

XFD488 goat anti-mouse IgG (H+L) *Cross Adsorbed*

Catalog Number: 16383

Unit Size: 1 mg

Product Details

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

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

Unit Details

Units	16383 (1 mg)
-------	--------------

Reconstitution Volume	1 mL ddH ₂ O
-----------------------	-------------------------

Chemical Properties

Appearance	Solid orange
------------	--------------

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

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

Spectral Properties

Excitation Wavelength	499 nm
-----------------------	--------

Emission Wavelength	520 nm
---------------------	--------

Applications

XFD488 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 488 (Alexa Fluor® is the trademark of Thermo Fisher). AAT Bioquest's iFluor® dyes are optimized for labeling proteins, in particular, antibodies. These dyes are bright, photostable and have minimal quenching on proteins. They can be well excited by the major laser lines of fluorescence instruments (e.g., 350, 405, 488, 555 and 633 nm). XFD488 goat anti-mouse IgG (H+L) conjugate has fluorescence excitation and emission maxima of ~491 nm and ~516 nm respectively. These spectral characteristics make them an excellent alternative to Alexa Fluor® 488 goat anti-mouse IgG (H+L) conjugate.