

## Product Information Sheet

Product Name:	Cy7® goat anti-rabbit IgG (H+L) *Cross Adsorbed*
Catalog Number:	16880
Unit Size:	1 mg
Concentration:	1 mg/mL
Formulation:	PBS, 2 mg/mL BSA
Reconstitution Volume:	1 mL ddH <sub>2</sub> O

### **Biological and Spectral Properties**

Stabilizer:	None
Appearance:	Dark 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-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 bovine IgG. The antibody is conjugated with Cy7® under optimal condition.
Excitation Wavelength:	749
Emission Wavelength:	776
Soluble in:	Water

**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.*