

**iFluor® 633 goat anti-mouse IgG (H+L) \*Cross Adsorbed\***

Catalog Number: 16558, 16782

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>16558 (200 ug)</b>     | <b>16782 (1 mg)</b>     |
|                       | 200 uL ddH <sub>2</sub> O | 1 mL ddH <sub>2</sub> O |

**Antibody Properties**

---

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

**Chemical Properties**

---

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

**Biological Properties**

---

|                       |                                                                                                                                                                                                     |
|-----------------------|-----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Stabilizer            | 2 mg/mL BSA                                                                                                                                                                                         |
| Appearance            | Solid                                                                                                                                                                                               |
| Preparation           | Goat anti-mouse IgG (H+L) is produced in goat with pooled total mouse IgG and affinity purified with mouse IgG coupled beads. The antibody is conjugated with iFluor® 633 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® 633 |
| Excitation Wavelength | 640 nm      |
| Emission Wavelength   | 654 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® 633 goat anti-mouse IgG (H+L) conjugate has fluorescence excitation and emission maxima of ~638 nm and ~655 nm respectively. These spectral characteristics make them an excellent alternative to Alexa Fluor® 633 goat anti-mouse IgG (H+L) conjugate (Alexa Fluor® is the trademark of Invitrogen).