody. This antibody was purified through affinity chromatography and conjugated to XFD350 (ex/em = 343/441 nm). XFD350 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 350 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 355 nm laser and 450/50 nm bandpass filter (for example, as in the BD Special Order LSRFortessa™ Cell Analyzer).