

XFD350 goat anti-mouse IgG (H+L) *Cross Adsorbed*

Catalog Number: 16380

Unit Size: 1 mg

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid
Molecular Weight	~150000
Soluble In	Water

Spectral Properties

Excitation Wavelength	343 nm
Emission Wavelength	441 nm

Applications

XFD350 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 350 (Alexa Fluor® is the trademark of Thermo Fisher). Our goat anti-mouse IgG whole antibodies have been cross-adsorbed against human IgG and human serum prior to conjugation to minimize cross-reactivity. This XFD350 labeled-goat anti-mouse IgG conjugate is prepared by the reaction of cross-adsorbed goat anti-mouse IgG whole antibody with XFD350 NHS ester, which has the same molecule as Alexa Fluor® 350 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.