

**iFluor® 594 goat anti-mouse IgG (H+L)**

Catalog Number: 16468, 16741

Unit Size: 200 µg, 1 mg

**Product Details**

---

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
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	200 µL ddH <sub>2</sub> O for 200 µg (#16468) and 1 mL ddH <sub>2</sub> O for 1 mg (#16741)
-----------------------	---

**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. 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	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® 594
Excitation Wavelength	587 nm
Emission Wavelength	603 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® 594 goat anti-mouse IgG (H+L) conjugate has fluorescence excitation and emission maxima of ~592 nm and ~614 nm respectively. These spectral characteristics make them an excellent alternative to Alexa Fluor® 594 goat anti-mouse IgG (H+L) conjugate (Alexa Fluor® is the trademark of Invitrogen).