

**iFluor™ 820 Goat Anti-human IgG (H+L)  
Antibody**Catalog number: 50136, 50137  
Unit size: 200 ug, 1 mg**Product Details**

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.
Expiration Date	12 months upon receiving
Concentration	1 mg/mL
Formulation	PBS, 2 mg/mL BSA

**Unit Details**

Unit	50136 (200 ug)	50137 (1 mg)
Reconstitution Volume	200 uL ddH <sub>2</sub> O	1 mL ddH <sub>2</sub> O

**Antibody Properties**

Species Reactivity	Human
Class	Secondary
Clonality	Polyclonal
Host	Goat

**Biological Properties**

Stabilizer	None
Appearance	Green solid
Preparation	Goat anti-human IgG (H+L) is produced in goat with pooled total human IgG, and affinity purified with human IgG coupled beads. The antibody is conjugated with iFluor™ 820 under optimal condition.
Application	Flow Cytometry (FACS), ELISA, HC, Western Blot
Soluble In	Water

**Spectral Properties**

Conjugate	iFluor™ 820
Excitation Wavelength	822 nm
Emission Wavelength	850 nm

## Applications

---

AAT Bioquest's anti-human secondary antibodies have well-characterized specificity for human immunoglobulins and are useful in the detection, sorting or purification of its specified target. This iFluor™ 820-labeled secondary antibody was prepared using AAT Bioquest's proprietary labeling technology. It demonstrated much brighter signal compared to the similar iFluor™ 820 goat anti-human IgG antibodies from other commercial sources, and thus can significantly increase assay sensitivities. Secondary antibodies offer increased versatility enabling users to use many detection systems (e.g. HRP, AP, fluorescence). They can also provide greater sensitivity through signal amplification as multiple secondary antibodies can bind to a single primary antibody. This antibody was purified through affinity chromatography and conjugated to iFluor™ 820 (ex/em = 822/850 nm).