

Product Information Sheet

Ordering Information

| | |
|---------------------|--|
| Product Number: | 16775 |
| Product Name: | iFluor™ 532 Goat Anti-mouse (H+L) IgG conjugate *Cross Adsorbed* |
| Unit Size: | 1 mg |
| Concentration: | 1 mg/mL |
| Formulation: | PBS, 2 mg/mL BSA |
| Stabilizer: | None |
| Appearance: | Red powder |
| Reconstitution: | In 1 mL ddH ₂ O |
| 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. |

Biological and Spectral Properties

| | |
|------------------------|--|
| Preparation: | Goat anti-mouse (H&L) IgG *Cross Adsorbed* 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™ 532 under optimal condition. |
| Excitation Wavelength: | 542 nm |
| Emission Wavelength: | 558 nm |
| Soluble in: | Aqueous buffers |

Applications: Immunofluorescence (IF), Flow Cytometry (FACS)

Recommended dilutions: For IF, the suggested staining concentration is at 75-750 ng/ml. For FACS, the suggested concentration is at 500 ng-5 µg/million cells in 1 mL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined.

**The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.*