

Product Information Sheet

Ordering Information

Product Number:	16782
Product Name:	iFluor™ 633 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:	Black 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™ 633 under optimal condition.
Excitation Wavelength:	638 nm
Emission Wavelength:	655 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.*