

## iFluor<sup>™</sup> 810 goat anti-rabbit IgG (H+L)

Catalog number: 48054, 48055 Unit size: 200 µg, 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	48054 (200 μg)	48055 (1 mg)
Reconstitution Volume	$200~\mu L~ddH_2O$	1 mL ddH <sub>2</sub> O
Antibody Properties		
Species Reactivity	Rabbit	
Class	Secondary	
Clonality	Polyclonal	
Host	Goat	
Chemical Properties		
Molecular Weight	~150000	
Biological Properties		
Stabilizer	None	
Appearance	Green solid	
Preparation	Goat anti-rabbit IgG (H+L) is produced in goat with pooled total rabbit IgG, and affinity purified with rabbit IgG coupled beads. The antibody is conjugated with iFluor™ 810 under optimal condition.	
Application	Flow Cytometry (FACS), ELISA, HC, Western Blot	
Soluble In	Water	
Spectral Properties		
Conjugate	iFluor™ 810	
Excitation Wavelength	811 nm	

## Applications

AAT Bioquest's iFluor<sup>™</sup> dyes are optimized for labeling proteins, in particular, antibodies. These dyes are bright, photostable, and have minimal quenching on proteins. They can be well excited by the major laser lines of fluorescence instruments (e.g., 350, 405, 488, 532-561, 633-647, and 808 nm). iFluor<sup>™</sup> 810 goat anti-rabbit IgG (H+L) conjugate has fluorescence excitation and emission maxima of 🛛 811 nm and 🖾 822 nm, respectively. These unique spectral characteristics makes iFluor<sup>™</sup> 810 goat anti-rabbit IgG (H+L) conjugate har stop anti-rabbit IgG (H+L) conjugates ideal for various NIR imaging applications, including Western blotting, ELISA, protein arrays, tissue section imaging, and *in vivo* imaging.