

# XFD647 Goat Anti-human IgG (H+L) Antibody

Catalog Number: 50168, 50169  
Unit Size: 200 ug, 1 mg

## Product Details

---

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	1 mg/mL
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 2 mg/mL BSA

## Unit Details

---

Reconstitution Volume	<b>50168 (200 ug)</b>	<b>50169 (1 mg)</b>
	200 uL ddH <sub>2</sub> O	1 mL ddH <sub>2</sub> O

## Antibody Properties

---

Species Reactivity	Human
Class	Secondary
Clonality	Polyclonal
Host	Goat

## Biological Properties

---

Stabilizer	2 mg/mL BSA
Appearance	Solid
Preparation	Goat anti-human IgG (H+L) is produced in goat with pooled total human IgG. The antibody is conjugated with XFD647 under optimal conditions.
Application	Flow Cytometry (FACS), IF, IHC, ELISA, WB
Recommended Dilutions	Suggested dilutions are only guidelines; users should titrate the product for their specific assay using appropriate controls

Application	Recommended dilution
Flow Cytometry (FACS)	1-5 µg/mL
IF	2 µg/mL

IHC	1-10 µg/mL
ELISA	100 ng/mL
WB	1-10 µg/mL

## Spectral Properties

---

Conjugate	Alexa Fluor® 647
Excitation Wavelength	650 nm
Emission Wavelength	671 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® 647-labeled secondary antibody was prepared using AAT Bioquest's proprietary labeling technology. It demonstrated much brighter signal compared to the similar Alexa Fluor® 647 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 XFD647 (ex/em = 650/671 nm). XFD647 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 647 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 640 nm laser and 670/14 nm bandpass filter (for example, as in the BD LSRFortessa™ Cell Analyzer).