

**iFluor® 350 goat anti-rabbit IgG (H+L)**

Catalog Number: 16600, 16795

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>16600 (200 ug)</b>     | <b>16795 (1 mg)</b>     |
|                       | 200 uL ddH <sub>2</sub> O | 1 mL ddH <sub>2</sub> O |

**Antibody Properties**

---

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

**Chemical Properties**

---

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

**Biological Properties**

---

|                       |                                                                                                                                                    |
|-----------------------|----------------------------------------------------------------------------------------------------------------------------------------------------|
| Stabilizer            | 2 mg/mL BSA                                                                                                                                        |
| Appearance            | Solid                                                                                                                                              |
| Preparation           | Goat anti-rabbit IgG (H+L) is produced in goat with pooled total rabbit IgG. The antibody is conjugated with iFluor® 350 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             | iFluor® 350 |
| Excitation Wavelength | 345 nm      |
| Emission Wavelength   | 450 nm      |

## Applications

---

AAT Bioquest's iFluor® dyes are optimized for labeling proteins, in particular, antibodies. These dyes are bright, photostable and have minimal quenching on proteins. They can be well excited by the major laser lines of fluorescence instruments (e.g., 350, 405, 488, 555 and 633 nm). iFluor® 350 goat anti-rabbit IgG (H+L) conjugate has fluorescence excitation and emission maxima of ~345 nm and ~442 nm respectively. These spectral characteristics make them an excellent alternative to Alexa Fluor® 350 goat anti-rabbit IgG (H+L) conjugate (Alexa Fluor® is the trademark of Invitrogen).