

**PE/AF 700 Goat Anti-human IgG (H+L)
Antibody *Cross Adsorbed***Catalog number: 50238, 50239
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 |

Unit Details

| | | |
|-----------------------|---------------------------|-------------------------|
| Unit | 50238 (200 ug) | 50239 (1 mg) |
| Reconstitution Volume | 200 uL ddH ₂ O | 1 mL ddH ₂ O |

Antibody Properties

| | |
|--------------------|------------|
| Species Reactivity | Human |
| Class | Secondary |
| Clonality | Polyclonal |
| Host | Goat |

Biological Properties

| | |
|-------------|--|
| Stabilizer | None |
| 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 PE/AF 700 under optimal condition. |
| Application | Flow Cytometry (FACS), ELISA, HC, Western Blot |
| Soluble In | Water |

Spectral Properties

| | |
|-----------------------|-----------|
| Conjugate | PE/AF 700 |
| Excitation Wavelength | 566 nm |
| Emission Wavelength | 721 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 PE/Alexa Fluor® 700-labeled secondary antibody was prepared using AAT Bioquest's proprietary labeling technology. It demonstrated much brighter signal compared to the similar PE/Alexa Fluor® 700 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 PE/Alexa Fluor® 700 (ex/em = 566/721 nm). It is compatible with the 561 nm laser and 710/50 nm bandpass filter (for example, as in the BD Special Order LSRFortessa™ Cell Analyzer).