

## PE/Texas Red® Goat Anti-Mouse IgG (H+L) Antibody

Catalog Number: 50312

Unit Size: 200 ug

### Product Details

---

|                    |   |
|--------------------|---|
| Storage Conditions | 2-8°C with minimized light exposure. Do not freeze.                         |
| Expiration Date    | 12 months upon receiving  |
| Concentration      | ~0.2 mg/mL  |
| Formulation        | Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide, 0.2% (w/v) BSA |

### Antibody Properties

---

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

### Biological Properties

---

|                       |  |
|-----------------------|--|
| Stabilizer            | 0.09% sodium azide, 0.2% (w/v) BSA   |
| Appearance            | Liquid   |
| Preparation           | Goat anti-mouse IgG (H+L) is produced in goat with pooled total mouse IgG. The antibody is conjugated with PE/Texas Red® under optimal conditions. |
| Application           | Flow Cytometry (FACS), IF, IHC   |
| 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           |

### Spectral Properties

---

|           |                              |
|-----------|------------------------------|
| Conjugate | PE (phycoerythrin)-Texas Red |
|-----------|------------------------------|

Excitation Wavelength      565 nm

Emission Wavelength      613 nm

## **Applications**

---

PE/Texas Red® Goat Anti-Mouse IgG (H+L) Antibody is a high-performance secondary antibody designed to bind both heavy and light chains of mouse immunoglobulin G (IgG). Conjugated to PE/Texas Red, this antibody delivers exceptionally bright red fluorescence with excitation/emission maxima of ~565/613 nm. The PE/Texas Red® fluorophore is prepared using AAT Bioquest's proprietary labeling technology to ensure superior signal intensity compared to similar antibodies from other sources. This antibody is ideal for flow cytometry, fluorescence microscopy, ELISA, and other immunoassays requiring sensitive and specific detection of mouse IgG. Its compatibility with 561 nm yellow-green laser lines and common bandpass filters (e.g., 613/18 nm) ensures seamless integration into standard fluorescence detection workflows.