

iFluor® 800 goat anti-mouse IgG (H+L)

Catalog Number: 48000, 48001

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 | 48000 (200 ug) | 48001 (1 mg) |
| | 200 uL ddH ₂ O | 1 mL ddH ₂ 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 iFluor® 800 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® 800 |
| Excitation Wavelength | 801 nm |
| Emission Wavelength | 820 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, 532-561, 633-647 and 808 nm). iFluor® 800 goat anti-mouse IgG (H+L) conjugate has fluorescence excitation and emission maxima of 801 nm and 820 nm, respectively. These spectral characteristics make them an excellent alternative to Alexa Fluor® 800 goat anti-mouse IgG (H+L) conjugate (Alexa Fluor® is a trademark of Invitrogen).