

iFluor® 546 goat anti-rabbit IgG (H+L) *Cross Adsorbed*

Catalog Number: 16688

Unit Size: 200 ug

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	200 uL 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® 546 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® 546
Excitation Wavelength	541 nm
Emission Wavelength	557 nm

Applications

iFluor® 546 is a bright orange fluorescent dye. iFluor® 546-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® 546 conjugates is pH-insensitive from pH 4 to pH 11. The iFluor® 546-labeled antibody conjugates can be well excited with either Nd:YAG laser (~532 nm) or Helium-Neon laser (~543 nm). iFluor® 546 family has the spectral properties essentially identical to those of Alexa Fluor® 546. Under the same conditions we tested, iFluor® 546 antibody conjugates are brighter and more photostable than the corresponding Alexa Fluor® 546. These spectral and labeling characteristics make the iFluor® 546 dye family a superior alternative to Alexa Fluor® 546. In addition, iFluor® 546 secondary antibody conjugates give higher signal/background ratios than the corresponding Alexa Fluor® 546-labeled conjugates.