

AF594-labeled goat anti-rabbit IgG (H+L)***Cross Adsorbed***

Catalog number: 16404

Unit size: 1 mg

Product Details

Storage Conditions	2-6°C and kept from light. To extend the shelf-life of this product, add an equal volume of glycerol to make a final concentration of approximately 50% glycerol and store at -20°C.
Expiration Date	12 months upon receiving
Concentration	1 mg/mL
Formulation	PBS, 2 mg/mL BSA

Unit Details

Unit	16404 (1 mg)
Reconstitution Volume	1 mL ddH ₂ O

Antibody Properties

Species Reactivity	Rabbit
Class	Secondary
Clonality	Polyclonal
Host	Goat

Chemical Properties

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

Biological Properties

Stabilizer	None
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 purified IgG has a minimal cross-reaction to human, horse, mouse and bovine IgG. The antibody is conjugated with Alexa Fluor™ 594 under optimal condition.
Application	Immunofluorescence (IF), Flow Cytometry (FACS)
Soluble In	Water

Spectral Properties

Conjugate	Alexa Fluor™ 594
-----------	------------------

Excitation Wavelength	590 nm
Emission Wavelength	618 nm

Applications

Our goat anti-rabbit IgG whole antibodies have been cross-adsorbed against human IgG and human serum prior to conjugation to minimize cross-reactivity. This AF488 labeled-goat anti-rabbit IgG conjugate is prepared by the reaction of cross-adsorbed goat anti-rabbit IgG whole antibody with AF594 NHS ester, the same molecule to Alexa Fluor® 488 NHS ester (Alexa Fluor is the trademark of ThermoFisher). Each conjugate has typically 4-6 fluorophores per IgG molecule. Fluorescent secondary antibody conjugates are useful in the detection, sorting, or purification of its specified target and ideal for fluorescence microscopy and confocal laser scanning microscopy, flow cytometry, and fluorescent western detection.