

## Product Information Sheet

### **Ordering Information**

Product Number:	16690
Product Name:	iFluor <sup>™</sup> 555 Goat Anti-rabbit (H+L) IgG conjugate *Cross Adsorbed*
Unit Size:	200 µg
Concentration:	1 mg/mL
Formulation:	PBS, 2 mg/mL BSA
Stabilizer:	None
Appearance:	Red powder
Reconstitution:	In 200 µL 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-rabbit (H&L) IgG is produced in goat with pooled total rabbit IgG, and affinity purified with rabbit IgG coupled beads. The purified IgG has a minimal cross-reaction to human, horse, mouse and swine IgG. The antibody is conjugated with iFluor <sup>™</sup> 555 under optimal condition.
Excitation Wavelength:	559 nm
Emission Wavelength:	569 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.*