

**Texas Red® goat anti-mouse IgG (H+L)**

Catalog Number: 16853

Unit Size: 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	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. The antibody is conjugated with Texas Red® 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	Texas Red®
Excitation Wavelength	586 nm
Emission Wavelength	603 nm

### Applications

---

AAT Bioquest's anti-mouse secondary antibodies are affinity-purified antibodies with well-characterized specificity for mouse immunoglobulins and are useful in the detection, sorting or purification of its specified target. This Texas Red-labeled secondary antibody was prepared using AAT Bioquest's proprietary labeling technology. It demonstrated much brighter signal compared to the similar Texas Red goat anti-mouse IgG antibodies from other commercial sources, 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.