

## iFluor™ 633 goat anti-rabbit IgG (H+L) \*Cross Adsorbed\*

Catalog number: 16704, 16835 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	16704 (200 ug)	16835 (1 mg)	
Reconstitution Volume	200 uL $ddH_2O$	1 mL ddH <sub>2</sub> O	
Antibody Properties			
Species Reactivity	Rabbit		
Class	Secondary		
Clonality	Polyclonal		
Host	Goat		
Chemical Properties			
Molecular Weight	~150000		
<b>Biological Properties</b>			
Stabilizer	None		
Appearance	Black 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 purified IgG has a minimal cross-reaction to human, horse, mouse and bovine IgG. The antibody is conjugated with iFluor™ 633 under optimal condition.		
Application	Immunofluorescence (IF), Flow Cytometry (FACS)		
Soluble In	Water		
Spectral Properties			
Conjugate	iFluor™ 633		

Excitation Wavelength	640 nm
Emission Wavelength	654 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> 633 goat anti-rabbit IgG (H+L) conjugate has fluorescence excitation and emission maxima of ~638 nm and ~655 nm respectively. These spectral characteristics make them an excellent alternative to Alexa Fluor<sup>®</sup> 633 goat anti-rabbit IgG (H+L) conjugate (Alexa Fluor<sup>®</sup> is the trademark of Invitrogen).