

iFluor® 790 goat anti-mouse IgG (H+L) *Cross Adsorbed*

Catalog Number: 16587, 16790

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	16587 (200 ug)	16790 (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® 790 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® 790
Excitation Wavelength	787 nm
Emission Wavelength	812 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® 790 goat anti-mouse IgG (H+L) conjugate has IR fluorescence excitation and emission maxima of ~780 nm and ~810 nm respectively. These spectral characteristics make them an excellent alternative to IRDye® 800 goat anti-mouse IgG (H+L) conjugate (IRDye® is the trademark of Li-COR).