

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

Catalog Number: 16541, 16777

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	<b>16541 (200 ug)</b>	<b>16777 (1 mg)</b>
	200 uL ddH <sub>2</sub> O	1 mL ddH <sub>2</sub> 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® 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.