

Product Information Sheet

Product Name:	iFluor™ 680 goat anti-mouse IgG (H+L) *Cross Adsorbed*	
Catalog Number:	16566	16784
Unit Size:	200 ug	1 mg
Concentration:	1 mg/mL	1 mg/mL
Formulation:	PBS, 2 mg/mL BSA	PBS, 2 mg/mL BSA
Reconstitution Volume:	200 uL ddH ₂ O	1 mL ddH ₂ O

Biological and Spectral Properties

Stabilizer:	None
Appearance:	Blue
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.
Preparation:	Goat anti-mouse (H&L) IgG 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™ 680 under optimal condition.
Excitation Wavelength:	682
Emission Wavelength:	701
Soluble in:	Water
<u>Applications:</u>	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.*