

iFluor[™] 405 goat anti-mouse IgG (H+L)

Catalog number: 16444, 16731 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	16444 (200 ug)	16731 (1 mg)	
Reconstitution Volume	200 uL ddH ₂ O	1 mL ddH ₂ O	
Antibody Properties			
Species Reactivity	Mouse		
Class	Secondary		
Clonality	Polyclonal		
Host	Goat		
Chemical Properties			
Molecular Weight	~150000		
Biological Properties			
Stabilizer	None		
Appearance	Light yellow solid		
Preparation	Goat anti-mouse IgG (H+L) is produced in goat with pooled total mouse IgG, and affinity purified with mouse IgG coupled beads. The antibody is conjugated with iFluor™ 405 under optimal condition.		
Application	Immunofluorescence (IF), Flow Cytometry (FACS)		
Soluble In	Water		
Spectral Properties			
Conjugate	iFluor™ 405	iFluor™ 405	
Excitation Wavelength	403 nm		

Applications

AAT Bioquest's iFluor[™] 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[™] 405 goat anti-mouse IgG (H+L) conjugate has fluorescence excitation and emission maxima of ~401 nm and ~420 nm respectively. These spectral characteristics make them an excellent alternative to Alexa Fluor[®] 405 goat anti-mouse IgG (H+L) conjugate (Alexa Fluor[®] is the trademark of Invitrogen).