

iFluor® 647 goat anti-rabbit IgG (H+L)

Catalog Number: 16642, 16809

Unit Size: 200 ug, 1 mg

Product Details

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

Chemical Properties

Appearance	Solid blue
Molecular Weight	~150000
Soluble In	Water

Spectral Properties

Excitation Wavelength	656 nm
Emission Wavelength	670 nm

Applications

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). iFluor® 647 goat anti-rabbit IgG (H+L) conjugate has fluorescence excitation and emission maxima of ~654 nm and ~674 nm respectively. These spectral characteristics make them an excellent alternative to Alexa Fluor®; 647 goat anti-rabbit IgG (H+L) conjugate (Alexa Fluor® is the trademark of Invitrogen).