

iFluor® 680 goat anti-rabbit IgG (H+L) *Cross Adsorbed*

Catalog Number: 16712, 16838

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	16712 (200 ug)	16838 (1 mg)
	200 uL ddH ₂ O	1 mL ddH ₂ O

Antibody Properties

Species Reactivity	Rabbit
Class	Secondary
Clonality	Polyclonal
Host	Goat

Chemical Properties

Molecular Weight	~150000
------------------	---------

Biological Properties

Stabilizer	2 mg/mL BSA
Appearance	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 antibody is conjugated with iFluor® 680 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® 680
Excitation Wavelength	684 nm
Emission Wavelength	701 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® 680 goat anti-rabbit IgG (H+L) conjugate has fluorescence excitation and emission maxima of ~682 nm and ~701 nm respectively. These spectral characteristics make them an excellent alternative to Alexa Fluor® 680 goat anti-rabbit IgG (H+L) conjugate (Alexa Fluor® is the trademark of Invitrogen).