

## iFluor® 594 Goat Anti-human IgG (H+L) Antibody \*Cross Adsorbed\*

Catalog Number: 50086, 50087

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>50086 (200 ug)</b>     | <b>50087 (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 and affinity purified with human IgG coupled beads. The antibody is conjugated with iFluor® 594 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

Flow Cytometry (FACS)

#### Recommended dilution

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             | iFluor™ 594 |
| Excitation Wavelength | 587 nm      |
| Emission Wavelength   | 603 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 iFluor® 594-labeled secondary antibody was prepared using AAT Bioquest's proprietary labeling technology. It demonstrated much brighter signal compared to the similar iFluor® 594 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 iFluor® 594 (ex/em = 588/604 nm). It is compatible with the 561 nm laser and 585/29 nm bandpass filter (for example, as in the BD FACSJazz™).