

iFluor® 860 goat anti-mouse IgG (H+L) *Cross Adsorbed*

Catalog Number: 48018, 48019

Unit Size: 200 ug, 1 mg

Product Details

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	1 mg/mL
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 2 mg/mL BSA

Unit Details

Reconstitution Volume	48018 (200 ug)	48019 (1 mg)
	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	2 mg/mL BSA
Appearance	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® 860 under optimal conditions.
Application	Flow Cytometry (FACS), IF, IHC, ELISA, WB
Recommended Dilutions	Suggested dilutions are only guidelines; users should titrate the product for their specific assay using appropriate controls

Application	Recommended dilution
Flow Cytometry (FACS)	1-5 µg/mL
IF	2 µg/mL
IHC	1-10 µg/mL
ELISA	100 ng/mL
WB	1-10 µg/mL

Spectral Properties

Conjugate	iFluor® 860
Excitation Wavelength	853 nm
Emission Wavelength	878 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, 532-561, 633-647, and 808 nm). iFluor® 860 goat anti-mouse IgG (H+L) conjugate has fluorescence excitation and emission maxima of 853 nm and 878 nm, respectively. These unique spectral characteristics makes iFluor® 860 goat anti-mouse IgG (H+L) conjugates ideal for various NIR imaging applications, including Western blotting, ELISA, protein arrays, tissue section imaging, and *in vivo* imaging.