

XFD594 goat anti-rabbit IgG (H+L) *Cross Adsorbed*

Catalog Number: 16404

Unit Size: 1 mg

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
--------------------	--

Expiration Date	12 months upon receiving
-----------------	--------------------------

Unit Details

Units	16404 (1 mg)
-------	--------------

Reconstitution Volume	0.2 mL ddH ₂ O
-----------------------	---------------------------

Chemical Properties

Appearance	Solid purple
------------	--------------

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

Soluble In	Water
------------	-------

Spectral Properties

Excitation Wavelength	590 nm
-----------------------	--------

Emission Wavelength	618 nm
---------------------	--------

Applications

XFD594 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 594 (Alexa Fluor® is the trademark of Thermo Fisher). 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 XFD594 labeled-goat anti-rabbit IgG conjugate is prepared by the reaction of cross-adsorbed goat anti-rabbit IgG whole antibody with XFD594 NHS ester, which has the same molecule as Alexa Fluor® 594 NHS ester. 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.