

iFluor™ 405 goat anti-mouse IgG (H+L) *Cross Adsorbed*

Catalog number: 16524, 16771 Unit size: 200 ug, 1 mg

| Product Details | | | |
|------------------------------|---|-------------------------|--|
| Storage Conditions | 2-6°C and kept from light. To extend the shelf-life of this product, add an equal volume of glycerol to make a final concentration of approximately 50% glycerol and store at -20°C. | | |
| Expiration Date | 12 months upon receiving | | |
| Concentration | 1 mg/mL | | |
| Formulation | PBS, 2 mg/mL BSA | | |
| Unit Details | | | |
| Unit | 16524 (200 ug) | 16771 (1 mg) | |
| Reconstitution Volume | 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 | None | | |
| Appearance | Light yellow 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 purified IgG has a minimal cross-reaction to human, horse, rabbit and bovine IgG. The antibody is conjugated with iFluor™ 405 under optimal condition. | | |
| Application | Immunofluorescence (IF), Flow Cytometry (FACS) | | |
| Soluble In | Water | | |
| Spectral Properties | | | |
| Conjugate | iFluor™ 405 | | |

| Excitation Wavelength | 403 nm |
|-----------------------|--------|
| Emission Wavelength | 427 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[™] 405 goat anti-mouse IgG (H+L) conjugate has fluorescence excitation and emission maxima of ~401 nm and ~420 nm respectively. These spectral characteristics make them an excellent alternative to Alexa Fluor[®] 405 goat anti-mouse IgG (H+L) conjugate (Alexa Fluor[®] is the trademark of Invitrogen).