

iFluor™ 555 goat anti-mouse IgG (H+L)

Catalog number: 16460, 16739

Unit size: 200 ug, 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 16460 (200 ug) 16739 (1 mg)

Reconstitution Volume 200 uL ddH₂O 1 mL ddH₂O

Antibody Properties

Species Reactivity Mouse

Class Secondary

Clonality Polyclonal

Host Goat

Chemical Properties

Molecular Weight ~150000

Biological Properties

Stabilizer None

Appearance Red solid

Preparation Goat anti-mouse IgG (H+L) is produced in goat with pooled total mouse IgG, and affinity purified

with mouse IgG coupled beads. The antibody is conjugated with iFluor™ 555 under optimal

condition.

Application Immunofluorescence (IF), Flow Cytometry (FACS)

Soluble In Water

Spectral Properties

Conjugate iFluor™ 555

Excitation Wavelength 557 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™ 555 goat anti-mouse IgG (H+L) conjugate has fluorescence excitation and emission maxima of ~559 nm and ~569 nm respectively. These spectral characteristics make them an excellent alternative to Alexa Fluor® 555 goat anti-mouse IgG (H+L) conjugate (Alexa Fluor® is the trademark of Invitrogen).