

## XFD488 Goat Anti-human IgG (H+L) Antibody

Catalog Number: 50156, 50157

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       |

### Chemical Properties

---

|                  |         |
|------------------|---------|
| Molecular Weight | ~150000 |
|------------------|---------|

### Biological Properties

---

|                       |                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                       |
|-----------------------|-------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Stabilizer            | None                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                  |
| Appearance            | Orange 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 antibody is conjugated with XFD488 under optimal condition.                                                                                                                                                                                                                                                                                                       |
| Soluble In            | Water                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                 |
| Application           | Flow Cytometry (FACS), ELISA, HC, Western Blot                                                                                                                                                                                                                                                                                                                                                                                                                                                        |
| Recommended Dilutions | <p>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.</p> <p><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></p> |

## Spectral Properties

---

|                     |        |
|---------------------|--------|
| Conjugate           | XFD488 |
| Excitation          |        |
| Wavelength          | 499 nm |
| Emission Wavelength | 520 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® 488-labeled secondary antibody was prepared using AAT Bioquest's proprietary labeling technology. It demonstrated much brighter signal compared to the similar Alexa Fluor® 488 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 XFD488 (ex/em = 499/520 nm). XFD488 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 488 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 488 nm laser and 530/40 nm bandpass filter (for example, as in the BD FACSJazz™).