

iFluor™ 350 goat anti-mouse IgG (H+L)
Cross AdsorbedCatalog number: 16520, 16770
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	16520 (200 ug)	16770 (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	Off-white 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 purified IgG has a minimal cross-reaction to human, horse, rabbit and bovine IgG. The antibody is conjugated with iFluor™ 350 under optimal condition.
Application	Immunofluorescence (IF), Flow Cytometry (FACS)
Soluble In	Water

Spectral Properties

Conjugate	iFluor™ 350
-----------	-------------

Excitation Wavelength	345 nm
Emission Wavelength	450 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™ 350 goat anti-mouse IgG (H+L) conjugate has fluorescence excitation and emission maxima of ~345 nm and ~442 nm respectively. These spectral characteristics make them an excellent alternative to Alexa Fluor® 350 goat anti-mouse IgG (H+L) conjugate (Alexa Fluor® is the trademark of Invitrogen).