

## iFluor™ 514 goat anti-rabbit IgG (H+L)

Catalog number: 16612, 16801 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	16612 (200 ug)	16801 (1 mg)
Reconstitution Volume	200 uL ddH <sub>2</sub> O	1 mL ddH <sub>2</sub> O
Antibody Properties		
Species Reactivity	Rabbit	
Class	Secondary	
Clonality	Polyclonal	
Host	Goat	
Chemical Properties		
Molecular Weight	~150000	
<b>Biological Properties</b>		
Stabilizer	None	
Appearance	Purple 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™ 514 under optimal condition.	
Application	Immunofluorescence (IF), Flow Cytometry (FACS)	
Soluble In	Water	
Spectral Properties		
Conjugate	iFluor™ 514	
Excitation Wavelength	528 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, 555 and 633 nm). iFluor<sup>™</sup> 514 goat anti-rabbit IgG (H+L) conjugate has fluorescence excitation and emission maxima of ~518 nm and ~542 nm respectively. These spectral characteristics make them an excellent alternative to Alexa Fluor<sup>®</sup> 514 goat anti-rabbit IgG (H+L) conjugate (Alexa Fluor<sup>®</sup> is the trademark of Invitrogen).