

iFluor® 568 goat anti-rabbit IgG (H+L) *Cross Adsorbed*

Catalog Number: 16692, 16832

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	16692 (200 ug)	16832 (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® 568 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® 568
Excitation Wavelength	568 nm
Emission Wavelength	587 nm

Applications

iFluor® 568 is a bright red fluorescent dye. iFluor® 568-labeled anti-IgG conjugates exhibit bright fluorescence signal and good photostability. Used for stable signal generation in imaging and flow cytometry, the fluorescence intensity of iFluor® 568 conjugates is pH-insensitive from pH 4 to pH 11. The iFluor® 568-labeled antibody conjugates can be well excited with Krypton ion laser (~568 nm). iFluor® 568 family has the spectral properties essentially identical to those of Alexa Fluor® 568. Under the same conditions we tested, iFluor® 568 antibody conjugates are brighter and more photostable than the corresponding Alexa Fluor® 568. These spectral and labeling characteristics make the iFluor® 568 dye family a superior alternative to Alexa Fluor® 568. In addition, iFluor® 568 secondary antibody conjugates give higher signal/background ratios than the corresponding Alexa Fluor® 568-labeled conjugates.