

## Product Information Sheet

### **Ordering Information**

Product Number:	16788
Product Name:	iFluor™ 750 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:	Dark blue powder
Reconstitution:	In 1 mL ddH <sub>2</sub> 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™ 750 under optimal condition.
Excitation Wavelength:	753 nm
Emission Wavelength:	779 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.*